



FLIPR® Penta

High-Throughput Cellular Screening System

Hardware and Software Version 5.0.1

User Guide

This document is provided to customers who have purchased Molecular Devices equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a non-transferrable license. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Each product is shipped with documentation stating specifications and other technical information. Molecular Devices products are warranted to meet the stated specifications. Molecular Devices makes no other warranties or representations express or implied, including but not limited to, the fitness of this product for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom. The sole obligation of Molecular Devices and the customer's sole remedy are limited to repair or replacement of the product in the event that the product fails to do as warranted.

For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: <http://www.moleculardevices.com/patents>

Product manufactured by Molecular Devices, LLC.
3860 N. First Street, San Jose, California, 95134, United States of America.
Molecular Devices, LLC is ISO 9001 registered.
©2020 Molecular Devices, LLC.
All rights reserved.



Contents

| | |
|--|------------|
| Safety Information | 7 |
| Warnings, Cautions, Notes, and Tips | 7 |
| Symbols on Instrument Labels | 8 |
| Electrical Safety | 9 |
| Chemical and Biological Safety | 9 |
| Moving Parts Safety | 10 |
| Ultraviolet (UV) Light Safety | 10 |
| Chapter 1: Introduction | 13 |
| Workflow | 14 |
| Before You Start | 14 |
| System Requirements | 14 |
| Software Installation | 16 |
| Powering On and Powering Off the System | 18 |
| Checking Instrument Status and Configuration | 20 |
| Chapter 2: System Hardware Overview | 23 |
| Five-Position Stage | 24 |
| Plate Support | 25 |
| Plate-Handling System | 26 |
| Liquid-Handling System | 29 |
| Optical System | 35 |
| Chiller | 45 |
| Observation Panel | 45 |
| Chapter 3: ScreenWorks Software Overview | 47 |
| Online Mode vs. Offline Mode | 47 |
| User Interface Overview | 48 |
| Using Protocols | 71 |
| Chapter 4: Calibrating and Signal Testing | 131 |
| Calibrating Optics | 131 |
| Running Signal Tests | 135 |

| | |
|--|------------|
| Chapter 5: Running an Experiment | 143 |
| Preparing Cells | 143 |
| Starting the System | 145 |
| Checking the System | 146 |
| Dye-Loading the Cells for Fluorescence Assays | 146 |
| Preparing a Source and Compound Plate | 147 |
| Setting Up an Assay Protocol | 149 |
| Optimizing the Optics and Fluid Dispensing | 150 |
| Starting the Assay Run | 155 |
| Chapter 6: Exporting Data | 157 |
| Exporting Time Sequence | 159 |
| Exporting Statistics | 160 |
| Exporting Group Statistics | 161 |
| Batch Exporting | 162 |
| Using ScreenWorks Peak Pro Software Version 2.0 | 165 |
| Interface Overview | 167 |
| Running Peak Pro 2.0 Analysis | 174 |
| Exporting Data With Peak Pro 2.0 Analysis | 180 |
| Chapter 8: Maintaining the Instrument | 185 |
| Washing Pipettor Tips and Pin Tools | 185 |
| Exchanging Hardware Components | 186 |
| Chapter 9: Troubleshooting | 209 |
| Instrument Status Colors | 209 |
| Troubleshooting Start-Up | 209 |
| General Troubleshooting | 211 |
| Troubleshooting the Pipettor | 214 |
| Troubleshooting the Optics | 218 |
| Troubleshooting the Yellow Plate | 222 |
| Troubleshooting the Tip Washer | 223 |
| Troubleshooting the Cell Reservoir | 224 |
| Troubleshooting Data | 225 |
| Troubleshooting Robotic Integration | 225 |
| Obtaining Support | 226 |

| | |
|---|------------|
| Appendix A: Technical Specifications | 227 |
| Appendix B: Consumables and Accessories | 231 |
| System Accessories | 231 |
| Plates | 234 |
| Assays | 237 |
| Recommended Assay Equipment and Supplies | 240 |
| Appendix C: FLIPR Penta Plate Dimensions | 241 |
| Appendix D: FLIPR Penta Instrument LED and Filter Combinations | 247 |
| Appendix E: Data Processing Algorithms | 249 |
| Hypothetical Experiment | 249 |
| Determining Crosstalk Correction | 250 |
| Determining Spatial Uniformity Correction | 251 |
| Determining Negative Control Correction | 252 |
| Determining Positive Control Scaling | 253 |
| Determining Subtract Bias | 254 |
| Determining Response Over Baseline | 255 |
| Appendix F: Using AquaMax Sterilant | 257 |
| Principle of Use | 257 |
| Contents | 257 |
| Materials Required but Not Provided | 257 |
| Storage | 258 |
| Reagent Preparation | 258 |
| Reagent Use | 258 |
| Warnings, Precautions and Limitations | 258 |
| Appendix G: Robotic Integration | 261 |
| Conventions | 261 |
| Interface Versioning | 261 |
| Instrument Overview | 262 |
| FLIPR Penta System Control Architecture | 271 |
| Setting Up Protocols for Remote Control Use | 278 |
| Command Set | 282 |
| Error Handling and Reporting | 302 |
| Remote Interface Revision History | 305 |

| | |
|--|------------|
| Barcode Specifications | 306 |
| Appendix H: Electromagnetic Compatibility | 309 |

Safety Information

The safety information section provides information on the safe use of the instrument. It includes the use of user-attention statements in this guide, a key to understanding the safety labels on the instrument, precautions to follow before operating the instrument, and precautions to follow while operating the instrument.

Read and observe all warnings, cautions, and instructions. Remember, the most important key to safety is to operate the instrument with care.



WARNING! If the instrument is used in a manner not specified by Molecular Devices, the protection provided by the equipment might be impaired.

Warnings, Cautions, Notes, and Tips

All warning symbols in the user guide are framed within a yellow triangle. An exclamation mark is used for most warnings. Other symbols can warn of other types of hazards such as biohazard, electrical, or laser safety warnings as are described in the text of the warning.

When warnings and cautions are displayed in this guide, be careful to follow the specific safety information related to them.

The following user-attention statements can be displayed in the text of Molecular Devices user documentation. Each statement implies a particular amount of observation or recommended procedure as described:



WARNING! A warning indicates a situation or operation that could cause personal injury if precautions are not followed. The warning symbol can vary depending on the warning. The definition of the symbol is included in the text of the warning.



CAUTION! A caution indicates a situation or operation that could cause damage to the instrument or loss of data if correct procedures are not followed.












Note: A note calls attention to significant information.





Tip: A tip provides useful information or a shortcut, but is not essential to the completion of a procedure.

Symbols on Instrument Labels

Each safety label found on the instrument contains an alert symbol that indicates the type of potential safety hazard related to the label. The following table lists the alert symbols that can be found on Molecular Devices instruments.

| Symbol | Indication |
|---|---|
|  | Indicates that the product documentation must be consulted. Indicates a warning for a situation or operation that could cause personal injury if precautions are not followed. There are specific details written next to the warning symbol. |
|  | Indicates a potential ultraviolet light hazard. |
|  | Indicates a potential pinch hazard. |
|  | Indicates a potential heat hazard. |
|  | Indicates power on. |
|  | Indicates power off. |
|  | Indicates the instrument manufacture date. |
|  | Indicates European technology conformity. |
|  | Indicates certification by TUV Rheinland for compliance with US and Canadian product safety standards. |

| Symbol | Indication |
|---|--|
|  | Indicates compliance with TÜV Rheinland product safety for Germany. |
|  | Indicates compliance with Chinese RoHS Pollution Control Requirements. |

Electrical Safety

To prevent electrically related injuries and property damage, inspect all electrical equipment before use and immediately report all electrical deficiencies. Contact Molecular Devices Technical Support to service equipment that requires the removal of covers or panels.

The instrument must be connected to a properly grounded power outlet to protect from the risk of electric shock. The main chassis of the instrument is grounded together with all related electrical components.

Do not remove the fixed covers, as there are no user-serviceable parts inside. All electrical work must be referred to Molecular Devices approved service personnel.

In the event of a liquid spillage into the main cavity of the instrument, disconnect the mains power supply before trying to clean up.

If the external covers on the instrument are removed, the power supply does not automatically stop.



WARNING! HIGH VOLTAGE Power off the instrument and disconnect the power cord before you do maintenance procedures that require removal of a panel or cover or disassembly of an interior instrument component.

Do not try to use the instrument until all covers are replaced.

To provide access for disconnecting power from the instrument, maintain a 66 cm (26 in.) minimum clearance area on the right side of the instrument.

To protect against fire hazard, replace the fuses only with the same type and rating as the original factory-installed fuses.

Chemical and Biological Safety

Normal operation of the instrument can involve the use of materials that are toxic, flammable, or otherwise biologically harmful. When using such materials, observe the following precautions:

- Handle infectious samples based on good laboratory procedures and methods to prevent the spread of disease.

- Observe all cautionary information printed on the original containers of solutions before their use.
- Dispose of all waste solutions based on the waste disposal procedures of your facility.
- Operate the instrument in accordance with the instructions outlined in this guide, and take all the required precautions when using pathological, toxic, or radioactive materials.
- Splashing of liquids can occur. When working with potentially hazardous liquids, take applicable safety precautions, such as wearing safety glasses and protective clothing.
- Observe the applicable cautionary procedures as defined by your safety officer when using hazardous materials.
- Observe the applicable cautionary procedures as defined by your safety officer when using flammable solvents in or near a powered-up instrument.
- Observe the applicable cautionary procedures as defined by your safety officer when using toxic, pathological, or radioactive materials.



WARNING! Never use the instrument in an environment where potentially damaging liquids or gases are present.

Moving Parts Safety

To prevent injury due to moving parts, observe the following:

- Never try to exchange labware, reagents, or tools while the instrument is operating.
- Never try to physically restrict the moving components of the instrument.
- Keep the interior of the instrument clear to prevent obstruction of the movement.

The motors use high-powered magnets. The linear drive units and encoders are delicate, so be very careful with them. To prevent serious damage to the instrument or its auxiliary parts, follow the preparation instructions in this guide before every process.

The instrument is equipped with an automatically locking door that locks whenever you run a process. The door prevents UV light from passing through during operation.

As a safety measure, if the door is open, an electromagnetic switch prevents the instrument from running. Never tamper with this switch, as it serves two purposes:

- It prevents the motors from running to reduce the potential of physical damage.
- It disables the UV light to prevent the risk of damage from UV radiation.

Motors and their related drives and cabling are sources of electromagnetic fields. Keep magnetic storage devices or strips, such as hard drives and credit cards, away from the instrument covers.



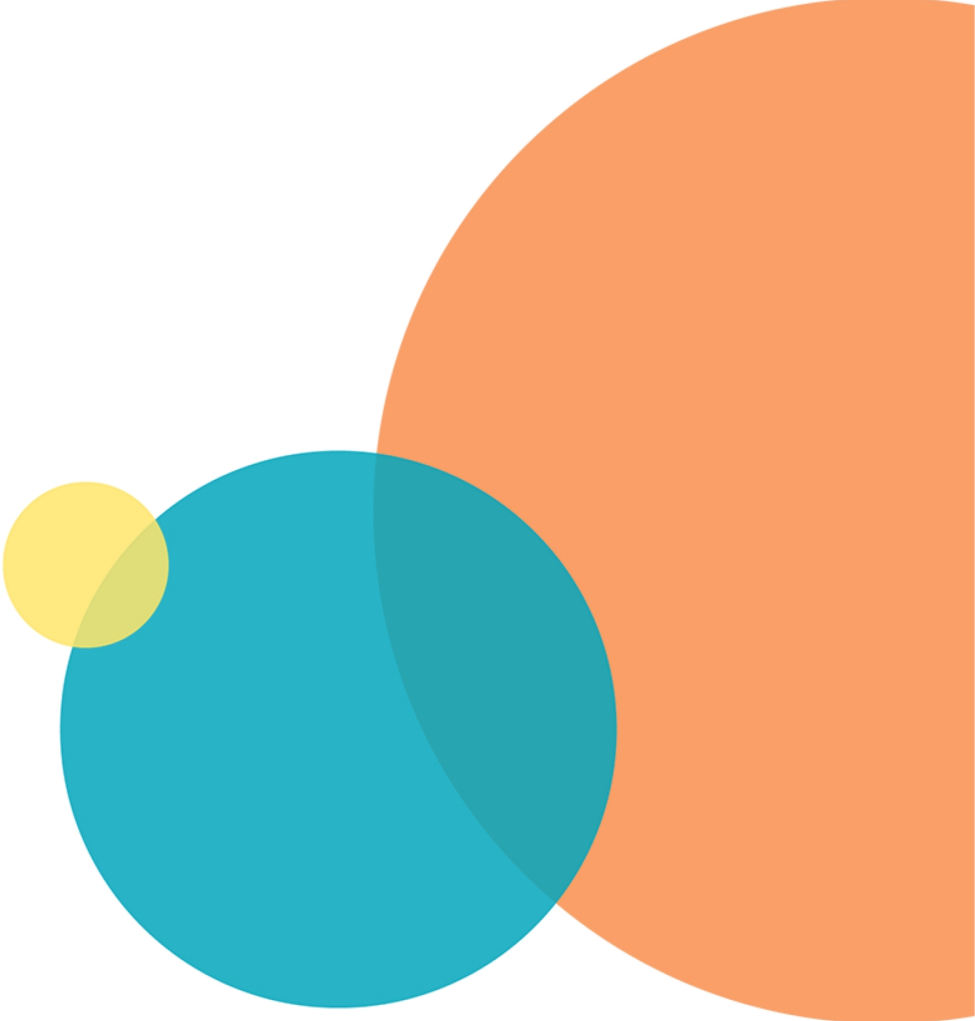
Note: Observe all warnings and cautions listed for all external devices attached to or in use during the operation of the instrument. See the applicable user guide for the operating and safety procedures of that device.

Ultraviolet (UV) Light Safety

The instrument is equipped with an automatically locking door that locks whenever you run a process. The door prevents UV light from passing through during operation.

As a safety measure, if the door is open, an electromagnetic switch prevents the instrument from running. Never tamper with this switch, as it serves two purposes:

- It prevents the motors from running to reduce the potential of physical damage.
- It disables the UV light to prevent the risk of damage from UV radiation.



The FLIPR® Penta High-Throughput Cellular Screening System provides an automated solution for identifying early leads in the drug discovery process and for evaluating drug efficacy and toxicity. With simultaneous pipette and read function, the system supports fast kinetic cellular assays. It can be quickly configured based on library size, detection mode, screening format, assay and target. A fully integrated solution from assay development to lead optimization.

The FLIPR Penta System includes:

- Simultaneous 96-well, 384-well or 1536-well liquid or cell transfer
- Expanded wavelength support
- User-configurable pipettors and optics
- FLIPR® Cycler, internal robotic plate handler option that works in conjunction with external robotic software
- Choice of camera:
 - EMCCD camera for fluorescence applications
 - HS EMCCD high-speed, high-sensitivity camera for fluorescence and luminescence
 - ICCD camera for fluorescence and luminescence
- Cell suspension option
- Compact platform with minimal facilities requirements
- ScreenWorks® Software and optional ScreenWorks® Peak Pro™ Software modules for analysis

This chapter includes the following introductory topics:

- [Workflow on page 14](#)
- [Before You Start on page 14](#)
- [System Requirements on page 14](#)
- [Software Installation on page 16](#)
- [Powering On and Powering Off the System on page 18](#)
- [Checking Instrument Status and Configuration on page 20](#)

Workflow

The FLIPR Penta High-Throughput Cellular Screening System uses the following sequence of processes:

1. Prepare the cells.
2. Power on the system.
3. Run a system check.
4. Load cell dye.
5. Prepare source plates.
6. Set up an assay protocol.
7. Run the experiment.
8. Analyze the data.

Before You Start



CAUTION! Before using the instrument, it is very important that you read and understand all the safety instructions. Then follow the procedures in [Powering On the System on page 18](#).

System Requirements

The following is an overview of electrical, physical and environmental requirements for the FLIPR Penta System.

Electrical

The FLIPR Penta System draws 12 A maximum and requires 100-240 VAC power source at 50-60 Hz. The system ships with a power cord appropriate for the receiving country. Additional shared outlets are required for computer and monitor. A power strip is acceptable for providing the additional outlets for the computer and monitor.

Minimum Lab Space

System dimensions are as follows:

- Without Cell Suspension Module or FLIPR Cyclor: approximately 39 inches wide x 27 inches deep x 70.25 inches high (99 cm wide x 69 cm deep x 1780.44 mm high).
- With Cell Suspension Module and FLIPR Cyclor: approximately 53 inches wide x 27 inches deep x 70.25 inches high (134.6 cm wide x 69 cm deep x 1780.44 mm high).



Note: To open the stage access door completely, you need more than 30 inches (76 cm) additional space in front of the instrument.

The external ScreenWorks Software host computer, is mounted on the right side of the instrument cabinet. The monitor, keyboard, and mouse are installed on an adjustable arm attached to the front right side of the instrument cabinet. Combined, they require an additional minimum amount of space of 40 inches (102 cm) wide x 40 inches (102 cm) deep for maneuverability.

For instruments with the ICCD camera, a chiller with dimensions 11 inches wide by 13 inches deep by 13 inches high (28 cm x 33 cm x 33 cm) is connected to the right side of the instrument with a 3-foot (91.4 cm) long tube. It can be placed anywhere within that 3-foot (91.4 cm) radius as long as there is access to the on/off button on the chiller.

A minimum 32 inch (81 cm) square footprint is required on the right side of the instrument for tip wash bottles.

The instrument cabinet should have servicing access space of 48 inches (122 cm) in front, 24 inches (61 cm) behind , and 10 inches (25.4 cm) on the left.

The FLIPR Penta System is designed with rolling castors so it can be readily moved to make necessary adjustments and perform maintenance. Leveling feet are also installed on the lower instrument chassis. These feet stabilize the instrument when integrated with a robot, but can also be used to establish a uniform instrument deck level in situations where the lab floor is not flat. When running an experiment, make sure the instrument feet are lowered and leveled.



WARNING! The FLIPR Penta System can weigh as much as 860 lbs (390 kg). Follow all necessary safety precautions and use proper moving techniques.

Computer System

The FLIPR Penta System uses two computers, an embedded computer and a host computer.

Embedded Computer

An embedded computer, supplied with the system in the lower chamber of the instrument cabinet, controls all of the FLIPR Penta System basic functions. These functions are initiated through the ScreenWorks Software installed on the host computer, and sent to the embedded computer to execute the function. This configuration allows separate data processing and instrument control to ensure maximum productivity during an experiment.

Host Computer

All normal user interaction with the FLIPR Penta System is done using the ScreenWorks Software that runs on the provided external host computer.

Table 1-1: ScreenWorks Software Computer Requirements

| Item | Description |
|------------------|--|
| CPU | 2.5 GHz, quad processor or faster |
| Operating system | Windows 10, 64-bit Windows 7, 32-bit and 64-bit older systems supported |
| Memory | 16 GB RAM 4 GB RAM minimum |
| Data Connection | 10/100 Ethernet port |
| Storage | 500 GB hard drive |

Software Installation

Your FLIPR Penta System is completely set up in your lab by your Molecular Devices representatives. The following software installation procedures are for when you might need to update the ScreenWorks Software System Control Software on your host computer.

You must have administrative privileges on the computer to install the ScreenWorks Software.

Installing the software includes the following procedures:

1. [Installing ScreenWorks Software on page 16](#)
2. [\(Optional\) Activating the ScreenWorks Peak Pro License on page 17](#)

Installing ScreenWorks Software

As of version 3.1, the installer automatically uninstalls the old software version as long as it is the same major release (3.1 to 3.2). If you are replacing ScreenWorks Software version 3.2 or version 4.x with version 5.0, manually uninstall the old software. See [Uninstalling ScreenWorks Software](#).

Use the following procedure to upgrade your ScreenWorks Software when needed.

To install the software:

1. Double-click the **ScreenWorks_5_0_x.exe** ScreenWorks Software installation file. A **Welcome to the ScreenWorks Setup Wizard** dialog is displayed.
2. Click **Next**.
3. In the **License Agreement** dialog, select **I accept the terms of the license agreement**, and click **Next**.

4. In the **Online/Offline Mode** dialog, designate the default mode in which you want the software to start.
 - In **Online**, ScreenWorks Software automatically looks for a connected instrument when the software is started.
 - In **Offline**, ScreenWorks Software does not automatically look for a connected instrument.

See [Online Mode vs. Offline Mode on page 47](#) for details.
5. Click **Next**.
6. In the **Destination Folder** dialog, the **Install ScreenWorks 5.0 to** field displays the default installation directory. To change the installation directory, click **Change**, navigate to the desired directory, then click **OK**.
7. Click **Next**.
8. In the **Select Program Folder** dialog, leave the displayed default **Program Folder** settings. Select **Anyone who uses this computer** to make ScreenWorks Software available to all users on the FLIPR Penta System host computer, then click **Next**.
9. In the **Configuring the ScreenWorks** installation dialog, if you want to make any changes, click **Back** to go to the previous screen, otherwise click **Next** to start the installation.
10. When the installation is complete, the **Completing the installation process** dialog appears. Click **Finish** to exit the wizard.

(Optional) Activating the ScreenWorks Peak Pro License

The ScreenWorks® Peak Pro™ Software functionality is license-protected. The software license activation enables the Peak Pro analysis functionality any time after the trial period expires. The following Peak Pro analysis licenses are available for purchase:

- ScreenWorks Peak Pro Software version 1.0—adds additional Kinetic Reduction Configuration types
- ScreenWorks Peak Pro Software version 2.0—adds a new Peak Pro 2.0 Analysis dialog and support for high-speed and high-sensitivity data



Note: There is a one time 14-day Peak Pro trial usage option in the Help menu. When the trial period expires, the Peak Pro 1.0 analysis functionality disappears, the Peak Pro 2.0 analysis functionality disables, and the Help menu trial option is inactive. The rest of the ScreenWorks Software version 5.0 remains functional.

To activate the ScreenWorks Peak Pro Software software license:

1. Start the ScreenWorks Software.
2. Click the **Help** tab.
3. Click **Software License**.
4. If you have not started the available 14-day Peak Pro trial, the option to start the 14-day Peak Pro trial, or enter the product key appears. Click **Yes**.

After the trial expires, only the option to provide a software license product key appears.

5. If you have internet connectivity, type the provided **Product Key** in the field and click **Activate Online**, and then follow the on-screen instructions.
6. If you do not have Internet connectivity, click **Activate Offline** and follow the on-screen instructions. Activate Offline requires the following:
 - Your product key
 - A separate computer with Internet connectivity
 - A USB drive for transferring files between the computers
7. Restart ScreenWorks Software to finish the license activation.

Uninstalling ScreenWorks Software

The following is the recommended method for removing ScreenWorks Software from a Windows-based computer because it also removes related information from the Windows Registry.

To uninstall ScreenWorks Software:

1. Open **Control Panel**.
2. Depending on your Windows operating system:
 - Windows 7—Click **Add or Remove Programs**
 - Windows 10—Click **Programs and Features**
3. In the list that displays, select **ScreenWorks**.
4. Click **Remove** or **Uninstall**.
5. Follow the on-screen instructions to finish uninstalling the software.

Powering On and Powering Off the System

The following procedures are for powering on and powering off the FLIPR Penta System, which includes connecting the ScreenWorks Software to the instrument.

Powering On the System

Before continuing, do as instructed in [Before You Start](#).

To power on the FLIPR Penta System:

1. Power on the computer and monitor.
2. Simultaneously press the **CTRL+ALT+DELETE** keys to launch the Windows operating system.
3. At the prompt type your password.



Note: The first time you log in after your system installation, the default password is *fliprtetra*.

Wait for the operating system to finish starting-up before continuing.

4. Power on the external chiller with the switch located on the left side of the chiller.

- Power on the FLIPR Penta System power switch located on the right side of the instrument.



CAUTION! The system goes through an initialization cycle to register all of the instrument components. This cycle is not complete until the green **Assay Finished (Unlock)** light on the upper door is the only light illuminated on the instrument status panel.

- Double-click on the  desktop icon to start the ScreenWorks Software.




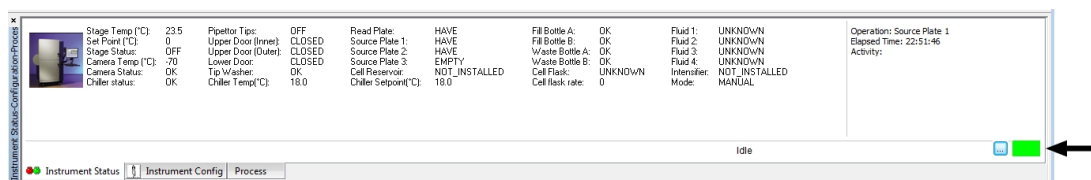
Tip: Do not repeatedly double-click the software icon. Starting the software can take several seconds.




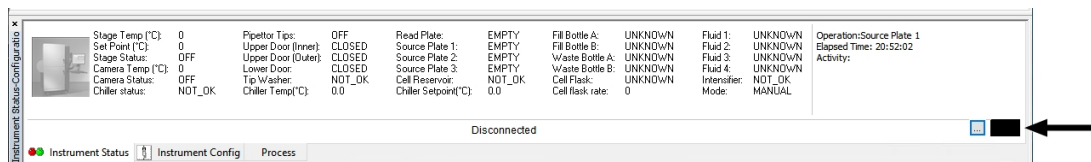
Note: The system is ready for use when the camera temperature is at operating temperature. The camera temperature and status display in the ScreenWorks Software Instrument Status panel. The camera operating temperature depends on the camera in your instrument. See [CCD Camera Options on page 35](#).

Connecting the Software to the Instrument

After powering on the instrument, when you start the ScreenWorks Software, it should automatically connect to the instrument. In the software, in the Instrument Status panel, in the lower right corner of the screen, a green connection status icon indicates the connection and in the toolbar, the  button displays.



If you see a black connection status icon, the software is disconnected from the instrument, and in the toolbar, the  button displays.



To connect to the instrument, click .

Powering Off the System and the Software

To power off the FLIPR Penta System:

1. At the end of an experiment run, wait for the **Assay Finished (Unlocked)** light on the upper door, Instrument Status Panel turns green, indicating the experiment is finished.
2. If the last protocol run did not remove the tips from the pipettor head, Molecular Devices recommends making sure that you remove them using the manual command. Failure to remove tips can result in an error on next start.
3. Exit ScreenWorks Software by selecting **File > Exit**.
4. Power off the computer and monitor.
5. Power off the FLIPR Penta instrument.
6. Power off the chiller.

Checking Instrument Status and Configuration

You can check instrument status on the front of the instrument and from within the ScreenWorks Software. You can only check the instrument configuration status from within the ScreenWorks Software. For more details see [System Status Panel](#) and [Software Status Tabs](#) on page 21.

System Status Panel

The system status panel, located next to the upper door handle, indicates if the door is locked or unlocked, depending on the instrument activity. It also includes an emergency Interrupt button to stop any running processes.

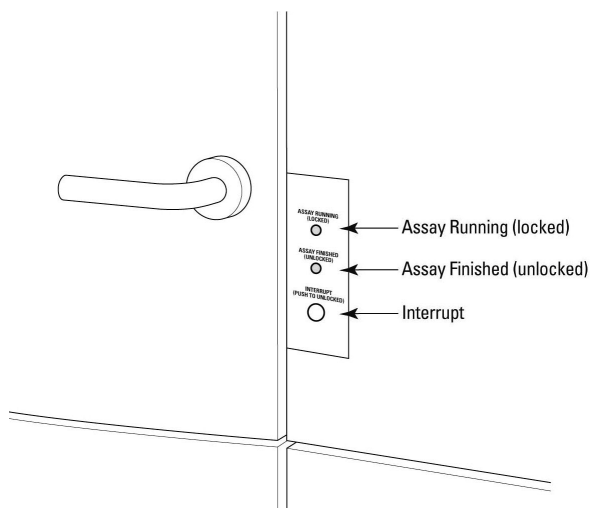


Figure 1-1: The System Status Panel

The panel has two lights and the Interrupt button. From the top of the panel these are:

- **Assay Running (Locked)**—Yellow light
The FLIPR Penta System is performing a task. The upper and lower doors are locked until the task finishes or is stopped using the Interrupt button.

- **Assay Finished (Unlocked)**—Green light
No tasks are being run and it is safe to open the upper and lower instrument doors.
- **Interrupt**—Flashing green light
A system override to stop all tasks, so you can access the instrument. When pressed, the light flashes until the system has reached a safe state to open the doors.



CAUTION! The Interrupt button immediately ends the experiment and should only be used in emergencies. Before you can use the instrument normally again, the system might need to be reinitialized from within the software by selecting **Instrument > Reset** from the menu.

Software Status Tabs

Within the ScreenWorks Software, the **Instrument Status-Configuration-Process** tab panel, located on the bottom of the main screen, is open automatically and reports the status of and settings for the FLIPR Penta System hardware; and it includes the processes used to create protocols. For more details, see [Instrument Status-Configuration-Process Panel on page 66](#).

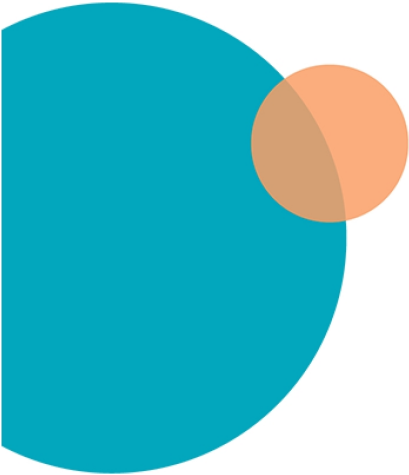
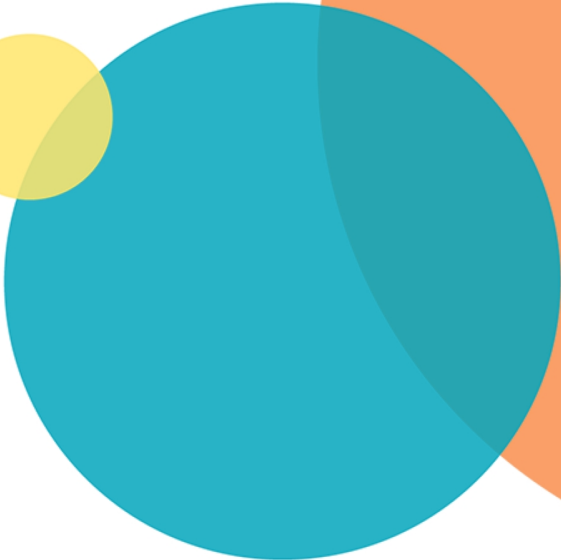
The screenshot displays the ScreenWorks software interface. The main window is titled 'ScreenWorks 5.0.0.146 - [Untitled1]'. The interface is divided into several sections:

- Menu Bar:** File, View, Instrument, Tools, Window, Help.
- Toolbar:** Contains various icons for file operations and instrument control.
- Main Display Area:**
 - Detail Graph:** A line graph showing 'Relative Fluorescence Units' on the y-axis (0 to 1) and 'Time (Seconds)' on the x-axis (0 to 1.0). The graph is currently empty.
 - Group Statistics Graph:** A table with columns for 'Group Name' and 'Well ID', and a 'Kinetic Reduction Value' column. It is currently empty.
- Bottom Status Bar:**
 - Instrument Status:** Active tab, showing a table of system parameters and their states.
 - Instrument Config:** Tab for configuration settings.
 - Process:** Tab for process management.

The 'Instrument Status' tab displays the following table:

| Parameter | Value/State | Parameter | Value/State | Parameter | Value/State | Parameter | Value/State |
|------------------|-------------|-------------------|-------------|--------------------|---------------|----------------|-------------|
| Stage Temp (°C) | 23.0 | Pipettor Top | ON | Read Plate | HAVE | Fill Bottle A | OK |
| Seal Temp (°C) | 0° | Upper Door (Inn) | CLOSED | Source Plate 1 | HAVE | Fill Bottle B | OK |
| Stage Temp (°C) | 0° | Upper Door (Out) | CLOSED | Source Plate 2 | HAVE | Wash Bottle A | OK |
| Camera Temp (°C) | 20° | Lower Door | CLOSED | Source Plate 3 | EMPTY | Wash Bottle B | OK |
| Camera Status | OK | 12° Heater | OK | Cell Reference | NOT_INSTALLED | Cell Plate | UNKNOWN |
| Chiller Status | OK | Chiller Temp (°C) | 17.9 | Cell Setpoint (°C) | 18.0 | Cell Wash Rate | 0 |

The bottom status bar shows 'Ready' on the left and 'NUM' on the right. A green light indicator is visible on the right side of the status bar.



Chapter 2: System Hardware Overview

The FLIPR Penta System consists of a cabinet 39 inches (965 mm) wide x 27 inches (686 mm) deep x 70 inches (1780 mm) high, with two compartments, accessed by manually opening front doors, top for wet components and bottom for dry components. Outside the cabinet are the wash system fluid bottles, optional Cell Suspension module, chiller, computer, monitor, keyboard, and mouse.

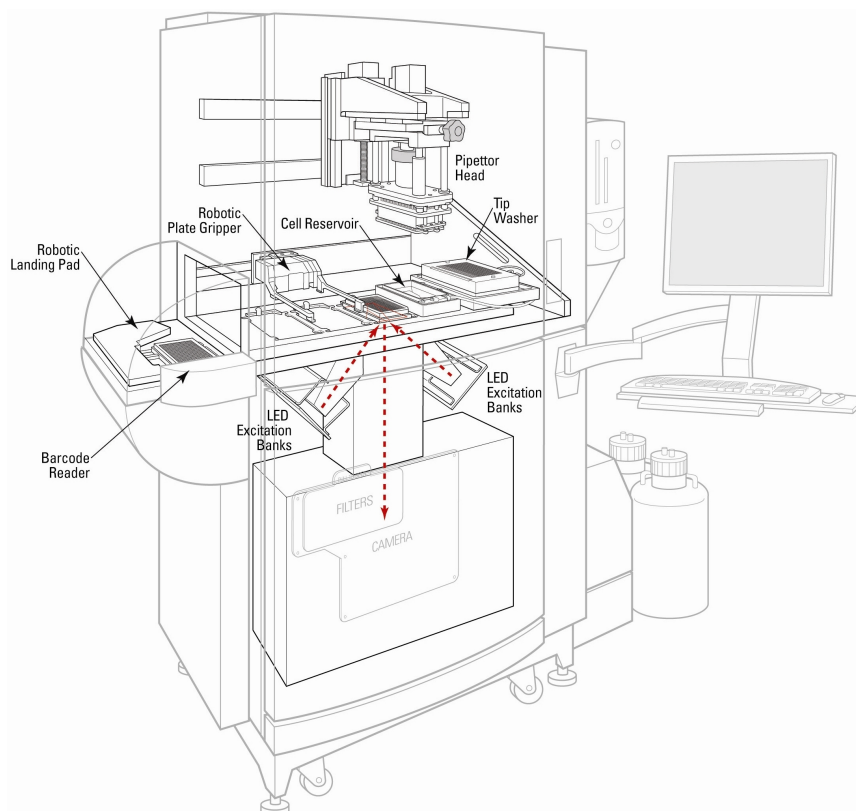


Figure 2-1: FLIPR Penta System With Optional Modules

For system component details, see the following topics:

- [Five-Position Stage on page 24](#)
- [Plate Support on page 25](#)
- [Plate-Handling System on page 26](#)
- [Liquid-Handling System on page 29](#)
- [Optical System on page 35](#)
- [Chiller on page 45](#)
- [Observation Panel on page 45](#)

Five-Position Stage

You run your experiment from the five-position stage in the upper compartment of the FLIPR Penta System. You can manually load plates through the upper compartment door before an experiment or robotically during the experiment using the optional FLIPR® Cycler internal plate handler.

The following are the five plate positions of the stage from left to right:

- Position 1—Tips and Source Plate 1
- Position 2—Source Plate 2
- Position 3—Read Plate
- Position 4—Cell Reservoir and Source Plate 3
- Position 5—Tip Washer

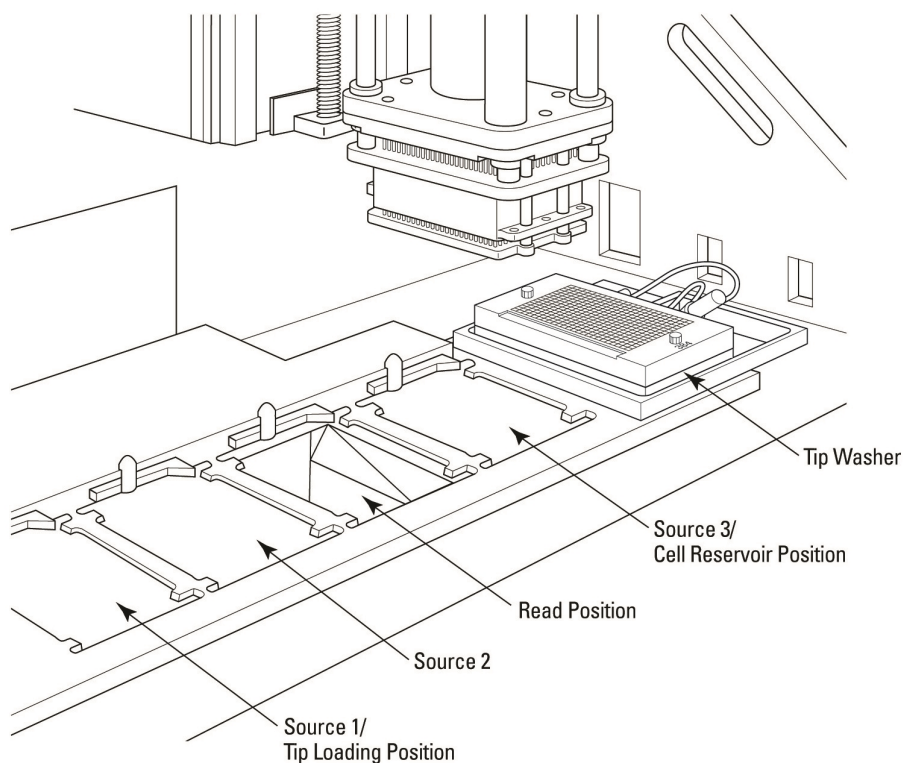


Figure 2-2: The Five-Position Stage

Positions 1, 2, and 4 support standard, low-volume, deep-well and reservoir source addition plates.

Position 1 is for tip loading and unloading, but this position can also be used as a source plate position after tips load and the tip box is removed.

Position 3 opens to the optics chamber below for excitation of fluorophores in read plate wells, the read position, and emission reading.

Position 4 is used as a source plate position when the Cell Reservoir, included with the optional Cell Suspension option, is not in use. A single Cell Reservoir is compatible with all of the FLIPR Penta System pipettor heads.

Positions 1 through 4 have a mechanical plate sensor to identify the presence of plates, tips or reservoirs.

Position 5 is used as a dedicated tip wash reservoir and should be configured to match respective pipettor heads, such as 96, 384, or 1536. Appropriate tip wash reservoirs are included in the purchase of a pipettor head. For more details, see [Washing Pipettor Tips and Pin Tools on page 185](#).

Robotic integration enables the optional FLIPR Cycler internal plate handler to automatically exchange in an experiment up to 12 source plates and tip racks and one read plate. See [Automating Plate Loading on page 26](#).

The plate indexers in Positions 1 through 4 registers plates and tip racks with well A1 in the lower left corner. The indexers also serve as mechanical sensors to detect the presence of plates or tips. If plates or tips are not sensed during a Manual Mode experiment, but requested by the software, the instrument stops and ends the experiment. During Remote Mode, the system notifies the SynchroMax™ ET plate handler robot or third-party plate handler robot that no plate or tip container is available and stops the instrument until it detects plates or tips. It is then the responsibility of the SynchroMax ET plate handler or third-party plate handler to deliver plates or tips to the system. See [SynchroMax ET External Plate Handler on page 28](#).



CAUTION! Sensors can only detect the presence of a plate or a tip container. Sensors cannot identify the type of plates or tips, or if the lids are on the plates or not. It is your responsibility to ensure that the correct plates and tips are loaded into position without lids.



Note: Molecular Devices no longer sells the SynchroMax ET plate handler robot. The FLIPR Penta System becomes a slave to whichever attached robot system is configured.

Regulating Stage Temperature

Plate stage positions 1, 2 and 4 for source plates, have temperature control. Stage temperature settings range from ambient 25°C (77°F) to 40°C (104°F). It can take approximately 15 minutes to reach equilibrium temperature.

Configure temperature regulation with the **Temperature Control** ON/OFF toggle command in the **Instrument > Manual Operation** menu or corresponding button.

Plate Support

FLIPR Penta System supports the use of 96-well, 384-well and 1536-well plates that conform to the ANSI standards submitted by the Society for Biomolecular Sciences.



Note: Only low-profile 1536-well plates are supported.

Black walled, clear-bottomed read plates provide an optimal imaging environment for fluorescence assays. These plates prevent signal diffraction while allowing excitation and signal access. Black walled, clear-bottomed plates or white walled, clear-bottomed plates can be used for luminescence assays.

For 96-well read plates, an optional slit-shaped mask can be used to minimize saturation and edge effects associated with these plates. Place the mask over the read position. See [Consumables and Accessories on page 231](#) for types of masks available.



Note: No mask is required for 1536-well and 384-well plates.

The 96- and 384-pipettor heads can be used with source or read plates with equal or one order higher well number. This is because the FLIPR Penta System can aspirate or dispense into quadrants of a plate. The following combinations are possible:

- The 96-pipettor head can be used with 96-well and 384-well plates.
- The 384-pipettor and pin tool head can be used with 384-well and 1536-well plates.
- The 1536-pipettor and pin tool head can be used only with 1536-well plates.

Deep-well plates or reservoirs can be substituted for standard well plates.

When compound is aspirated or delivered to a plate with a greater number of wells than the pipettor head, the quadrant number (1 to 4) must be entered in the protocol configuration in ScreenWorks Software for each dispense.

Many of the compatible source and read plates are available through Molecular Devices. See [Consumables and Accessories on page 231](#).

For more compatible plate details, see [FLIPR Penta Plate Dimensions on page 241](#).

Plate-Handling System

The plate handling system includes the following methods for placing plates on the five-position stage to run your experiments:

- [Loading Plates Manually on page 26](#)
- [Automating Plate Loading on page 26](#)

Loading Plates Manually

In Manual Mode, before running an experiment, position all of the assay components in the five-position stage by hand in the upper compartment. No additional plate or tip changes can be made after the experiment starts.



Note: Keep the top compartment door closed during normal system operation. All system functions stop when the door is open.

Automating Plate Loading



Note: Molecular Devices no longer offers the SynchroMax ET external plate handler robot. The following information is for reference only.

Support for automated plate delivery with the FLIPR Penta System is facilitated using optional robot systems.

To increase the number of plates you can use in an experiment, and minimize personnel requirements, use the optional FLIPR Cyclor internal plate handler coupled with the optional external SynchroMax ET plate loader robot or a third-party plate loader robot, for example, stacker system or robotic arm. One read and up to 12 source plates and tip racks can be shuttled in and out of the FLIPR Penta System in one experiment using automated delivery.

When using automated delivery the SynchroMax ET plate loader or third-party plate loader delivers plates to, and picks them up from, the landing pad on the outside left of the instrument, from where the FLIPR Cyclor plate handler shuttles them in and out of the plate stage compartment. The shuttle door over the landing pad opens and closes to maintain a light-tight environment within the compartment.

During automated operation the SynchroMax ET plate loader or third-party plate loader controls the FLIPR Penta System by sending instructions to load protocols, run experiments, and retrieve plates from the landing pad. These commands are executed immediately upon receipt by the instrument. Persisting instrument settings cannot be made from the remote controlling program—these must be configured in ScreenWorks Software before control is passed to the plate-delivery system.

To start Remote Mode:

- To pass control to the plate-delivery software, in ScreenWorks Software, select **Instrument > Set Remote Mode**.

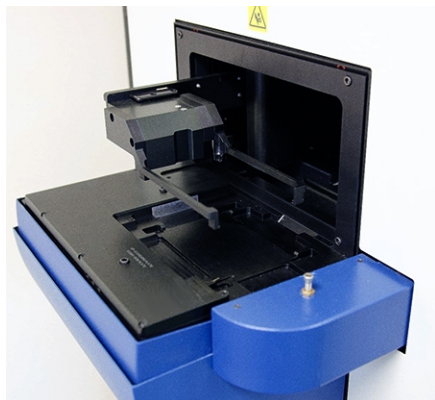
To stop Remote Mode:

- The FLIPR Penta System stays in remote control until you select **Instrument > Set Manual Mode**.

The third-party plate loader software communicates with the FLIPR Penta System computer through the serial communication port using TCP/IP. SynchroMax ET software is installed with ScreenWorks Software on the FLIPR Penta System computer so it is able to communicate directly with the instrument. See [Robotic Integration on page 261](#) for remote control syntax.

FLIPR Cyclor Internal Plate Handler

The optional FLIPR Cyclor plate handler system is a plate gripper that runs along the back wall of the upper read compartment, giving it access to Positions 1 through 4 in the five-position stage and the landing pad on the outer left-side of the cabinet.



The FLIPR Cycler plate handler shifts source plates and tip containers between these locations when under the control of the external SynchroMax ET plate holder robot or a third-party plate handler.

The FLIPR Cycler plate handler carries standard, low volume and deep well 96-well, 384-well and 1536-well plates that conform to proposed ANSI standards submitted by the Society for Biomolecular Sciences. In addition, the FLIPR Cycler plate handler can move Molecular Devices qualified tips.



Note: Reservoirs can be used during robotic integration, however the FLIPR Cycler plate handler is not able to move reservoir. All reservoirs must be loaded manually before running an experiment, including the Cell Reservoir.



CAUTION! While the system is compatible with plates that conform to proposed ANSI standards submitted by the Society for Biomolecular Sciences, some plates may not be handled as reliably by the FLIPR Cycler plate handler because of their low plate weight.

During robotic integration, Molecular Devices recommends evaluating plates and tips handling by the FLIPR Cycler plate handler for handling reliability before starting a screen. The FLIPR Penta System is only validated for use with the Molecular Devices tips. We are not able to guarantee performance or troubleshoot any instrument issues with any other tips.

The upper and lower door of the FLIPR Penta System must remain closed for the duration of the experiment. Plates are transported in and out of the instrument only through the FLIPR Cycler plate handler robotic landing pad door.



WARNING! Do not place your fingers in the FLIPR Cycler shuttle door as this can cause injury.

SynchroMax ET External Plate Handler



Note: The SynchroMax ET plate handler robot is no longer an available option for new purchases. Other robotic systems are still supported.

The SynchroMax ET robot is a six-stack plate handler that use to be available as an optional purchase with the FLIPR Penta System. It delivers plates to and from the landing pad, integrating with the FLIPR Cyclor internal plate handler, which ferries the plates to and from their appropriate locations in the five-position stage. The configuration interface of the SynchroMax ET software is opened directly from within ScreenWorks Software.

Liquid-Handling System

The liquid-handling system includes all the following components involved in moving liquids to and from the five-position stage:

- [Standard Pipettor Head](#)
- [Pin Tool Head on page 30](#)
- [Pipettor Tip and Pin Tool Washer on page 31](#)
- [Cell Suspension on page 32](#)

Standard Pipettor Head

Standard pipettor heads are available in 1536-tip, 384-tip and 96-tip formats.

The 384-pipettor head and 96-pipettor head both use disposable plastic tips. In contrast, the 1536-pipettor head uses a stainless steel tip block with a disposable 1536-tip gasket.

You can wash the plastic tips, replace them between each compound addition, or at the end of an experiment. The 1536-tip block is washed at specified times.

Pipettor operations are controlled from within ScreenWorks Software protocols, or some operations, for example, loading tips can be done individually directly through commands in the Instrument menu. Connectors on the back of the pipettor head identify the head format as 1536-tips, 384-tips, or 96-tips, so ScreenWorks Software only offers valid plate formats and pipetting parameters for protocol setup.

The standard pipettor head uses air displacement to control aspiration and dispense speed and volume. Configure the volume of compound to be transferred in the software. You can also configure drawing compound from multiple source plates to dispense into one destination plate or to aspirate from one plate and dispense to multiple well plates or quadrants.

The instrument also has 384-pin and 1536-pin pipettor heads available with pin tool and specific aspirating, dispensing, and wash protocols. These protocols activate only when the appropriate head is installed.

Table 2-1: FLIPR Penta Fluidics Specifications

| Wells | Pipettor Head Range | Pipetting Increments | Precision |
|-------|---------------------|----------------------|---------------|
| 96 | 5–200 µL | 1 µL | 3% CV @ 75 µL |
| 384 | 1–25 µL | 0.5 µL | 4% CV @ 25 µL |
| 1536 | 0.5–3 µL | 0.1 µL | 6% CV @ 3 µL |

The 96-pipettor head and 384-pipettor head displace air in the disposable pipette tips.

In the 1536-pipettor head, a plunger for each of the 1536 tips presses against an elastic gasket seated on the tip block. When the plungers move down, they create an initial seal between the gasket and tip-block. After creating the seal, further plunger movement causes air displacement in the tip-block.

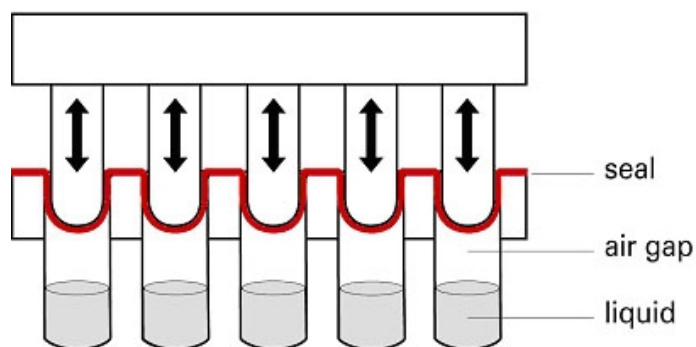
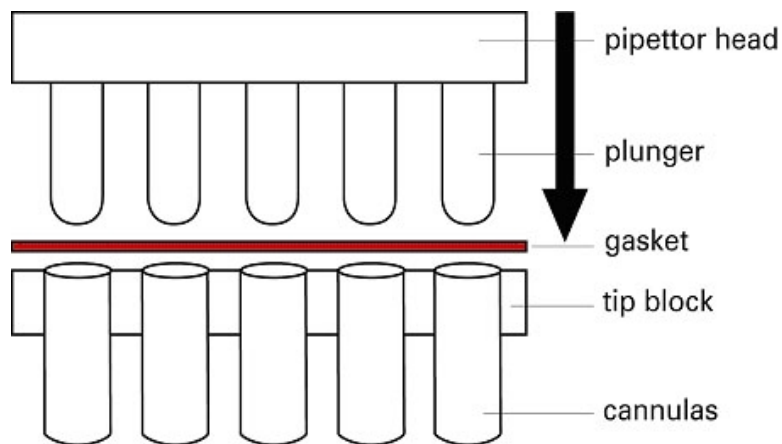


Figure 2-3: Seal creation in the 1536-pipettor head

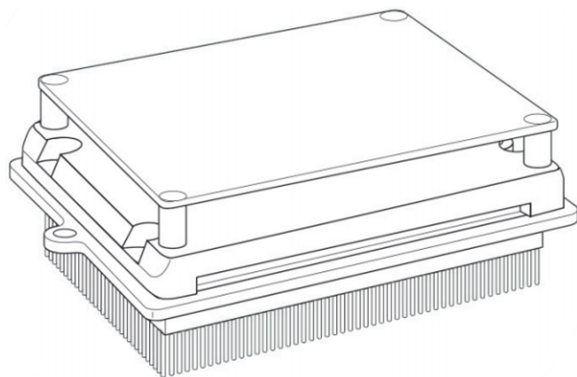


CAUTION! Performance depends on tip-block and gasket seating and can be compromised if the seal breaks.

Use only Molecular Devices recommended tip-blocks and gaskets to ensure the highest accuracy and to reduce the possibility of damaging the pipettor and tips. For recommendations, see [Consumables and Accessories on page 231](#).

Pin Tool Head

Pin tools are blocks of solid or slotted pins, where the pins replace the hollow tips used with a standard pipettor. The pins use capillary action to pick up and transfer liquid from one plate to another. Their ability to accurately and reliably transfer compounds in nanoliter volumes allows you to supply test compounds in 100% DMSO solution, removing the need to prepare intermediate dilution plates.



The volume that each pin picks up is determined by the size of the pin (and, if slotted, the size of the slot) and the withdrawal speed of the pin from the liquid—a faster removal speed leaves more liquid on the pin.

Pins for the 384 pin tool are supplied in four sizes, giving a total range (across all these sizes) from 84 nL to 656 nL. The 1536 pin tool has seven different pin sizes, giving a total range from 19 nL to 117 nL.

Each pin size has a specified volume range that it carries:

- The lowest reported volume is for a tip removal speed of 7.8 mm/s.
- The highest reported volume is for a tip removal speed of 57.0 mm/s.

You specify the precise pick up volumes and tip removal speeds during your assay development.

The FLIPR Penta System can be fitted with 384- or 1536-pin tool heads. Pin tools themselves, in the appropriate 384 or 1536 format, can be easily and rapidly replaced to change the pin size. See [Exchanging Pin Tools on page 195](#).

All the pin tools used with the FLIPR Penta System are available with a hydrophobic and lipophobic coating to prevent or reduce the nonspecific binding of proteins and lipids to the pins.

In order to ensure uniform compound pick-up across the entire pin tool, pins can be configured to ‘float’ in source plate wells. Individual pins are not rigidly attached to the pin block, having a small amount of vertical movement up into the block. When set to float, the pin head moves down very low so that all pins sit on the bottom of the well and push up a little into the block. This ensures that all pins are equally immersed in their wells, for example, sitting on the bottom. This outcome could not be guaranteed if the pins were rigidly fixed to the block, given that plate bottoms are often not completely flat.

Pipettor Tip and Pin Tool Washer

The washer consists of a reservoir top of the selected pipettor format, mounted over a wash basin. The wash basin is connected to two solvent-supply carboys and two waste carboys located on the floor beneath the computer monitor. A basin beneath the tip washer base drains to a waste carboy, to safely remove any solvent that overflows from the reservoir.

Wash solution fills the reservoir for a calibrated amount of time. Solvent is then drained from the reservoir after each wash cycle. Up to five wash cycles can be configured within a single Wash Tips process in your experiment protocol. See [Wash Cell Reservoir Process on page 124](#), and [Creating New Protocols on page 124](#).

For pipettor tips, set volume of solvent is drawn up, optionally held for a time, and then expelled, up to 20 times.

For pins, vertical motion of the tip block is used to agitate the wash solvent around the pins.

The option is available to wash tips or pins in up to two solutions before reusing the tip washer. When additional wash solutions are required, pipettor tips or pins can be washed in a boat or reservoir, located in one of the source plate positions and using a protocol Mix Fluid process configuration. See [Mix Fluid Process on page 113](#).



WARNING! High volumes of volatile, flammable solvents in the reading chamber can cause explosive conditions. Use of 100% isopropanol, etc., in the tip washer is particularly discouraged without additional ventilation. Consult your facilities expert to determine the appropriate ventilation to avoid explosive conditions.

Pin tools are supplied with blotting stations that can be loaded into one of the plate positions on the stage, such as Position 4 (Source 3).

Blot pin steps can be configured in the protocol to remove fluid from the pins, for example, following pin washing.



Note: A waste bottle sensor override (P/N 0700-0827) is available in the FLIPR Penta System accessory kit to bypass the waste sensor and dispose waste in containers other than the dedicated waste carboys.

To exchange the wash reservoir top, see [Exchanging Pipettor Tips and Pin Tool Wash Reservoir Top on page 191](#).

Cell Suspension

The Cell Suspension module is optional. The first installation of the cell reservoir is done by a Molecular Devices certified field service engineer.

The Cell Suspension module consists of two components:

- The Cell Suspension module ([Figure 2-4](#)) located externally on the right-side of the instrument.
- The Cell Reservoir ([Figure 2-6](#)) installs in Position 4 (Source 3) in the 5-position stage. See [Five-Position Stage on page 24](#).

The Cell Suspension module ([Figure 2-4](#)) consists of a shelf with a magnetic motor mounted underneath it, a cell flask with a magnetic stirrer, up to four fluid bottles for automated cleaning, and a removable cover for keeping cells in a dark environment.

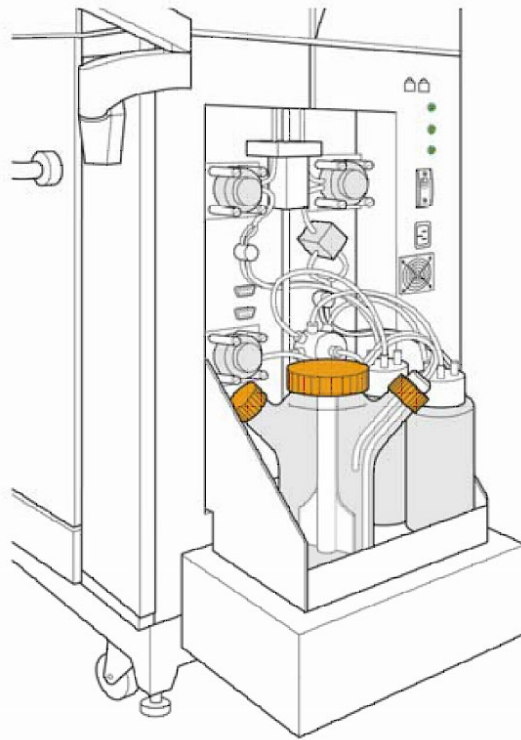


Figure 2-4: External Cell Suspension Module

The spinner flask contains a stirrer, which is driven by a magnetically coupled motor mounted in the lower part of the Cell Suspension module. The Cell Suspension stir speed can be set in a protocol or with manual instrument controls. Stir speed of 5 equals approximately 1 revolution per second. From the uniform suspension in the spinner flask cells are pumped into the Cell Reservoir, where the pipettor head in 96, 384, or 1536 format removes the appropriate amount and cells are automatically pumped to a specified destination. A protocol in ScreenWorks Software controls the stir speed and source/destination of cell suspension activity.

The cell valve selects the source for filling, or the destination for draining. There are 8 valve positions: flask fill for filling the reservoir, flask drain for draining the reservoir without causing air bubbles, Waste Bottle A, Waste Bottle B, and Fluid 1 to Fluid 4.

Fluid 1 to Fluid 4 are user specifiable, and can be cleaning solutions, water or buffer. You can choose to pump cells back into the cell flask or to any other fluid bottles. Bottles for Fluid 1 and Fluid 2 are automatically included with the Cell Suspension option.

Figure 2-5 demonstrates how the system is connected and shows all the possible combinations for protocol development.

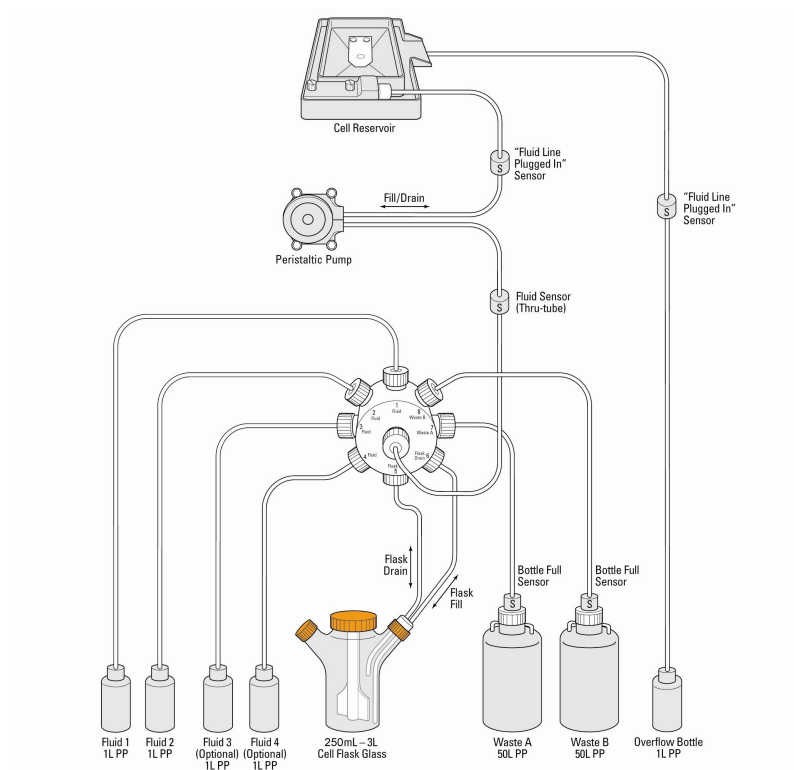


Figure 2-5: Cell Suspension Module Connections

The Cell Reservoir is a special plate type that has one fluid line used for both input and output, and an electronic plate ID that is part of the electrical/fluidic connector, which identifies the reservoir. In the case of an overflow, an overflow trough catches excess fluid and pipes it to the tip washer overflow trough, which directs it to the overflow bottle.

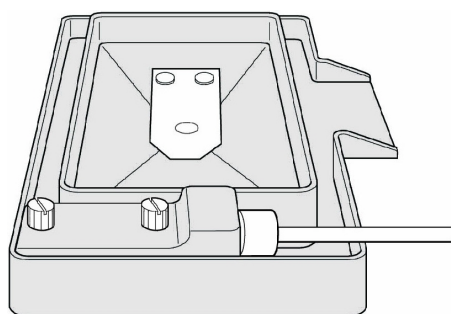


Figure 2-6: Cell Reservoir

When in use, the Cell Reservoir is filled from any of the bottles in the external Cell Suspension module by a pump with adjustable speed and direction.

Wash the Cell Reservoir by either adding the Wash Reservoir process to the protocol (see [Creating New Protocols on page 124](#)), selecting **Wash Reservoir** in the **Instrument > Manual Operation** menu, or manually removing the reservoir and autoclaving it.

Optical System

The FLIPR Penta System optics are housed in the bottom compartment of the main cabinet. In fluorescence assays, where excitation is required, light from light-emitting diodes (LEDs) is directed at the base of the read plate exposed in position 3 in the 5-position stage above. Light emitted from the plate travels down through emission filters before being captured in the CCD camera.

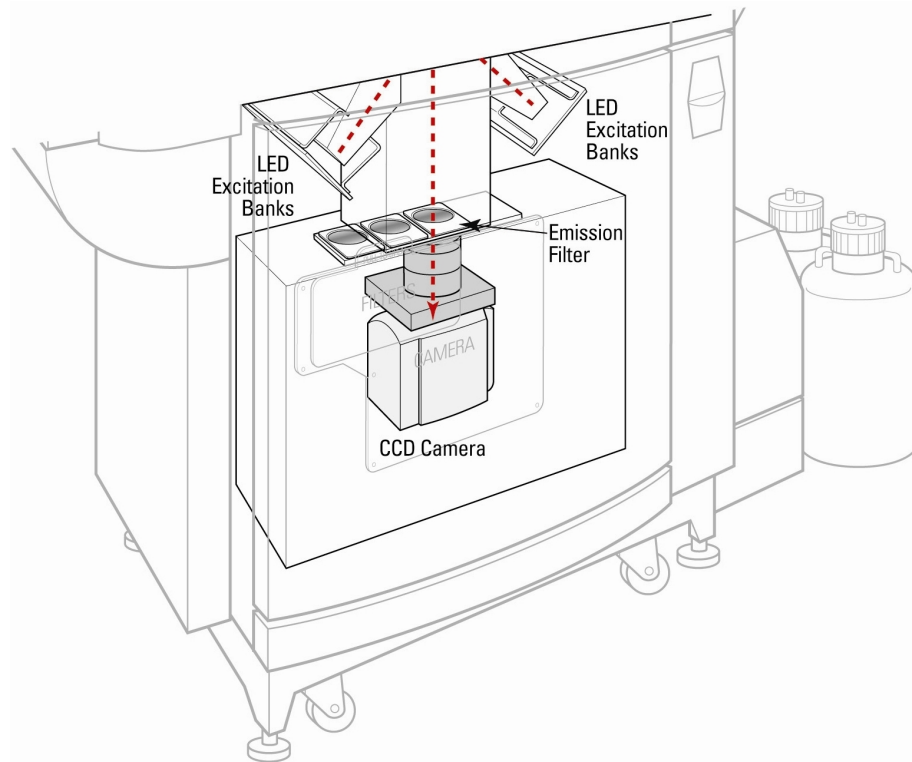


Figure 2-7: System Optics

CCD Camera Options

Three camera options are available for purchase with the FLIPR Penta System:

- EMCCD (Electron Multiplying CCD)—camera for fluorescence only experiments
- HS EMCCD (High-Speed, High-Sensitivity Electron Multiplying CCD)—camera for both high-speed, high-sensitivity, fluorescence, and luminescence assays
- ICCD (Intensified CCD)—camera for both fluorescence and luminescence assays

EMCCD Camera

The camera is located directly beneath the read plate on the five-position stage. The camera is thermoelectrically cooled and requires about five minutes to reach its operating temperature of -70°C (-94°F) $\pm 2^{\circ}\text{C}$.



CAUTION! The instrument does not let you use the camera before it has reached proper operating temperature. Check the camera temperature on the Instrument Status panel before starting an experiment.

The camera uses frame-transfer technology that shifts pixel values from the exposed portion of the chip to the data processing portion. This method allows a high acquisition rate and eliminates the possibility of camera shutter failure.

The camera is an integrating-type detector using temporal integration to build up the signal-to-noise ratio. Depending on the intensity of the emitted light (reliant on dye efficiency and LED power), it may be necessary to vary camera exposures. This prevents the measured fluorescence signal from being dominated by detector noise.

Images are taken of the bottom of the entire plate for a time specified in the ScreenWorks Software protocol; exposure time can be set from 0.05 to 30 seconds. In ratiometric experiments for example, where two or more wavelengths are measured, the number of images captured increases so that an image is taken for each wavelength at the specified rate. From each image a relative light value is calculated for each well. The FLIPR Penta System reports relative fluorescence units (RFUs) in a range from zero to approximately 12,000.

The light intensity detected by each pixel on the CCD chip can be amplified using the Gain setting in ScreenWorks Software (**Settings Process**). This parameter has a range from one to 240. Amplification is exponential, with increments increasing as you go higher up the range. Fluorescence assays typically use a gain of 80 whereas luminescence assays should use 200 as a starting point. Gain optimization should be done during assay development to determine optimal conditions for your screen.

When a luminescence experiment follows a fluorescence experiment we recommend that you cycle the camera temperature to eliminate ghost images that may have been created during the fluorescence assay. Select **Cycle Camera Temperature** from the **Instrument > Manual Operation** menu to choose this option. The camera warms up to room temperature to release ghost images, before cooling back down to approximately -70°C (-94°F).



Note: Luminescence readings are available with the EMCCD camera option, however camera sensitivity is not optimal for this type of experiment.

HS EMCCD Camera

The camera is located directly beneath the read plate on the five-position stage. The camera is thermoelectrically cooled and requires about five minutes to reach its operating temperature range between -70°C (-94°F) $\pm 2^{\circ}\text{C}$.



CAUTION! The instrument does not let you use the camera before it has reached proper operating temperature. Check the camera temperature on the Instrument Status panel before starting an experiment.

The camera uses frame-transfer technology that shifts pixel values from the exposed portion of the chip to the data processing portion. This method allows a high acquisition rate.

The camera is an integrating-type detector using temporal integration to build up the signal-to-noise ratio. Depending on the intensity of the emitted light, which is reliant on dye efficiency and LED power, it might be necessary to use shorter or longer camera exposures. This prevents the measured fluorescence signal from being dominated by detector noise.

Images are taken of the bottom of the entire plate for a time specified in the ScreenWorks Software protocol. There are a variety of exposure settings available depending on the mode in which the camera is run. In ratiometric experiments for example, where two or more wavelengths are measured, the number of images captured increases so that an image is taken for each wavelength at the specified rate. From each image a relative light value is calculated for each well. The FLIPR Penta System reports relative fluorescence units (RFU) in a range from zero to approximately 64,000 RFU.

The light intensity detected by each pixel on the CCD chip can be amplified using the **Gain** setting in ScreenWorks Software (**Settings Process**). This parameter for the Fluorescence Reading Mode has a range from 1.5 to 128. In the Camera Mode > Normal, Fluorescence assays typically use a gain of 8 for bright calcium assays depending on the exposure time and LED Power. This needs to be optimized for each different assay. Luminescence assays should use a gain of 1.5 to 500 as a starting point. Run gain optimization tests during assay development to determine optimal conditions for your screen.



Tip: When using the Camera Mode > HighSpeed, for optimal performance, use a Camera Gain of 20 or higher and adjust the Exposure Time and Excitation Intensity for your experiment.

When a luminescence experiment follows a fluorescence experiment, cycle the camera temperature to eliminate ghost images that can be created during the fluorescence assay. Select **Instrument > Manual Operation > Cycle Camera Temperature** to choose this option. The camera warms up to room temperature to release ghost images, before cooling back down to approximately -70°C (-94°F).

When using the HighSpeed camera mode or the Sensitivity camera mode, the camera behaves differently than in the Normal camera mode. The HighSpeed and the Sensitivity camera mode use 2x2 pixel binning in the camera, which combines 4 pixels into one. This is done to both increase sensitivity to smaller signals, as well as for speed. The effect of overexposure in the camera sensor when using this binning mode is different than for a non-binned exposure, due to the design of the sensor. For more details see [Determining Protocol Saturation Levels for the HS EMCCD Camera on page 139](#).

Stabilizing the HS EMCCD Camera

When you use the HS EMCCD camera, the optional stabilization process can be run before your experiment starts to avoid camera signal drift.

The Stabilizing camera process takes 50 seconds to complete.



Tip: You might be prompted to run camera stabilization, but if it is not important for your protocol, you can skip it and run your experiment immediately.

If your biology is time-sensitive you can ready the instrument before running your experiment by selecting **Instrument > Manual Operation > Pre-Stabilize**.

If your biology is not time-sensitive, within your protocol settings, you can select **Auto Stabilize** to start before your experiment begins. See [Edit Read Mode on page 73](#).



Note: When **Auto Stabilize** is saved as part of your protocol, if the camera is not already stabilized and your protocol has multiple reads, **Auto Stabilize** only runs the first time.

Your instrument maintains the camera stability for a limited time. Look in the Instrument Status tab to view the stabilization timer countdown. As long as your protocol is unchanged, and the instrument door is closed, camera stability is sustained for two hours between experiment runs. The two hour expiration countdown restarts after every experiment run.

When you open the instrument door, you have up to 15 minutes with the instrument door open before you have to run the camera stabilization process again. Within the 15 minute time limit, after you close the door, the timer returns to the two hour expiration countdown.

Understanding HS EMCCD Camera Protocol Saturation

When setting up a protocol, the read exposure settings must be set to avoid oversaturation on the HS EMCCD camera sensor. If well signals hit saturation, then they cannot be quantitatively compared. The camera provides a saturation warning when 5% of well values on the plate exceed 90% of the camera sensor dynamic range. So you want to keep your maximum signals below this level. Typically, the HS EMCCD camera threshold is about 45000 counts; but it can vary a few thousand counts between cameras. Use the Protocol Signal Test dialog to determine your camera saturation level and to adjust your exposure accordingly for your biology.

As the exposure increases past the saturation threshold, you get a saturation warning, and the raw image in the Protocol Signal Test dialog shows the wells becoming bright white, as before with Normal camera mode. See [Figure 2-8](#) and [Figure 2-9](#).

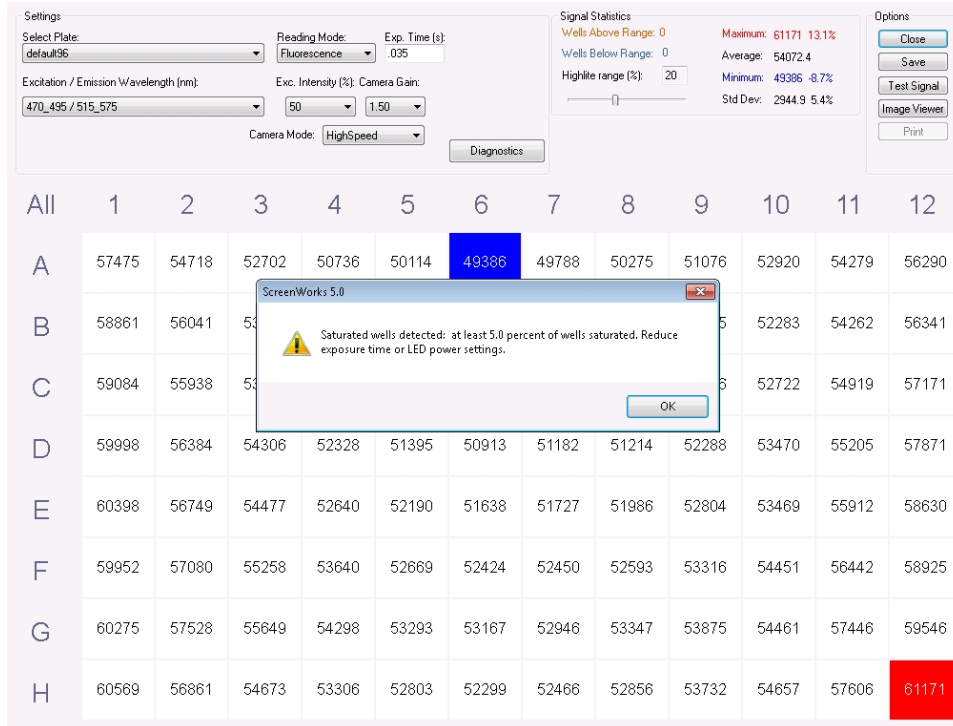


Figure 2-8: HighSpeed Mode Signal Test Saturation With Warning

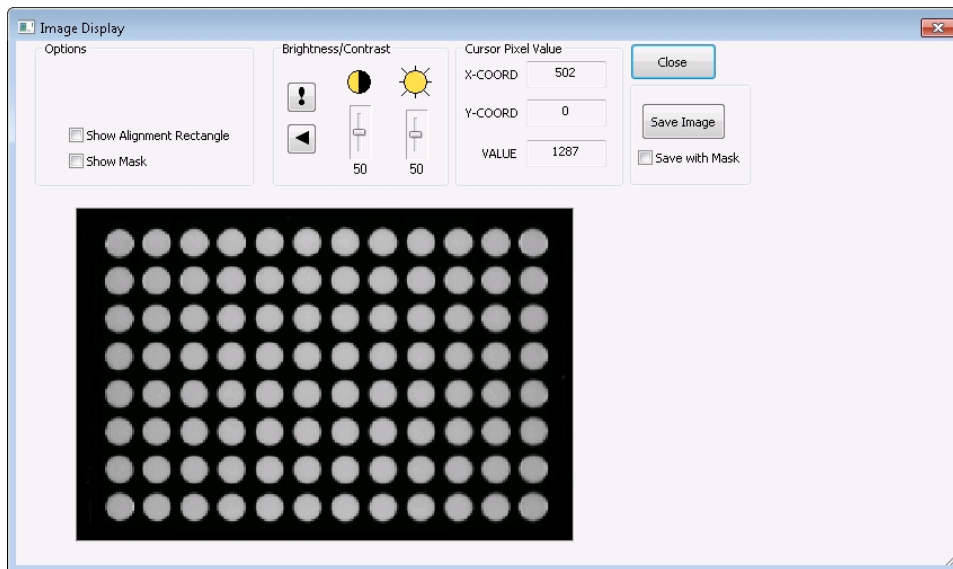


Figure 2-9: Image Display of HighSpeed Mode Signal Test Saturation

For a short range of over exposure the saturation warning works; but as the exposure increases further, the brightest pixel values start to get darker in the center of the well areas, as shown in Figure 2-10.

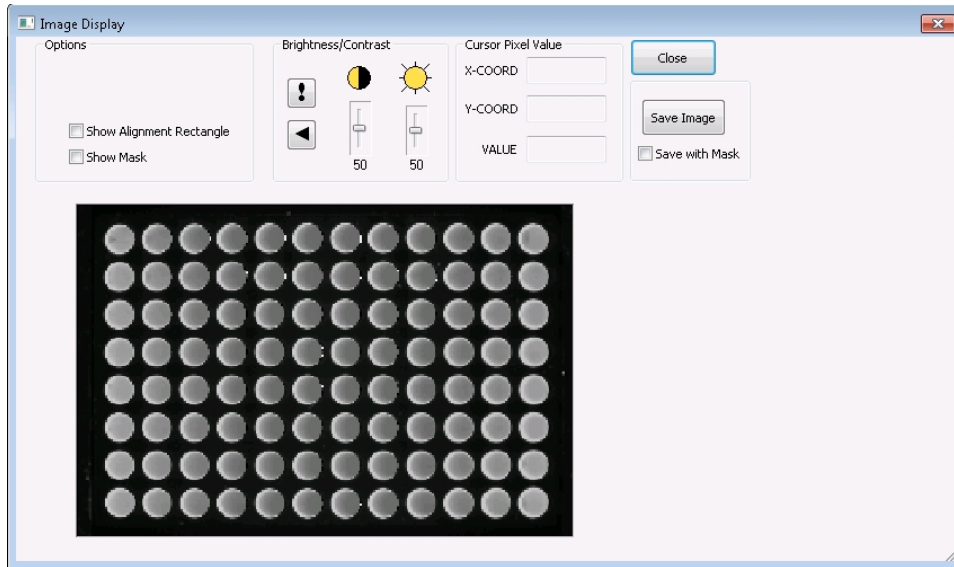


Figure 2-10: Image Display of HighSpeed Mode Signal Test Saturation Reversing

As the brightest pixels get darker, the pixel values drop lower, instead of staying at the maximum saturation value. This reversal effect happens when the binned image pixels are overexposed. The overall average for the image decreases, and after about a 4x overexposure increase, no saturation warning appears. As the exposure increases further, the wells start to bloom and flood across the image as before, except that the center of the streaks are darker, and no saturation warning appears, as shown in [Figure 2-11](#).

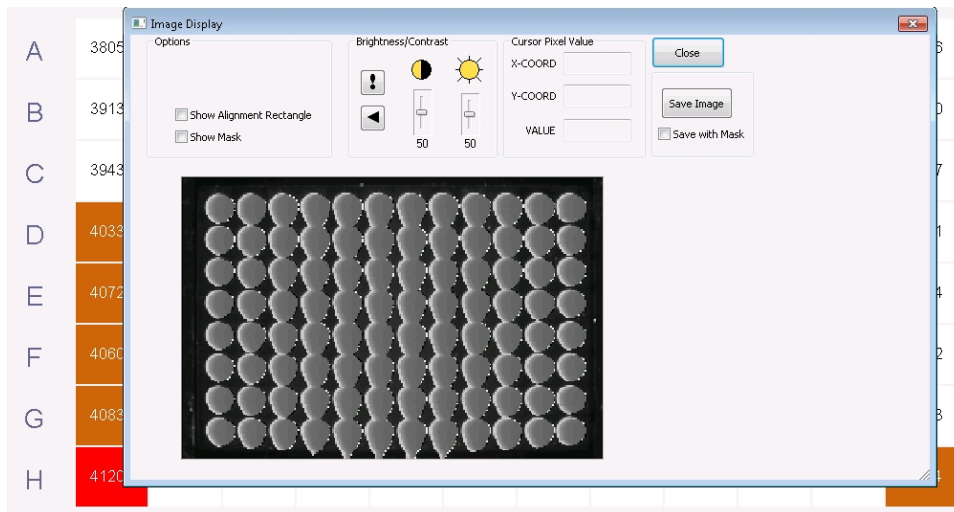


Figure 2-11: Image Display of HighSpeed Mode Signal Test Blooming Without Saturation Warning

This is a problem for detecting when the image measurements are saturated, and the system cannot give a reliable saturation warning for grossly overexposed well signals when using HighSpeed camera mode.

For adjusting the exposures for your specific biology, see [Determining Protocol Saturation Levels for the HS EMCCD Camera on page 139](#).

ICCD Camera

The camera is mounted directly beneath the read plate on the five-position stage, but at a slightly different height than the EMCCD cameras. This camera is thermoelectrically cooled and requires about five minutes to reach its operating temperature range between -20°C (-4°F) $\pm 5^{\circ}\text{C}$.



CAUTION! The instrument does not let you use the camera before it has reached proper operating temperature. Check the camera temperature on the Instrument Status panel before starting an experiment.

Similar to the Standard EMCCD camera, the camera also uses frame-transfer technology that shifts pixel values from the exposed portion of the chip to the data processing portion. In the ICCD camera, the signal is amplified in the intensifier before reaching the CCD chip. Using this method there is less noise and therefore the camera is significantly more sensitive than the Standard EMCCD camera.

When using the ICCD camera, images are taken of the bottom of the plate, amplified in the intensifier, and transferred through a fiber optic taper to the CCD chip. The Gain control of the intensifier allows for a bright signal from fluorescence as well as a dim signal from luminescence to be enhanced accordingly, so as to provide the best signal and not saturate the CCD chip below. In Fluorescence mode the Gain is preset to 2000, whereas in Luminescence mode, the Gain defaults to 280,000, but can be lowered if the luminescence assay is very bright.

This camera has a Gate Open % feature, adjustable in Fluorescence mode only. This feature controls how long the intensifier is on for each of the frames collected during the exposure time. This controls the signal intensity of the assay.

The recommended value range for the ICCD camera is 40,000 to 50,000 (at maximum signal). This camera does not require temperature cycling when changing between Fluorescence and Aequorin modes.

All Cameras

When calculated, relative luminescence units (RLU) and relative fluorescence units (RFU) are displayed in real time in the ScreenWorks Software **Analysis** process window (within the limits of computer processing speed). Data for one wavelength (for example, **Read** mode), for all 1536-wells, 384-wells or 96-wells of the read plate, display in the Multi-Well Graph on the right side of this view. The **Detail Graph** can be populated by selecting wells in the Multi-well graph. When an experiment is completed, post-assay analysis can be done via ScreenWorks Software, or data can be exported.

In normal operation images are discarded after relative luminescence units (RLU) and relative fluorescence units (RFU) values are measured. However, for quality control purposes, users can define the number of images per **Read with Transfer Fluid** step. Up to 100 images per experiment can be retained. These images can be useful for troubleshooting problems, such as cells lifting from well bottoms during compound addition. Images are saved as *.tif files with the same name and to the same directory as the data file. They can be viewed by clicking the **Image** button in the Analysis process page when the resulting data file is open in ScreenWorks Software.



Note: The robustness of an assay is not dependent on the size of the signal. It is better determined by the signal-to-noise ratio. A commonly used calculation for determining assay robustness is the Z' factor equation.

LED Modules

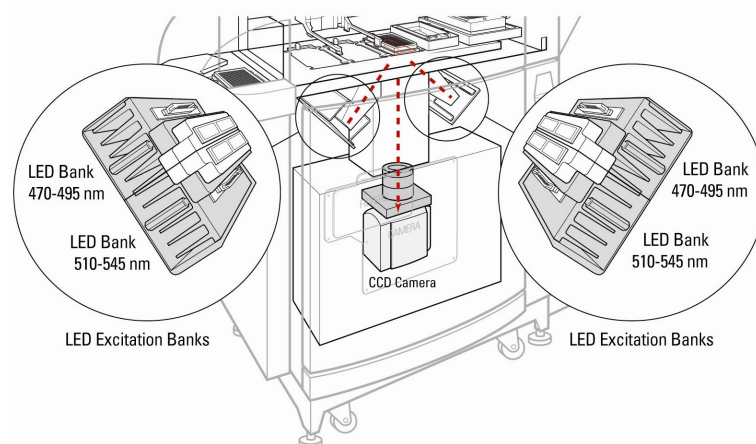


Figure 2-12: LED Configuration

The FLIPR Penta System has a total of four LED banks providing illumination for plate reading. The LED banks pulse on, two at a time, only when an image is to be captured, protecting cells from possible dye photo-bleaching.

The LED banks are divided between two modules, one on either side of the read plate position above. LED banks in corresponding positions in either module are paired. The paired LED banks pulse simultaneously during an experiment, so that light strikes the read plate base from two directions, helping to ensure that the entire plate is maximally illuminated.

Typically, the two LED bank pairs are set up with LEDs of different wavelengths, for example, one pair might have LEDs of range 470 nm to 495 nm, while the other pair might be 510 nm to 545 nm. Ratiometric experiments can be set up to use both of these wavelengths, in which case the paired banks fire alternately.



Note: Despite the two excitation wavelengths firing alternately, output data files show time points for each as occurring simultaneously.

Deflectors around the LED banks direct all light from the LEDs through excitation bandpass filters that further refine the wavelength. The light is then funneled into light pipes that focus it onto the base of the read plate.

The FLIPR Penta System does not operate without the full complement of LED banks installed, however blank LED banks can be used for one pair if only one excitation wavelength is available. You can change the LED banks, if needed. See [Exchanging LED Modules on page 198](#).

Unless you buy additional LED banks, your FLIPR Penta System includes one set of default calcium LED banks (470 nm–495 nm) and one set of blank LED banks.

Use the ScreenWorks Software Protocol builder Settings screen to configure the LEDs for your experiment. See [Setup Read Mode on page 73](#).

No LED warmup time is required before running an experiment. Start time is only dependent on the time it takes for the camera to cool down and for the stage to heat up, if the option is used.

The LEDs are air-cooled by fans, however the light output varies slightly as they heat up. To help with the heat transfer, a piece of foam is inserted in the back of the LED bank on each side. Also to counteract the temperature change, a temporal correction is automatically applied to the LED feedback circuit to normalize the system.

Flat field calibration is automatically applied to the read plate to adjust for non-uniformity of illumination across the plate. To manually calibrate the system, see [Calibrating Optics on page 131](#).



WARNING! Do not look into LED banks when turned on, especially at intensities over 30% or in the UV spectrum, unless you are viewing them through the observation panel. If light is seen escaping the instrument when the LED modules are turned on, shut down immediately and call Molecular Devices Technical Support.

Emission Filters

A three-position filter slider holding up to three 60 mm diameter interference filters is located in front of the CCD camera. The slider can be alternated in front of the camera to separate out the emission band of the dye being used. These filters can be used with a single excitation wavelength or can be paired up with additional emission filters and LEDs in a ratiometric experiment, for example, as excitation LEDs alternate between two different wavelengths, filters change at the same time so that each image taken by the camera matches the right emission filter with the excitation LED bank. The most common FLIPR Penta System configuration is a LED excitation wavelength of 470–495 nm with a 515–575 nm bandpass emission filter. For luminescence experiments it is also possible to run without the filter.

Emission LEDs and emission filters are user-changeable in approximately 5 minutes. See [Exchanging Filters on page 201](#). When installed, filters are mechanically sensed and the filter configuration can be viewed in the ScreenWorks Software, in the **Instrument Configuration** panel. If they are used in an experiment, the instrument prompts you to calibrate a new emission filter to correspond to an LED module.



Note: If the needed filter is not available, a custom filter cassette (P/N 5018841) is available for purchase for use with filters created by an outside vendor. When installed, these filters appear in ScreenWorks Software as Custom 1, Custom 2 and Custom 3.

Fluorescence Mode

In fluorescence mode the system LEDs illuminate the bottom of a 1536-well, 384-well, or 96-well ‘read’ plate containing cells loaded with fluorescent dye, and measure the fluorescence in each well. By taking a sequence of measurements in conjunction with compound application, changes in fluorescence emission characteristics due to the change in ion concentration, for example, Ca²⁺, TI⁺, or Membrane Potential.

Light-emitting diodes (LEDs) in the FLIPR Penta System produce light at distinct wavelength ranges to excite the fluorescent dye that has been added to the cells in the read plate wells. The entire well plate bottom is illuminated. Fluorescent light emitted by the dye—again, for the entire plate—passes through an emission filter before being captured in a CCD camera, high-speed, high-sensitivity HS EMCCD, EMCCD or ICCD. Fluorescence is measured from each well independently, and converted into a numerical value. The FLIPR Penta System can be configured with two LED banks and three emission filters, allowing the software configuration of up to four excitation/emission wavelength combinations (‘read modes’). Up to four different fluorescence effects can be measured within a single experiment.

Luminescence Mode

The FLIPR Penta System also provides luminescence mode with the HS EMCCD camera option and the ICCD camera option. The instrument has a light-tight enclosure to operate in luminescence mode. A specialized high-sensitivity HS EMCCD camera can be installed in place of the standard EMCCD camera to detect signal at both high-speed fluorescence or Aequorin (Luminescence) mode. The camera is mounted directly beneath the read plate, so images are taken of the entire bottom of the plate. For cell suspension experiments, an integrated Cell Reservoir allows uniform cell suspension to be pipetted directly into the read plate. From the 3 filter positions available on FLIPR Penta System, it is recommended to have one open position, so no filter is used during the luminescence assay.

As with fluorescence, luminescence is measured from each well independently, and converted into a numerical value.

Chiller

The chiller sits outside of the FLIPR Penta instrument and is connected with a cable and tubing. It uses a special provided cooling liquid, and the instrument embedded computer and firmware control it.



Figure 2-13: The Chiller

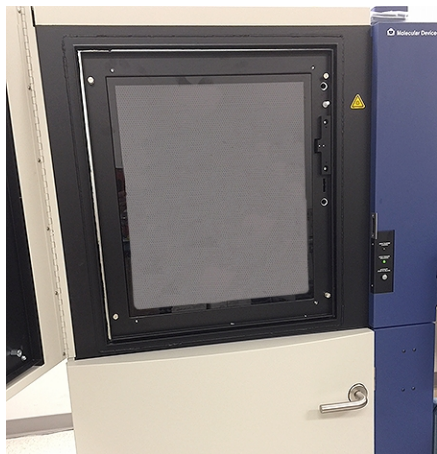
Because the FLIPR Penta System is a light-tight instrument, there is very limited airflow inside the instrument enclosure. The external chiller provides the operating environment requirement for all of the available camera options.

Observation Panel

Use the observation panel while testing or troubleshooting the FLIPR Penta System to view hardware movements in the upper top compartment.

Under normal operating conditions, the upper door must be closed to run an experiment, and keep light out of the chamber. When the observation panel is mounted onto the chamber, the door can be left open, allowing you to view movements of the pipettor and FLIPR Cycler plate handler. Normal instrument control comes from the ScreenWorks Software. In Remote mode SynchroMax ET plate handling, or third-party plate handling software controls the instrument.

To mount the panel, remove it from the door and attach it with the four captive thumbscrews to the top compartment frame.



When not in use, the observation panel attaches to the inside of the upper door.



To acquire quality data, re-attach the observation panel to the inner door before running an experiment.



CAUTION! If pretending to run in luminescence mode with the ICCD camera, DO NOT touch the white door switches. Room light damages the Intensifier. The door switches detect the open door to protect the camera.

Only use the observation panel to view internal pipettor movement or FLIPR Cycler movement.





CAUTION! Do not use it during experiments when data is being acquired.

This chapter provides an overview of the ScreenWorks Software functionality, including the following topics:

- [Online Mode vs. Offline Mode](#)
- [User Interface Overview on page 48](#)
- [Menu Bar on page 50](#)
- [Experiment Window on page 65](#)

Online Mode vs. Offline Mode

The software has two startup modes:

-  Online—the software is connected to the instrument.
-  Offline—the software is disconnected from the instrument.

The default start-up mode is determined during software installation in the **Online/Offline** dialog. Regardless of the startup mode, pipettor head and tip washer type must always match.



Note: To be able to select either mode while generating protocols, you must install the software in Offline mode.

When the software is open, you can switch modes by selecting from the **Instrument** menu, **Go Online** or **Go Offline**.



Note: Switching modes after the software is opened does not change the software startup mode chosen at the time of installation. To change the default startup mode, the ScreenWorks Software must be reinstalled.

Online Mode

When started in Online mode, ScreenWorks Software checks for instrument connections. If no connections are sensed, you are notified. You can then either check the connections and attempt to connect again, or choose to run the software in Offline mode.

Create new protocols in Online mode.

When ScreenWorks Software starts in Online mode and connects to the instrument, the default installation configuration file is overwritten using the current instrument settings and plate library information.

If you are working in Online mode and then switch to Offline mode, the instrument setup configuration is remembered as the last setting.

Offline Mode

When ScreenWorks Software starts and no instrument connection registers, you are notified and can then either check the connections and attempt to connect again in Online mode, or choose to run the software in Offline mode.

When ScreenWorks Software starts in Offline mode, you can configure the following hardware options:

- Camera Type
- Pipettor Format

Protocols created in Offline mode with hardware settings that do not match current hardware settings are flagged. You must change the hardware settings to match those in the protocol in order to run it.

User Interface Overview

The ScreenWorks Software main screen includes menu and toolbar across the top, and a status bar at the bottom. The main working area in the middle has two sections:

- Experiment window—the central proportion of the main window, for protocol configuration, data viewing and analysis.
- Instrument Status-Configuration-Process panel—at the bottom of the main work area, for displaying instrument information and creating protocols. This panel contains the Instrument Status tab, Instrument Config tab, and the Process tab.

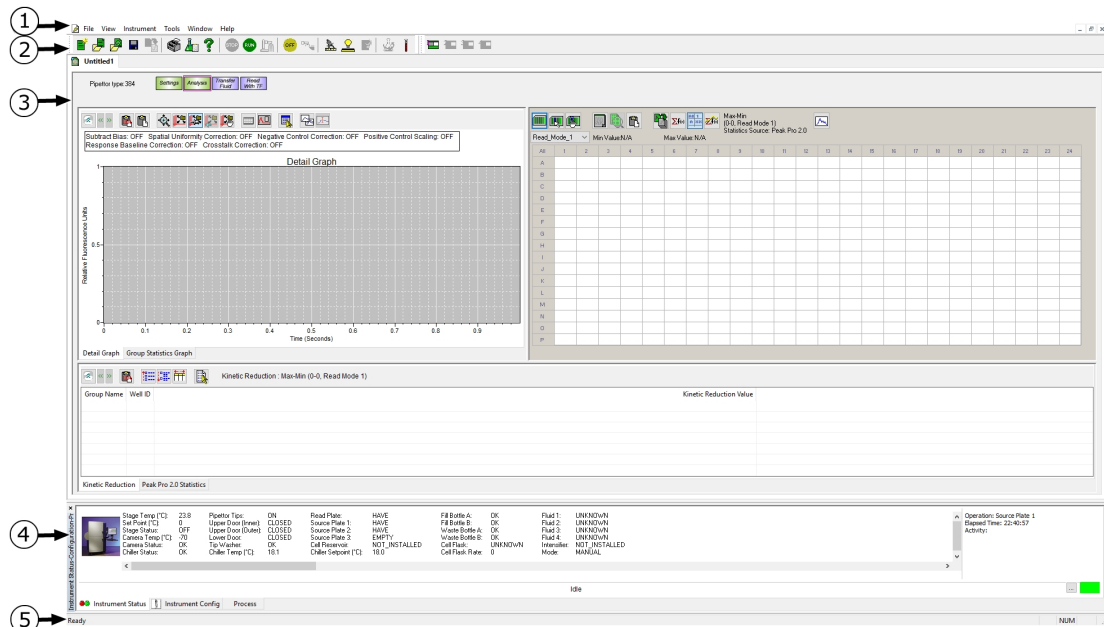


Table 3-1: Main Window Layout

| Item | Description |
|------|-------------|
| 1 | Menu Bar |


Table 3-1: Main Window Layout (continued)

| Item | Description |
|------|---|
| 2 | Toolbar |
| 3 | Experiment Window |
| 4 | Instrument Status-Configuration-Process Panel |
| 5 | Status Bar |

Menu Bar

The menu bar contains the following six sub-menus that group together related functions:

- File Menu
- View Menu on page 52
- Instrument Menu on page 53
- Tools Menu on page 58
- Window Menu on page 62
- Help Menu on page 62

 **Tip:** Some menu options can be opened with a keyboard shortcut using the **ALT** key to underline the letter in each menu title that is used to open the menu; for example, when you click **ALT** the **I** in the **Instrument** menu is underlined. Click the **I** key and the **Instrument** menu opens.

File Menu

The **File** menu enables you to select the following options:

Table 3-2: File Menu Options

| Option | Description | Keyboard Shortcut |
|--------|---|------------------------|
| New | Opens a new protocol (*.fmp) in the Experiment window, with default settings for one read with fluid transfer. New protocols are named <i>Untitled[n]</i> , where <i>n</i> is a number. | Ctrl+N |
| Open | Opens an Open File dialog to open a saved Protocol (*.fmp), Data (*.fmd) or Image (*.png) file, as selected in the submenu. By default, the dialog opens in the folder set in Tools > Set Default Directories . Note: Only data files created with ScreenWorks Software can be opened in ScreenWorks Software. Data files from previous FLIPR Tetra Systems (versions 1.X through 2.X) cannot be opened. | Ctrl+O Ctrl+Shift+O |
| Close | Closes the currently active protocol or data file displayed in the Experiment window. If modifications have been made to the file, you are prompted to save the modifications. | Alt+F,C |
| Save | Saves the currently active protocol or data file, displayed in foreground of the Experiment window. Also saves if the displayed file is a default protocol that is untitled, the Save As dialog appears so that you can name the file. | Ctrl+S |

Table 3-2: File Menu Options (continued)

| Option | Description | Keyboard Shortcut |
|----------------------|--|-------------------|
| Save As | Opens the Save As dialog and allows you to save a protocol or data file under a new name or format (for example, save a data file as a protocol). Also allows you to save a protocol or data file as an earlier version of the ScreenWorks Software. When changing formats, select the file type from the Save as type drop-down list. | Alt+F,S |
| Save All Files | Saves all of the opened protocol and data files. If the default protocol file was opened, the Save As dialog is displayed so that you can name the files. | |
| Close All Files | Closes all of the opened protocol and data files. If any of the files had changes since they were opened or the default protocol file was opened, the Save or Save As dialog is displayed so that you can name the files. | |
| Export | Opens the Export File or Batch Export dialog, as selected in the submenu, to manually export data from the open data file, or other data files on disk; see Exporting Data on page 157 . | |
| Page Setup | Opens the Page Setup dialog to configure the printer and print settings for the document. | Alt+F,G |
| Print Preview Report | Displays how the document looks before printing. | Alt+F,V |
| Print Report | The Print Report dialog allows you to select the graphs or reports to print from the current data file. Having made the selection you can open Print Preview Report to check before printing, or open the Print dialog to print. | Crtl+P |
| 1–6 Data Files | Lists the six most recently opened data files, with the most recently opened at the top. | |
| 7–10 Protocol Files | Lists the four most recently opened protocol files, with the most recently opened at the top. | |
| Exit | Closes ScreenWorks Software. If you have unsaved data, you are prompted to save it before closing. | Alt+F,X |

Saving Data Files as Protocol Files

When the active file is a protocol file, you can add, remove, or change processes in the file, then save the amended protocol and run it. Protocol information stored in data files cannot be edited, nor used to run a new experiment, however it is possible to extract this information to a new protocol file. Experiments run using this file have exactly the same steps as the steps used to create the data file.

To store a data file as a protocol file, select **Save As**, assign a protocol name, select **.fmp* as **File Type** and click **Save**. The stored file is stripped of all data and only associated protocol information is stored in the protocol file.



Note: Saved changes only affect the protocol. The data file from which the protocol was derived remains intact.

View Menu

The **View** menu options enable you to select the following options:

The keyboard shortcut is **ALT+V**.

Table 3-3: View Menu Options

| Option | Description |
|--------------------|---|
| Experiment Setup | Displays the Experiment Setup view of a protocol or data file, showing processes and associated dialogs in the Experiment window. |
| Experiment Summary | Displays in the Experiment window the Experiment Summary for the current protocol or data file in the Experiment window. To return to the Experiment Setup view, select View > Experiment Setup . |
| Save Layout | Saves the proportions displayed in the Experiment window (including Multi-Well and Detail Graphs as well as Group Statistics window) and Instrument Status panel to one of four layouts. These layouts can be toggled between when the appropriate Restore Layout selection is made. |
| Restore Layout | Restores the screen layout to one of four proportions saved in the Saved Layout selection. |
| Instrument Status | Toggles the Instrument Status panel in and out of view. See Instrument Status Tab on page 67 . |
| Toolbar | Toggles the Toolbar in and out of view. See Toolbar on page 62 . |
| Status Bar | Toggles the Status Bar at the bottom of the main window in and out of view. See Status Bar on page 71 . |

Instrument Menu

The **Instrument** menu commands enable you to access instrument communication and manual dialogs.

Table 3-4: Instrument Menu Commands

| Options | Description |
|-------------------|--|
| Go Online/Offline | <p>Switches the instrument between Online mode and Offline mode. See Online Mode vs. Offline Mode on page 47.</p> <p>Go Online—Displays when ScreenWorks Software is offline. Click to connect the software to the instrument, giving ScreenWorks Software control of the instrument.</p> <p>Go Offline—Displays when ScreenWorks Software is connected to the instrument. Click to disconnect the software from the instrument.</p> |
| Run Experiment | Instructs ScreenWorks Software to start an experiment using the uppermost protocol file in the experiment. |
| Stop Experiment | <p>Instructs ScreenWorks Software to stop the experiment.</p> <p>Note: Stop Experiment should only be used in emergencies to halt an experiment. If used, an instrument reset may be required.</p> |
| Manual Operation | <p>Commands in this submenu control specific hardware operations.</p> <p>Note: The commands are disabled if the associated hardware is not available.</p> <p>Load Tips—Instructs the pipettor to load tips or pin tool from the stage Position 1.</p> <p>Unload Tips—Instructs the pipettor to unload tips or pin tool to the stage Position 1. Unload tips to a tip rack, otherwise, make sure there is a container at stage Position 1 to receive tips.</p> <p>Wash Tips—Instructs the pipettor to wash tips or pin tool. A dialog opens for you to configure the wash.</p> <p>Drain Wash Basin—Enables you to drain the wash basin to the waste bottles when there is an instrument error.</p> <p>Yellow Plate Signal Test—Coordinates the reading of the stage Position 3 read plate for the current protocol and displays the numerical results in the Signal Test dialog. For more details, see Signal Test Dialog Options on page 55.</p> <p>Protocol Signal Test—Coordinates the reading of the stage Position 3 read plate and to display the numerical results in the Signal Test dialog. Instrument settings can be saved to the open protocol. For more details, see Signal Test Dialog Options on page 55.</p> <p>Change Head—Instructs the pipettor head to move over stage Position 3 to be exchanged to a new head format. See Exchanging Pipettor or Pin Tool Head on page 186.</p> |

Table 3-4: Instrument Menu Commands (continued)

| Options | Description |
|---------------------------------|--|
| Manual Operation (continued) | <p>Cycle Camera Temperature—Cycles the camera temperature when you want to run a low-fluorescence or luminescence experiment immediately after running a high-fluorescence experiment. The temperature cycle runs for about 15 minutes, warming to room temperature, 20°C to 25°C (68°F to 77°F), and then cooling to operating temperature: EMCCD camera: -70°C (-94°F) ± 2°C HS EMCCD camera: -70°C (-94°F) ± 2°C ICCD camera: -20°C (-4°F) ± 5°C</p> <p>Temperature Control—Turns heating on and off for stage source plate Positions 1, 2, and 3, and enables you to specify the heating temperature.</p> <p>Cell Flask Stirring Control—Enables you to set the stir speed rate for the cell flask.</p> <p>Wash Cell Reservoir—Enables Cell Reservoir washing by selecting the Fluid Source, Fluid Destination, Fill Speed, Drain Speed, along with the number of Wash Cycles and Hold Time. It also enables you to pre-coat the tubes, which is recommended for the first run with any cells in suspension.</p> <p>Drain Cell Reservoir—Enables you to drain the Cell Reservoir to a specified destination for manually pouring cells into the Cell Reservoir, or in case of an instrument error.</p> <p>Note: Resetting the instrument automatically drains the Cell Reservoir to the waste bottle.</p> |
| Set Remote/ Manual Mode | <p>Switches the instrument between manual and remote modes.</p> <p>Set Remote Mode—Enables you to integrate a third-party robotics system with the FLIPR Penta System. See Robotic Integration on page 261.</p> <p>ScreenWorks Software automatically goes into Remote Mode when you open SynchroMax Automation.</p> <p>Set Manual Mode—Disables any robotics control, and switches instrument communication to the ScreenWorks Software.</p> |
| SynchroMax Automation | <p>Opens the SynchroMax dialog for configuring plate handling with the SynchroMax ET. See SynchroMax™ Automation on page 57.</p> |
| Reset | <p>Reinitializes the system to clear any fatal errors as designated in red at the bottom of the Instrument Status panel.</p> |
| Clear Error | <p>Clears minor system errors indicated in yellow in the Instrument Status panel. See Instrument Status Colors on page 209.</p> |
| Calibration | <p>Opens the Calibration dialog where Flat Field Calibration Calibrations can be done. See Calibrating Optics on page 131.</p> |
| Refresh Configuration | <p>Refreshes the instrument configuration after you change hardware settings.</p> |

Signal Test Dialog Options

The signal test has two functions:

- Checking the state of the overall system: This test is typically run using the yellow test plate with the respective plate format for your assay. Open the dialog with the **Yellow Plate Signal Test** option in the main menu. Outlined in the table below are the default settings used when performing the yellow plate signal test. These settings cannot be saved with a protocol for later use.
- Checking initial fluorescence of a plate before running an assay: This test is typically run to evaluate the assay plate before running an experiment. Open the dialog with the **Protocol Signal Test** option in the main menu. Settings outlined in the table below reflect the Normal camera mode and default to the settings in the currently open protocol when running the protocol signal test. These settings can be saved in the open protocol for later use.

Table 3-5: Yellow Plate Signal Test and Protocol Signal Test Setting Options

| Options | Description |
|--------------------------------|--|
| Select Plate | Choose the plate type from the drop-down list. |
| Excitation/Emission Wavelength | Select the appropriate excitation/emission wavelength pair for the signal test from the drop-down list. Note: Only calibrated excitation/emission wavelength pairs are displayed in the drop-down list. |
| Reading Mode | Select Fluorescence or Luminescence from the list. |
| Camera Gain | Select the camera gain for the signal test from the list. For the 470–495/515–575 nm excitation/emission pair when reading a yellow test plate: Standard EMCCD Camera is 50 ICCD camera is 2000 HS EMCCD camera is 8 when in Camera Mode > Normal, and is 20 when in Camera Mode > HighSpeed |
| Camera Mode | In Reading Mode > Florescence: Normal, Sensitivity, and HighSpeed for the HS EMCCD camera only In Reading Mode > Luminescence: Normal and Sensitivity |
| Gate Open | This option is available only for the ICCD camera. Typical Gate Open values for Fluorescence are around 6% but can be set to 100%. Luminescence gate is always 100%, and not user-adjustable. |

Table 3-5: Yellow Plate Signal Test and Protocol Signal Test Setting Options (continued)

| Options | Description |
|----------------------|--|
| Excitation Intensity | Select the LED intensity for the signal test from the drop-down list. Excitation intensity is scaled as a percentage of the total LED output (0–100%). For the 470–495/515–575 nm excitation/emission pair when reading a yellow test plate: Standard EMCCD Camera is 80 ICCD camera is 30 HS EMCCD camera is 50 |
| Exposure Time | Enter the amount of time (in seconds) to keep the camera shutter open during the Signal Test. For the 470–495/515–575 nm excitation/emission pair when reading a yellow test plate: Standard EMCCD Camera is 0.4 seconds ICCD camera is 0.53 HS EMCCD camera is 0.005 |
| Highlight Range | Highlights well values that lie within the set statistical range. Set the statistical range by using the slider. |
| Wells Above Range | Displays the number of wells above the statistical range as determined by the Highlight Range. |
| Wells Below Range | Displays the number of wells below the statistical range determined by the Highlight Range. |
| Maximum | Displays the largest value on the signal test plate. |
| Average | Displays the average value on the signal test plate. |
| Minimum | Displays the smallest value on the signal test plate. |
| Std. Dev | Displays the standard deviation of the signal test plate. |
| Test Signal | When clicked, initiates a new signal test. |
| Save | When clicked, saves the signal test as an ASCII text file (*.sig). |
| Print | When clicked, prints the signal test. |
| Image Viewer | Displays the CCD image showing where relative fluorescence units (RFU) values were derived for the signal test. Show > Hide Mask —Shows or hides the mask used in the test, indicating the pixel area used to derive the RLU value for each well. View Image —Refreshes signal test image. |
| Diagnostics | For Technical Support use only. |

For procedures, see [Running Signal Tests on page 135](#).

SynchroMax™ Automation



Note: Molecular Devices no longer offers the SynchroMax ET external plate handler robot. The following information is for reference only.

The **SynchroMax Automation** option is only enabled when you have a SynchroMax ET external plate handler installed.

When you open the **SynchroMax** dialog from **Instrument > SynchroMax Automation**, ScreenWorks Software automatically goes into **Remote Mode**, giving control of the instrument to the SynchroMax ET software. ScreenWorks Software returns to **Manual Mode** when the **SynchroMax** dialog closes.

To run a series of experiments with the SynchroMax ET plate handler, first open the protocol file that you want to use in ScreenWorks Software. Only one protocol is used in a run for as many times as the SynchroMax ET plate handler delivers plates.

Having prepared the protocol, open the **SynchroMax** dialog and select a **Stack Layout Template**. These files specify, for each plate location in the five-position stage, how many plates load to that position, from which stack, and where they are moved after use. A number of **Stack Layout Templates** are supplied—select the one that fits your assay.

When a **Stack Layout Template** is open in the dialog, the graphic in the dialog displays on which stacks the plates load. You can use this as a guide for loading the plates. All plates in a single stack should be of the same type.

If there are any stage positions that use the same plate throughout the duration of the run, these stage positions should be loaded manually before the run begins. Stage positions that have plates brought to them during the run must be empty at the start of the run.

Before selecting **Run**, make sure all output racks are empty prior to beginning the experiment. When plates are in position and the FLIPR Penta System ready, click **Run** in the **SynchroMax** dialog. After checking that the plate configuration is compatible with the protocol **Settings** configuration, the instrument runs until all plates have been used. During this time the **Run** button changes to **Stop**, so you can stop the run before it completes if necessary.



Note: **Stop** should only be used in emergencies to halt an experiment. If used, an instrument reset may be required.

When the SynchroMax ET plate handler is active, the system clears all plates at the end of an experiment. The SynchroMax plate handler remembers only those plates it loaded during the experiment and checks if any of those plates remain when the experiment completes. If so, they are removed. A reagent reservoir can be used in a source position within the FLIPR Penta System as long as it is manually loaded and the SynchroMax template does not include the loading or removal of plates to that position.

Unless you need to run another set of experiments, click **Done** to close the dialog and return ScreenWorks Software to **Manual Mode**.

Tools Menu

| Options | Description |
|-------------------------|---|
| Set Default Directories | <p>Opens the Set Default Directories dialog to designate directories in which to store protocol and data files, signal tests, group templates and to which export files are written.</p> <p>The Open > Protocol, Data and Image File commands in the File menu open to the folders set here, as do the File > Export dialogs. In addition, Group Templates *.fmg can be exported and imported from the default folder defined in the directory.</p> <p>The default protocol directory is: C:\Documents and Settings\[your_user_name]\My Documents\Molecular Devices\ScreenWorks\MyProtocols [MyData, MySignalTests, or MyGroupTemplates]</p> <p>The default data and export directories are: C:\Documents and Settings\[your_user_name]\My Documents\Molecular Devices\ScreenWorks\MyData \</p> <p>The default signal test directory is: C:\Documents and Settings\[your_user_name]\My Documents\Molecular Devices\ScreenWorks\MySignalTests</p> <p>The default group template directory is: C:\Documents and Settings\[your_user_name]\My Documents\Molecular Devices\ScreenWorks\MyGroupTemplates</p> <p>Note: It is recommended that you save all protocol and data files on the local hard drive to ensure instrument function or data is not lost if your server fails during an experiment.</p> |
| Plate Library | Opens the Plate Process Definition dialog, which lists current plate definitions and allows you to add additional plate definitions to the system. See Plate Process Definition on page 59 for details. |
| Open Error Log | Opens an error log generated by the FLIPR Penta System. This feature is for technical support and requires a password. |
| Save Error Log | Save error logs. Logs are saved as *.fel (FLIPR Penta System Error Log) encrypted files. The files can be forwarded to Molecular Devices Technical Support. |
| Assay Log | <p>For data files only, opens a dialog reporting when the protocol steps were applied.</p> <p>Note: This log can be accessed by selecting CTRL+SHIFT+A.</p> |

Plate Process Definition

All plates currently in the system are listed in the **Plate Process Definition** dialog, opened from **Tools > Plate Library**. They are categorized first into read and source plates and then by well-number format.

A description of commands in the Plate Process Definition and Define Camera Parameters, both accessed via **Tools > Plate Library** is provided below. For instructions on adding a new read or source plate, see [Define Basic Plate Parameters Dialog on page 60](#). The Plate Library is camera-specific, so if any plates are added with one camera type, they are unavailable after the camera type changes.

The **Plate Process Definition** dialog is accessible only to administrators, and is password protected. The default user name is *fliprtetra* and password is *flipr*.

The following fifteen default plates are included with the system software:

- Default 96
- Default 96 small volume
- Default 96 no slit mask
- Default 96 boat
- Default 384
- Default 384 small volume 5X3 mask
- Default 384 small volume 3X3 mask
- Default 384 boat
- Default 384 blot
- Default 1536
- Default 1536 boat
- Default 1536 blot
- Default Cell Reservoir 96
- Default Cell Reservoir 384
- Default Cell Reservoir 1536

Table 3-6: Plate Process Definition Dialog Options

| Options | Description |
|--------------|--|
| Close | Closes the Plate Process Definition dialog. |
| Open Plate | Opens the Define Basic Plate Parameters dialog for the plate selected in the tree view. |
| Copy Plate | Opens the Define Basic Plate Parameters dialog allowing you to add a new plate. See Define Basic Plate Parameters Dialog on page 60 . Note: When copying a plate, make sure that you select a plate with a similar well format, as the number of wells and plate mask are transferred to the new plate. |
| Delete Plate | Deletes the plate displayed in the Plate Process Definition field. |
| Collapse All | Collapses the Plate Process Definition tree. |
| Expand All | Displays all items in the Plate Process Definition tree. |

Define Basic Plate Parameters Dialog

This dialog opens when you select **Open Plate** or **Copy Plate** in the **Plate Process Definition** dialog. Use the dialog to establish the critical dimensions of the plates you are using.



Note: Dimensions of default plates cannot be modified. However, the plate mask can be realigned.

Table 3-7: Define Basic Plate Parameters Dialog Options

| Options | Description |
|------------|--|
| Plate Name | Enter the name you want assigned to the new plate (maximum number of characters is 50). |
| Plate Type | Select Source Plate or Read Plate for the new plate. Source Plate —Creates a plate that does not include a plate mask. These plates are typically used for compound storage. Read Plate —Use for all plates that are read in the read position and require a plate mask. |

Table 3-7: Define Basic Plate Parameters Dialog Options (continued)

| Options | Description |
|----------------------|---|
| Plate Specifications | <p>Area for entry of physical plate dimensions. Refer to the diagram in the upper right corner of the dialog.</p> <p>For best results, obtain the plate dimensions from the plate manufacturer. See FLIPR Penta Plate Dimensions on page 241.</p> <p>Note: Molecular Devices is not responsible for instrument malfunctions if plate specifications are not correct.</p> <p>Rows—Enter the number of rows on the plate. Columns—Enter the number of columns on the plate. Well Shape—Select the appropriate shape from the drop-down list. Well Volume (μL)—Enter the maximum well volume in microliters. X (mm)—Enter the distance (in mm) from the left side of the plate to the center of well A1. Y (mm)—Enter the distance (in mm) from the top of the plate to the center of well A1. Bottom (mm)—Enter the distance (in mm) from the bottom of the plate skirt to the inside of the well. Top (mm)—Enter the distance (in mm) from the bottom of the plate skirt to the top of the well. Well Offset (mm)—Enter the distance (in mm) from the center of one well to the center of adjacent well.</p> |

Define Camera Parameters Dialog

When **Read Plate** is identified as the **Plate Type** in the **Define Basic Plate Parameters** dialog, the **Finish** button is enabled. Before clicking **Finish**, place a read plate with 10^{-8} M fluorescein in the read position. When selected, the instrument reads the plate and defines a plate mask for the plate definition created. If a plate is in position, but no mask can be defined, the plate definition is saved as a source plate.

Window Menu

The **Window** menu options allow you to change the way the Experiment window is viewed. The keyboard shortcut is **ALT+W**.

Table 3-8: Window Menu Options

| Options | Description | Keyboard Shortcuts |
|-----------------|--|--------------------|
| Cascade | Aligns the open windows so that they are overlapped and staggered with the title bars visible. | Alt+W,C |
| Tile | Divides the screen into as many segments as there are windows and aligns them so that they are displayed side-by-side. Each file occupies a segment and is visible. However, each of the file images is scaled according to the number of files being displayed. | Alt+W,T |
| 1–10 Data Files | Lists the open files (up to 10) | |

Help Menu

The **Help** menu options provide access to user guides and information about the software. The keyboard shortcut is **ALT+H**.

Table 3-9: Help Menu Options

| Options | Description |
|-------------------------------------|---|
| FLIPR Penta System User Guide (PDF) | Opens a PDF version of this manual that is appropriate for the version of software installed. |
| Release Notes (PDF) | Opens a PDF version of the <i>ScreenWorks Software Release Notes</i> . |
| Protocol Guide (PDF) | Opens a PDF guide that provides detailed information, procedures, and optimization guidelines for running common types of experiments. |
| Software License | Opens the Software License dialog, which enables ScreenWorks Peak Pro Software activation. |
| Contact Us | Links to the Molecular Devices website contact page. |
| About ScreenWorks | Opens the About ScreenWorks dialog, which displays the version numbers for ScreenWorks Software, the Firmware EC, Firmware Motion, and Remote Interface. |

Toolbar

The toolbar, beneath the menu bar, contains tool button shortcuts for a number of main menu commands.

The toolbar can be hidden or shown from the **View > Toolbar** toggle command.

For more detailed descriptions of the buttons, see [Menu Bar on page 50](#).

Table 3-10: Toolbar Buttons


























| Button | Name | Description |
|---|-----------------------|--|
|  | New Document | Opens a new protocol file (*.fmp) with default settings for an assay with a single read with fluid transfer. Files are named <i>Untitled[n]</i> where <i>n</i> is a number. |
|  | Open Protocol File | Opens the Open File dialog to browse and open protocol (*.fmp) files. |
|  | Open Data File | Opens the Open File dialog to browse and open data (*.fmd) or image files. |
|  | Save File | Opens the Save File dialog that allows you to save the current file in the desired location. |
|  | Export File | Opens the Export File dialog that allows you to manually export the current file in the desired location. |
|  | Print | Opens the Print dialog. |
|  | Experiment Summary | Displays protocol contents (process steps and correction) for the current protocol or data file. This button toggles with Experiment Setup . |
|  | Experiment Setup | Opens the Experiment Setup window to edit the current protocol, view or analyze data in a data file. This button toggles with Experiment Summary . |
|  | Help | Opens various documentation in PDF format. |
|  | Stop | Stops the experiment currently running. |
|  | Run | Starts the protocol selected in the Experiment window. |
|  | SynchroMax Automation | Opens the SynchroMax™ Automation window to select the desired template to run the SynchroMax ET plate handler. |
|  | Online Mode | This button is displayed when System Control Software is in Offline Mode . Click to connect to the instrument and go into Online Mode. This button toggles with Offline Mode . |

Table 3-10: Toolbar Buttons (continued)

| Button | Name | Description |
|---|---------------------------------|---|
|  | Offline Mode | This button is displayed when ScreenWorks Software is in Online Mode. Click to disconnect from the instrument and go into Offline Mode . This button toggles with Online Mode . |
|  | Remote Mode | Instructs ScreenWorks Software to disable the manual connection and to only receive Remote commands from a third-party robot. This button toggles with Manual Mode . |
|  | Manual Mode | Instructs ScreenWorks Software to disable the remote connection and to only receive Manual commands from the ScreenWorks Software user interface. This button toggles with Remote Mode . |
|  | Calibration | Opens the Calibration dialog where Flat Field calibrations can be performed. |
|  | Yellow Plate Signal Test | Opens the Yellow Plate Signal Test dialog to display the numerical results using the Yellow Plate. |
|  | Protocol Signal Test | Opens the Protocol Signal Test dialog to display the numerical results before running an experiment. Settings defined in this signal test can be saved to the protocol *.fmp file. |
|  | Set Spinner Flask Stirring Rate | Opens the Spinner Flask Control dialog where the stirring rate can be set. |
|  | Set Stage Temperature | Opens the Set Temperature dialog where the stage temperature can be set in degrees Centigrade, or disabled. |
|  | Restore Layout 1 | Restores the Experiment window layout to those defined as Save To Layout 1 in the View menu. |
|  | Restore Layout 2 | Restores the Experiment window layout to those defined as Save To Layout 2 in the View menu. |
|  | Restore Layout 3 | Restores the Experiment window layout to those defined as Save To Layout 3 in the View menu. |
|  | Restore Layout 4 | Restores the Experiment window layout to those defined as Save To Layout 4 in the View menu. |

Experiment Window

The Experiment window is the main working area in ScreenWorks Software for either:

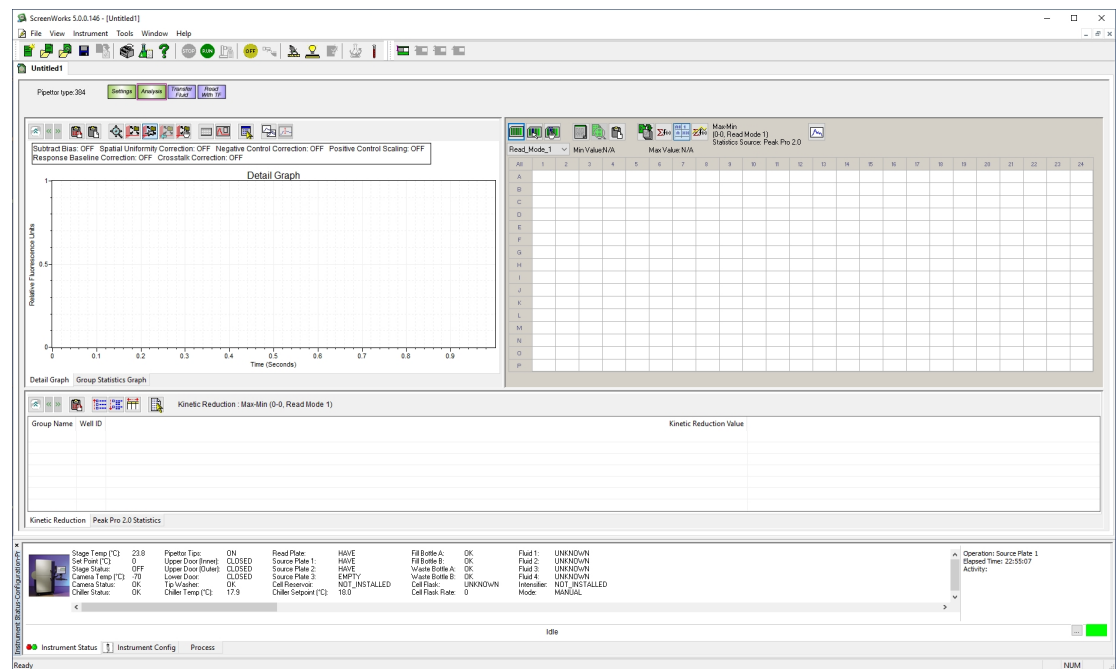
- [Experiment Setup on page 65](#)
- [Experiment Summary on page 66](#)

The Experiment window can have one protocol file and multiple data files open simultaneously. Use the options in the Windows menu to view these file panes individually, cascaded, or tiled. See [Window Menu on page 62](#).

Only the selected open file pane is active.

Experiment Setup

By default, ScreenWorks Software opens the **View > Experiment Setup** screen in the Experiment window.

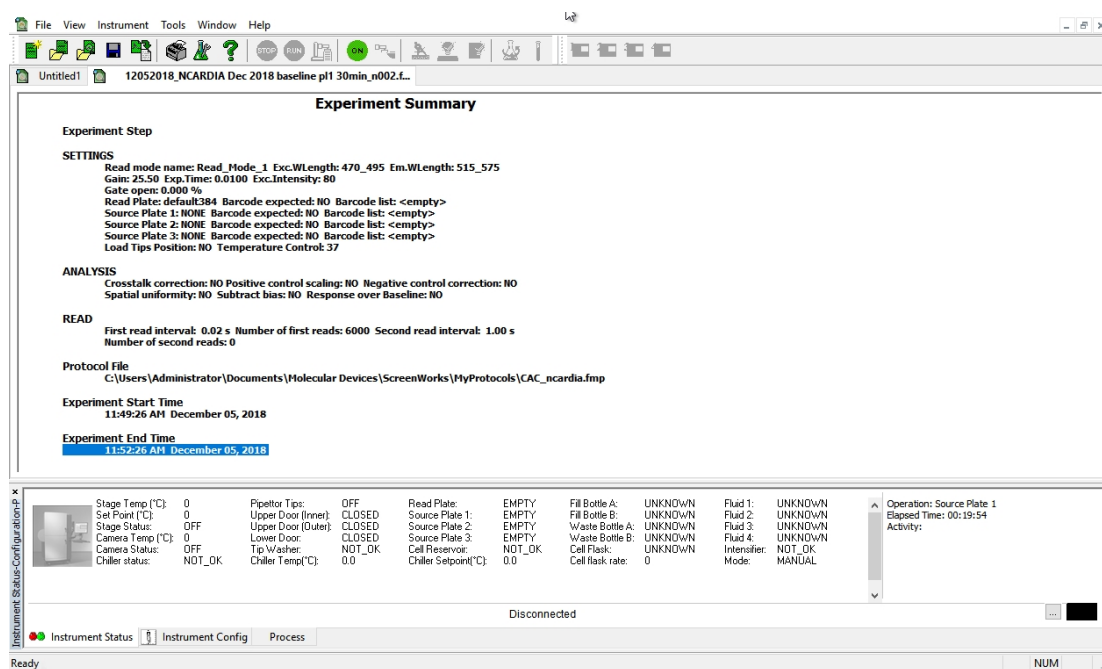


A new protocol file displays as a tab named **Untitled1**, and contains four default processes. See [Understanding Processes on page 72](#). This default protocol file can be used to create a new protocol file, closed, or ignored to work with an existing protocol file or data file.

Protocol (*.fmp) and data (*.fmd) files both show the process steps they use at the top of the window, ordered from left to right. Clicking on the process icons opens the related settings dialog. A data file is a protocol file that also includes acquired data in the Analysis process dialog.

Experiment Summary

The Experiment Summary is a single, text-based protocol summary screen.



To view the Experiment Summary page:

- Select **View > Experiment Summary**.

Instrument Status-Configuration-Process Panel

By default the **Instrument Status-Configuration-Process** panel is open and docked to the bottom of the main screen. See [User Interface Overview on page 48](#).

The following tabs are in the panel:

- [Instrument Status Tab on page 67](#)
- [Instrument Configuration Tab on page 69](#)
- [Process Tab on page 70](#)

To undock and relocate the panel:

1. Click and drag on the title bar on the left of the panel.
2. Drag the panel to where you want it.

To hide the panel, do one of the following:

- Click the **x** in the title bar of the dialog.
- Select the **View > Instrument Status** menu option.

To show the panel when hidden, select the **View > Instrument Status** menu option.

Instrument Status Tab

The current status of the system hardware components report in the **Instrument Status** tab panel.

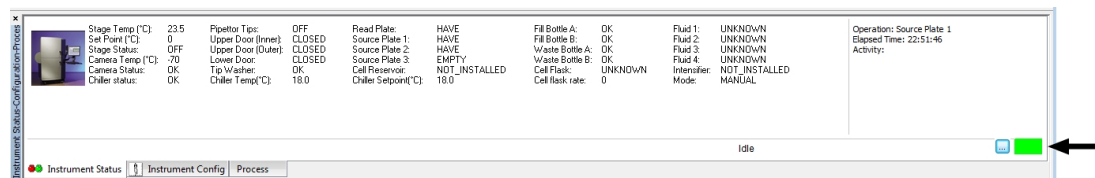


Figure 3-1: Example Instrument Status Tab Panel With Online Status Indicator

Table 3-11: Instrument Status Options

| Options | Description |
|--------------------|--|
| Stage Temp (°C) | Displays what the stage temperature is currently. |
| Set Point | Displays the stage temperature setting. |
| Stage Status | Displays whether the heated stage is turned on or off. |
| Camera Temp (°C) | Displays the camera temperature. Operating temperature for the camera depends on the installed camera. EMCCD camera: -70°C (-94°F) ± 2°C HS EMCCD camera: -70°C (-94°F) ± 2°C ICCD camera: -20°C (-4°F) ± 5°C |
| Camera Status | Indicates whether the camera is turned on or off. |
| Chiller status | Indicates whether the chiller is available or not. |
| Pipettor Tips | Reports when tips are loaded on the pipettor head. |
| Upper Door (Inner) | Reports whether the inner-upper door (observation panel) is open or closed. Note: The system runs as long as the inner door is closed, however data might not be valid if the outer door is open. |
| Upper Door (Outer) | Reports whether the outer-upper door is open or closed. |
| Lower Door | Reports whether the lower door is open or closed. |
| Tip Washer | Indicates the status of the tip washer. |
| Chiller Temp (°C) | Reports the current temperature in the chiller. |
| Read Plate | Reports when a plate is present in Position 3 (Read Plate position). |
| Source Plate 1 | Reports when a plate is present in Position 1 (Source Plate 1 or Tip Loading position). |
| Source Plate 2 | Reports when a plate is present in Position 2 (Source Plate 2). |

Table 3-11: Instrument Status Options (continued)

| Options | Description |
|-----------------------|--|
| Source Plate 3 | Reports when a plate is present in Position 4 (Source Plate 3 or Cell Reservoir). |
| Cell Reservoir | Reports if Installed or Not_Installed . |
| Chiller Setpoint (°C) | Reports the set point from the chiller. |
| Fill Bottle A | Reports when bottle A is empty of wash solution. |
| Fill Bottle B | Reports when bottle B is empty of wash solution. |
| Waste Bottle A | Reports when waste bottle A is full. |
| Waste Bottle B | Reports when waste bottle B is full. |
| Cell Flask | Reports the last known state of the stir Cell Flask. At start the state is <i>Unknown</i> until the Cell Flask is used. |
| Cell flask rate | Reports the set stir rate of the cell flask. If the Cell Suspension option is installed and the stir rate is 0, an exclamation sign displays. |
| Fluid 1 | Reports the last known state of the Fluid 1 bottle. At the start the status is <i>Unknown</i> until that Fluid is used. |
| Fluid 2 | Reports the last known state of the Fluid 2 bottle. At the start the status is <i>Unknown</i> until that Fluid is used. |
| Fluid 3 | Reports the last known state of the Fluid 3 bottle. At the start the status is <i>Unknown</i> until that Fluid is used. |
| Fluid 4 | Reports the last known state of the Fluid 4 bottle. At the start the status is <i>Unknown</i> until that Fluid is used. |
| Intensifier | Reports if Installed or Not_Installed . |
| Mode | Reports if ScreenWorks Software is in Manual or Remote mode. |

Identify Connection Status

To view the connection status:

- In the bottom-right corner of the panel, identify the color-coded connection status icon, which reports the connection status between ScreenWorks Software and the instrument. Green means connected, and black means disconnected. For more details, see [Instrument Status Colors on page 209](#).

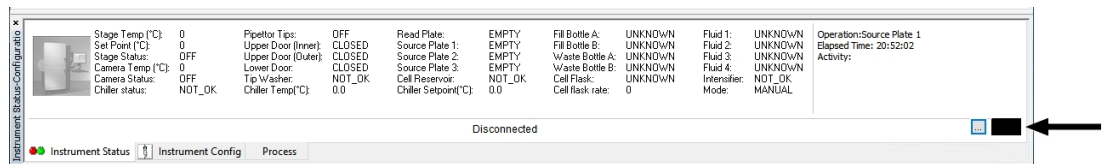



Figure 3-2: Example Instrument Status Tab Panel With Offline Status Indicator

View Messages and Faults

To view status messages and faults:

- Next to the color-coded connection status icon, click  **Instrument Status** to see a list of the last thousand messages in the Instrument Status History dialog.

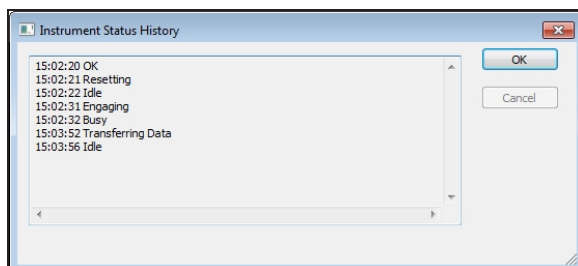


Figure 3-3: Example Instrument Status History Dialog



Tip: If needed, the dialog text can be copied to the clipboard to paste elsewhere.

Instrument Configuration Tab

The **Instrument Configuration** tab indicates the current instrument configuration of the LED banks, emission filters, pipettor head, and optional FLIPR Cycler.

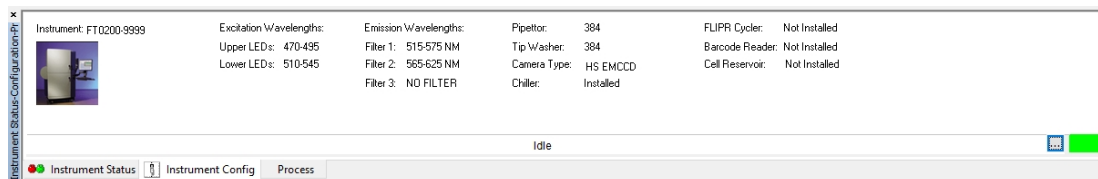


Figure 3-4: Example Online Instrument Config Tab Panel

If the system is Offline, you can configure these settings to define protocols. For camera types details, see the *CCD Camera Options* section of the *FLIPR Penta High-Throughput Cellular Screening System User Guide*. See [Online Mode vs. Offline Mode on page 47](#).



CAUTION! If a protocol created offline does not match the Instrument Configuration when opened online, the protocol does not run until the configuration of the protocol and instrument match.

Table 3-12: Instrument Configuration Options

| Options | Description |
|------------------------|--|
| Excitation Wavelengths | Displays the excitation wavelengths installed on the system. Upper LEDs —Displays the wavelength range of the top set of LED banks in the LED modules. Lower LEDs —Displays the wavelength range of the lower set of LED banks in the LED modules. |

Table 3-12: Instrument Configuration Options (continued)

| Options | Description |
|----------------------|---|
| Emission Wavelengths | Displays the emission filter wavelengths installed on the system. Up to three filters can be installed at the same time. |
| Pipettor | Displays the type of pipettor head (96, 384, 1536, 384 pin tool, or 1536 pin tool) installed on the system. Note: The pipettor head format must agree with the tip wash reservoir format. A warning is issued if these are different. |
| Tip Washer | Displays the type of tip washer (96, 384 or 1536) installed on the system. |
| Camera Type | Select from EMCCD, HS EMCCD, or ICCD camera. |
| Chiller | Reports when a chiller is installed. |
| FLIPR Cycler | Reports when the FLIPR Cycler is installed. |
| Barcode Reader | Reports when a barcode reader is installed. |
| Cell Reservoir | Use the check box to indicate whether or not the Cell Reservoir is installed. |

Process Tab

The Process tab is used to create new protocols and edit existing protocols.



Figure 3-5: Example Online Process Tab Panel Docked

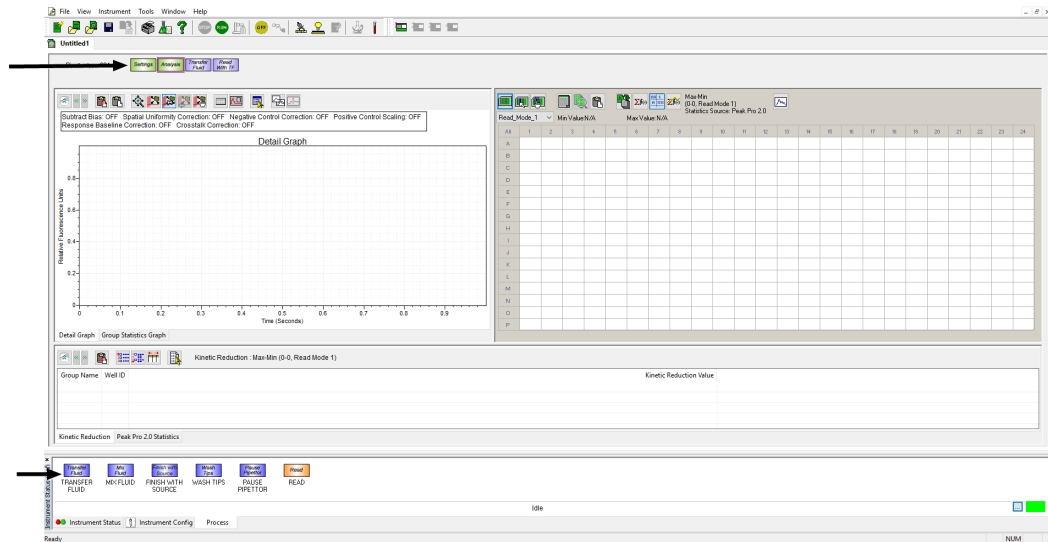
The Process tab panel displays processes that can be incorporated into your experiment protocols. See [Using Protocols on page 71](#). The color of the process icon depends on the function in the protocol. See [Protocol Process Icon Colors on page 71](#).



Note: The available processes vary according to the instrument configuration. For example, different process opts appear when the instrument is configured to use a pin tool instead of a pipettor.

For details about each process, see [Understanding Processes on page 72](#).

Protocol Process Icon Colors



Green—indicates that the process is required to run an experiment and cannot be deleted.

Blue—indicates that the process involves liquid handling. These steps are always run series with each other, and never run simultaneously.

Orange—indicates that the process runs simultaneously with any liquid handling process, if possible.

Purple—indicates linked processes, such as **Mix with TF** and **Read with TF**, which are connected to the **Transfer Fluid** process.

Status Bar

The status bar, across the bottom of the main window, provides tool tips for commands in the main menu. When you open a menu from the menu bar and place the cursor over a command, a description of the command is displayed in the status bar.

The status bar can be hidden or shown from the **View > Status Bar** toggle command.

Using Protocols

The protocols you create determine how the system functions while running an experiment. Protocols consist of a series of process configurations. As you create your experiment protocol, you configure each process. See [Creating New Protocols on page 124](#). Process options are located in the Process tab in the Instrument Status-Configuration-Process panel. See [Process Tab on page 70](#).

*** Tip:** Refer to the *FLIPR Penta High-Throughput Cellular Screening System Protocol Guide* for specific assay protocols.

To save a data file as a protocol file, see [Saving Data Files as Protocol Files on page 128](#).

Understanding Processes

The following provides details for using the available Process options used to create your experiment protocols.

Settings Process

Settings is the first process in all protocols. Use it to define read modes, plate positions, and data directories. You can also configure automatic file naming, and set plate temperature.

Select the green **Settings** icon in the Experiment window to view the **Settings** page where the settings are made.

Pipettor type:384

Setup Read Mode:

| Read Mode Name | Exc. Wlength(nm) | Em. Wlength(nm) | Gain | Exp. Time(s) | Exc. Intensity(%) | Gate Open(%) |
|---|------------------|-----------------|------|--------------|-------------------|--------------|
| <input checked="" type="checkbox"/> Read_Mode_1 | 470_495 | 515_575 | 1.00 | 0.0500 | 50 | N/A |
| <input type="checkbox"/> Read_Mode_2 | 470_495 | 515_575 | 1.00 | 0.0500 | 50 | N/A |
| <input type="checkbox"/> Read_Mode_3 | 470_495 | 515_575 | 1.00 | 0.0500 | 50 | N/A |
| <input type="checkbox"/> Read_Mode_4 | 470_495 | 515_575 | 1.00 | 0.0500 | 50 | N/A |

Assign Plate to Position:

| Position | Plate Name | Barcode Expect | Bar Code List | Flask speed |
|----------------|------------|----------------|---------------|-------------|
| Read Plate | default384 | NO | | N. |
| Source Plate 1 | default384 | NO | | N. |
| Source Plate 2 | N0NF | NO | | N. |

Load Tips Position

Data File Name:

Include Date User defined name:

Include Bar Code Example: MMDDYYYY_n000.fmd

Temperature Control (C):

Use Manual

Folders:

Data folder:

Auto Export folder:

Setup Read Mode

The **Setup Read Mode** table at the top of the **Settings** page displays settings for up to four read modes that the FLIPR Penta System can run in a single experiment.

Setup Read Mode:

| Read Mode Name | Exc. Wlength(nm) | Em. Wlength(nm) | Gain | Exp. Time(s) | Exc. Intensity(%) | Gate Open(%) |
|---|------------------|-----------------|------|--------------|-------------------|--------------|
| <input checked="" type="checkbox"/> Read_Mode_1 | 470_495 | 515_575 | 1.00 | 0.0500 | 50 | N/A |
| <input type="checkbox"/> Read_Mode_2 | 470_495 | 515_575 | 1.00 | 0.0500 | 50 | N/A |
| <input type="checkbox"/> Read_Mode_3 | 470_495 | 515_575 | 1.00 | 0.0500 | 50 | N/A |
| <input type="checkbox"/> Read_Mode_4 | 470_495 | 515_575 | 1.00 | 0.0500 | 50 | N/A |

Each read mode defines:

- Reading Mode
- Excitation and Emission Wavelength
- Excitation Intensity
- Camera Mode (EMCCD camera and HS EMCCD camera)
- Exposure Time
- Camera Gain
- Gate Open (ICCD camera only)
- Auto Stabilize (HS EMCCD camera only)

For protocols that use multiple read modes, the instrument alternates between the selected read modes for each time point, outputting distinct sets of readings, one for each mode. ScreenWorks Software automatically calculates the ratio of the readings from each mode for each time point.

Protocol files that use read modes not supported by your instrument configuration display the invalid read mode rows in red.

To enable a read mode:

- In the **Setup Read Mode** table, select a **Read Mode Name** check box.

To disable a read mode:

- In the **Setup Read Mode** table, deselect a **Read Mode Name** check box.

Edit Read Mode

To edit read mode settings:

1. Select a **Read Mode Name** row and click **Edit Mode**.
Alternatively, double-click a **Read Mode Name** row.
2. In the **Edit read mode** dialog, change the following settings options as needed, then click **OK** to finish.



Note: Available options depend on the camera installed in your instrument.

Table 3-13: Edit Read Mode Dialog Options

| Options | Description |
|--------------------------------|--|
| Read Mode Name | Enter a name for the selected read mode. |
| Reading Mode | Select Fluorescence or Luminescence . |
| Excitation/Emission Wavelength | Select a needed excitation LED and emission filter wavelength pair from the list. NONE instructs the instrument to not turn on any LEDs or use an empty emission filter position. These are typically used in luminescence experiments. Note: Only calibrated wavelength pairs are available when the software is in Online mode. See Calibrating the Optics on page 132 . |
| Excitation Intensity | Select a value from the list to regulate the intensity of light emitted by the LEDs for a given Fluorescence read. The range is 20% to 100%. |
| Camera Mode | Select HighSpeed , Sensitivity , or Normal to specify your camera performance preference. Note: This option is only available with the HS EMCCD camera. See Understanding HS EMCCD Camera Protocol Saturation on page 38 . |
| Exposure Time | Enter a time in seconds to regulate the time that light is collected and measured by the camera. A longer exposure increases signal intensity. Exposure time affects the read interval for data collection. For example, in Camera Mode > Normal, an exposure time of 0.4 seconds added to a required 0.1 seconds camera integration time makes the highest update frequency 0.5 seconds. Depending on the camera, the range is 0.01 seconds to 30 seconds. For example, using the HS EMCCD camera in Camera Mode > HighSpeed, the range is 0.001 seconds to 30 seconds. |
| Camera Gain | Select a value from the list to regulate the amplification of the camera power. Increasing the camera gain increases the signal. Ranges depend on the camera in use: EMCCD camera: 1–240 HS EMCCD camera: Reading Mode >Fluorescence : Camera Mode > Normal: 1-128 or Camera Mode > HighSpeed or Sensitivity: 20-128. Luminescence mode: 1–500 ICCD camera: Fluorescence read mode: preset to 2000. Luminescence mode: 2000–280,000 |
| Auto Stabilize | Automatically adjusts the camera for 50 seconds for optimal functionality before the protocol starts. Note: This option is only available with the HS EMCCD camera. See Stabilizing the HS EMCCD Camera on page 38 . |
| Gate Open | Select the percentage of each frame captured to further regulate the signal intensity. Note: This option is only available with the ICCD camera in Fluorescence mode. |



Note: Differences in **Camera Gain** or **Exposure Time** between read modes significantly increase the minimum update time. For fastest update time, use the same gain and exposure time for all read modes.

Assign Plate to Position

Current plate assignments are reported in the **Assign Plate to Position** table. The pipettor head must have a plate type assignment for each of the positions to be used in your experiment.

Assign Plate to Position:

| Position | Plate Name | Barcode Expect | Bar Code List | Flask speed |
|----------------|------------|----------------|---------------|-------------|
| Read Plate | default384 | NO | | N/A |
| Source Plate 1 | default384 | NO | | N/A |
| Source Plate 2 | NONE | NO | | N/A |

Load Tips Position Edit Plate...

For each plate position, the table lists:

- **Plate Name**—the type of plate, as configured in **Tools > Plate Library**.
- **Bar Code Expect**—Whether or not you want the system to read a bar code from the plate. Bar code numbers can optionally be incorporated into output data file names.
- **Bar Code List**—in data files only, the bar codes of plates used in the experiment, in the given plate position.

To change these settings, click **Edit Plate**.

If Position 1 is to be used only for tips, and not to hold a source plate, select the **Load Tips Position** check box. This removes Source Plate 1 from the assignment table.

If the Cell Reservoir is installed in the system, Source 3 is automatically dedicated to the head format, such as **defaultcellres384**.

Edit Plate

To change a plate position assignment:

1. Select a **Position** row and click **Edit Plate**.
Alternatively, double-click a **Position** row.
2. In the **Change Plate** dialog, change the following settings options as needed, then click **OK** to finish.

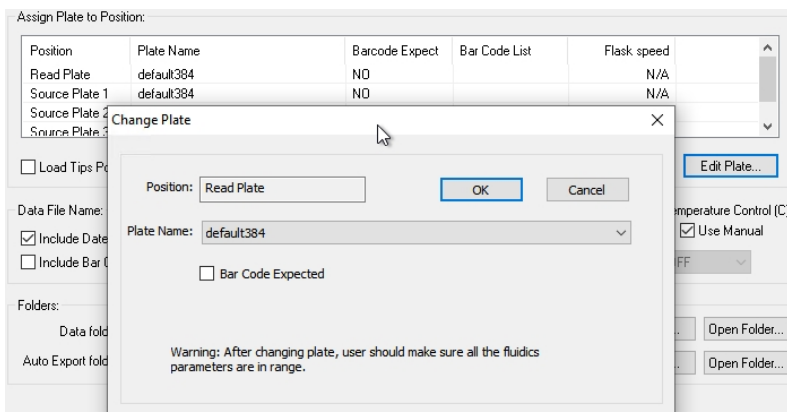


Table 3-14: Change Plate Dialog Options

| Options | Descriptions |
|--------------------------|---|
| Position | Displays the plate position you are editing. |
| Plate Name | Select the plate type from the drop-down list. The list contains default plates and any additional plates that have been added using Tools > Plate Library . Note: Only plates with the same number of wells or one order of complexity higher than the pipettor head format are displayed. |
| Bar Code Expected | Check this option if using a bar code for the plate. The bar code number can be incorporated into the file names of data files generated by the protocol. In Manual Mode , you are prompted to enter the bar codes prior to starting an experiment. In Remote Mode the bar code number is automatically scanned on the FLIPR Cyclor landing pad or is passed to ScreenWorks by the third-party plate handler. |
| Cell Flask Spinning Rate | When the Cell Reservoir from the Cell Suspension option is installed, spinning rate can be set here or Use Manual can be selected and the manual settings apply. |



CAUTION! Each plate position has a mechanical sensor to detect the presence of a plate, tip rack or boat in each position. However, the instrument does not distinguish the type of apparatus placed in a position.

Data File Name

Use the **Data File Name** options to configure a file-naming protocol for data files created in experiments. The file-naming protocol applies to ***.fmd** (data) and ***.png** (image) files.

Data File Name:

Include Date User defined name:

Include Bar Code Example: MMDDYYYY_rXXX.fmd

Table 3-15: Date File Name Options

| Options | Description |
|--------------|---|
| Include Date | Check to include the date in the file name. It is checked by default. Note: To ensure that data files are not overwritten, we recommend that this option remains checked. |

Table 3-15: Date File Name Options (continued)

| Options | Description |
|-------------------|--|
| Include Bar Code | When selected, the bar codes of the first five plates used in the experiment are included in the file names. If there are fewer than five bar codes, then only those available are used. Note: Bar Code Expected must be selected for the desired plate position in order for the bar code to be incorporated. |
| User Defined Name | Enter a string of characters (up to 25) to be added to the file name. |

The base structure for file names is as follows:

Date_UserDefinedName_Barcode1_Barcode2_Barcode3_Barcode4_Barcode5_NNN.fmd

where NNN is an integer from 1 to 999. The integer value NNN starts at 001 and increments for each data file generated with the same base data file name.

The date uses format: MMDDYYYY, where M = month, D = day, and Y = year.

If a software crash or other error occurs during a run that interferes with normal instrument operation, then a data file is automatically stored as:

InterruptedExperiment_Date_Time_NNN.fmd

Folders

Use the **Folders** fields to specify into which directories your data and export files save.

- **Change**—allows you to select a different directory.
- **Open Folder**—opens the Windows Explorer dialog directly to the directory designated.

By default, data, image and export files save in the following location:

C:\Documents and Settings\[your user name]\My Documents\Molecular Devices\ScreenWorks\MyData

Temperature Control

The **Temperature Control** field allows you to set the plate-stage heaters to a needed temperature for an experiment. Temperature range is from ambient 25°C (77°F) to 40°C (104°F).

Select **Use Manual** to control the heaters manually from ScreenWorks Software at the time the experiment is run.

If at the start of an experiment, the stage temperature is different from the selected value, a message displays. You can wait until the stage reaches the needed temperature or run the experiment anyway.



Note: When the temperature is manually turned on through the **Instrument** menu, a setting of “0” in the protocol does not affect the temperature setting. To turn off the temperature, manually select **Instrument > Manual Operations > Temperature Control Off**.

Analysis Process

The **Analysis** process is included in all protocols and is always the second process after **Settings**.

- [Viewing Data](#)
- [Grouping, see page 86](#)
- [Corrections, see page 91](#)
- [Exporting Data, see page 157](#)

Viewing Data

Acquired data can be viewed in ScreenWorks Software on the main **Analysis** process page, in the **Multi-Well Graph**, or in larger **Detail Graph**. The options are described in this section.

Multi-Well Graph

The major part of the **Analysis** process page consists of a grid representing the wells of the read plate used in the protocol or data file (for example, with 96-wells, 384-wells, or 1536-wells). This is called the **Multi-Well Graph**. For protocol files this graph is empty, as no data have been acquired, but in data files each cell of the graph has a trace of one of the measurements taken from the corresponding read plate well. The maximum and minimum relative light units values of the displayed data are reported above the Multi-Well Graph.

The Multi-Well Graph displays, for each well represented, one trace of relative light units versus time. If two or more read modes were recorded, select the mode to view from the display list box above the graph. To see the ratio of two read modes, enable **Ratiometric Options** in the **Configure Corrections > Corrections** dialog, and select **Ratio** in the display list.

The **Read Mode** defined in the Multi-Well Graph is also applied to any traces displayed in the **Detail Graph**.

The traces displayed in the Multi-Well Graph can be adjusted by the application of different data-correction options available in the **Correction** dialog, for example, the traces can be scaled relative to the average positive control response. See [Corrections on page 91](#).

If groups have been defined for the assay, wells in the Multi-Well Graph are color-coded to represent these. See [Grouping on page 86](#). In addition, there is the option to include the kinetic reduction value below each well trace in the Multi-Well Graph.

Buttons, just above the Multi-Well Graph, open dialogs that offer various data analysis options; some relevant to protocol set up, some to already-acquired data in data files, and some to both. Multi-Well Graph buttons and descriptions include:

Table 3-16: Multi-Well Graph Button Options













| Buttons | Name | Description |
|---|----------------------|---|
|  | Configure Groups | Opens the Grouping dialog to classify wells in the plate into groups (for example, positive and negative controls). Groups created in protocol files are transferred to all the data files generated by the protocol. In data files, new groups can be created or existing groups edited. See Grouping on page 86 . |
|  | Select Groups | Opens the Select Groups dialog to classify which groups are quickly displayed in the Detail Graph from a predefined list based on the groups defined in the Grouping dialog. Note: Hold down the SHIFT or CTRL key to select multiple groups. |
|  | Group Selection Mode | Button toggles between Group Selection Mode and Well Selection Mode . During Group Selection Mode , wells selected within a predefined group displayed in the Multi-Well Graph cause all wells within those groups to be displayed in the Detail Graph . |
|  | Well Selection Mode | Button toggles between Group Selection Mode and Well Selection Mode . Well Selection Mode only displays wells selected in the Multi-Well Graph on the Detail Graph with their respective group colors assigned. |
|  | Notes | In protocol files, use this dialog to write comments to be stored with all data files generated by the protocol. In the data files, view the comments (now read-only) in the same dialog. |
|  | Images | For data files only, where images have been saved during an experiment (by checking Save Images in a Read with TF step) these can be viewed by clicking the Images button. |
|  | Copy | Copies data in the Multi-Well graph to the clipboard so it can be pasted in a different program such as Microsoft Word. |

Table 3-16: Multi-Well Graph Button Options (continued)

| Buttons | Name | Description |
|---|------------------------------|---|
|  | Configure Auto-Export | <p>Available for protocol files, this button opens the Auto-Export dialog allowing configuration of statistics and time sequence files to be automatically created whenever the protocol is run. This button also works to export data after a data file has been created.</p> <p>For data files the button changes to Export, and opens the File Export dialog, also opened from the File > Export command. See Exporting Data on page 157.</p> |
|  | Configure Correction | <p>Opens the Corrections dialog to apply various corrections to modify data display. There is also the option to view ratiometric data. Settings made here in protocol files affect how data are viewed when the protocol is run, but since all raw data are stored in data files, these options all remain available for acquired data as well. See Corrections on page 91.</p> |
|  | Hide Kinetic Reduction Value | Hides the kinetic reduction values that are displayed in the Multi-Well Graph. |
|  | Show Kinetic Reduction Value | Shows the kinetic reduction value in each well of the Multi-Well Graph. |
|  | Configure Kinetic Reduction | <p>Opens the Kinetic Reduction Configuration dialog to define the parameters used to define the kinetic reduction.</p> <p>Reduction Type—Defines the reduction to be applied to the kinetic data traces displayed in the Multi-Well Graph. See Kinetic Reduction Types on page 81.</p> <p>Start Read—Define the first read to be used to determine the kinetic reduction.</p> <p>End Read—Define the last read used to determine the kinetic reduction.</p> <p>Read Mode—Select the read mode to apply the kinetic reduction to.</p> |

Kinetic Reduction Types

The **Reduction Type** field is located in the **Kinetic Reduction Configuration** dialog. Available options depend on the software version in use. A separate license for ScreenWorks® Peak Pro™ Software Software Version 1.0 adds advanced peak detection and characterization measurements to the field of standard measurement options (see [Figure 3-6](#)). This functionality is optimal for relatively fast events that have been acquired with long sampling intervals. Contact your Molecular Devices Sales Representative for details.



CAUTION! High speed camera data is not compatible with Peak Pro 1.0. You must use ScreenWorks Peak Pro Software Software Version 2.0 for the high speed camera data analysis. See [Using ScreenWorks Peak Pro Software Version 2.0 on page 165](#).

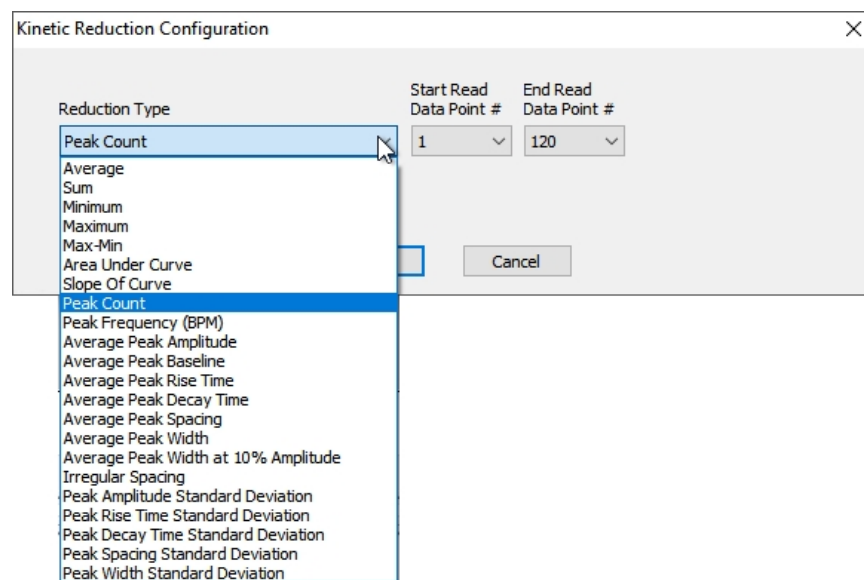


Figure 3-6: Kinetic Reduction Types Including Additional ScreenWorks Peak Pro 1.0 Reductions

Standard measurement options and definitions include:

- **Average**—Numerical average of relative light units counts of the selected reads.
- **Sum**—Numerical sum of relative light units counts of the selected reads.
- **Minimum**—Lowest detected count (a single number) of all the selected reads.
- **Maximum**—Highest detected count (a single number) of all the selected reads.
- **Maximum-Minimum**—Result of subtracting the minimum count (a single number) from the maximum count (a single number).
- **Area Under Curve**—Numerical calculation of the area under the curve for the selected reads.
- **Slope of Curve**—Result of calculating the slope between two selected reads.

(With optional license) ScreenWorks Peak Pro Software Software Version 1.0 measurement options and definitions include:

- **Peak Count**—Number of peaks or action potentials detected within the read data that meet the specifications in the Configure Peak Detection dialog settings.
- **Peak Frequency (BPM)**—The Beats-Per-Minute frequency of the peaks detected based on the inverse of the peak temporal spacing. At least two peaks are required for this measurement.
- **Average Peak Width**—Average peak width measurement in seconds from the 50% peak amplitude from baseline.
- **Peak Width Standard Deviation**—Standard deviation of average peak widths.
- **Average Peak Amplitude**—The average peak amplitude relative to the Average Peak Baseline measurement.
- **Peak Amplitude Standard Deviation**—Standard deviation of peak amplitudes.
- **Average Peak Baseline**—Average baseline amplitude measurement from the base of each peak detected.
- **Average Peak Spacing**—Average peak spacing in seconds. At least two peaks are required for this measurement.
- **Peak Spacing Standard Deviation**—Standard deviation of peak spacing.
- **Average Peak Rise Time**—Average time measured on the rising edge for each peak. Equal to the time between the 10% and 90% peak amplitudes.
- **Rise Time Standard Deviation**—Standard deviation of rise times.
- **Average Peak Decay Time**—Average time measured on the falling edge for each peak. Equal to the time between the 90% and 10% peak amplitudes.
- **Peak Decay Time Standard Deviation**—Standard deviation of decay times.
- **Average Peak Width at 10% Amplitude**—Average peak width in seconds at the 10% peak amplitude from baseline.
- **Irregular Spacing**—Identifies wells that have irregular spacing between peaks. Wells with regular peak temporal spacings are marked as **OK**. Wells with missing peaks are marked as **MISS**. Wells with extra peaks are marked as **EXTRA**. Wells with peak spacing characteristics of both missing and extra are marked as **IRREG**.

The **Configure Peak Detection** button activates when a Peak Pro measurement is selected from the **Reduction Type** field (see [Figure 3-7](#)). The **Configure Peak Detection** dialog allows the Peak Pro measurement settings to be adjusted and optimized (see [Figure 3-8](#)).

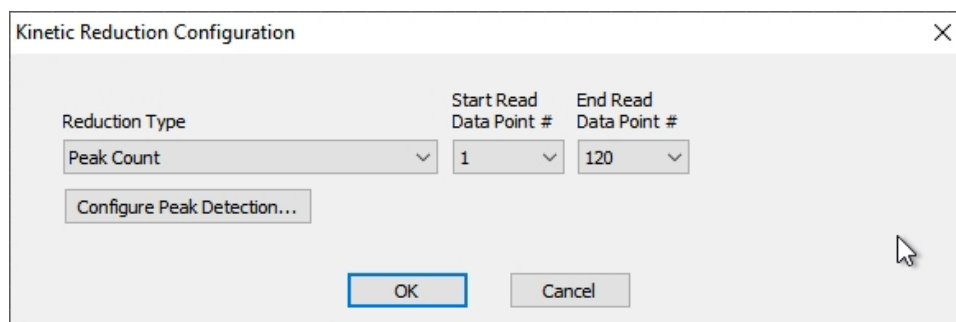
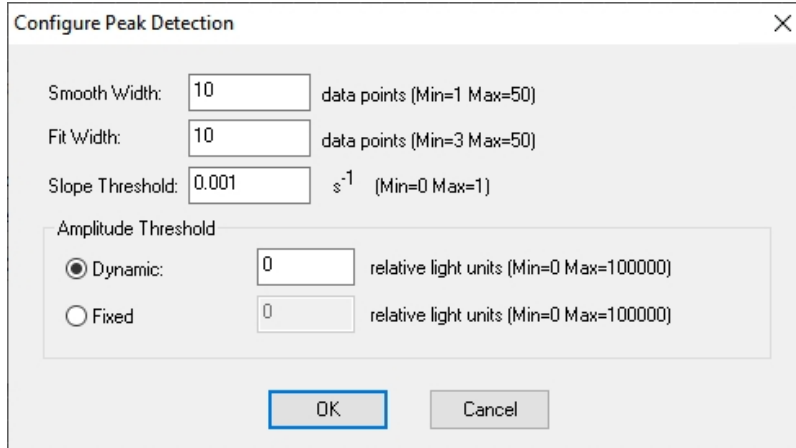


Figure 3-7: Configure Peak Detection Button Activation

 **Tip:** Start with default settings.

Configure Peak Detection dialog options include:

- **Smooth Width:** Running average to reduce noise, set between 5 to 10.
- **Fit Width:** Number of data points that are fit for each peak, set between 5 to 15.
- **Slope Threshold:** Distinguishes between cardio beats and noise, set to 0.001.
- **Amplitude Threshold Dynamic:** Value setting is relative to the data average.
- **Amplitude Threshold Fixed:** Value setting is relative to zero.



Configure Peak Detection

Smooth Width: data points (Min=1 Max=50)

Fit Width: data points (Min=3 Max=50)

Slope Threshold: s⁻¹ (Min=0 Max=1)

Amplitude Threshold

Dynamic: relative light units (Min=0 Max=100000)

Fixed: relative light units (Min=0 Max=100000)

OK Cancel

Figure 3-8: Configure Peak Detection Dialog

Detail Graph

All the read modes for wells selected in the multi-well graph can be displayed enlarged in a detail graph. You can select individual or multiple wells to view, or select by group.



Note: When the FLIPR Penta System is configured for two or more read modes, it cycles through each mode alternating between them for each reading. Although readings from each mode within a cycle occur at different times, they are represented as occurring simultaneously.

The data displayed in the **Detail Graph** have the same corrections applied as the data displayed in the Multi-Well Graph. The **Detail Graph** reports these settings at the top of the graph window:

- Subtract Bias
- Spatial Uniformity Correction
- Negative Control Correction
- Positive Control Scaling
- Response Over Baseline
- Crosstalk Correction

These correspond directly to options in the **Correction** dialog, opened from the **Analysis** process page. To change these settings, change them in the **Correction** dialog. This changes all data in the **Multi-Well Graph** and **Detail Graph** to reflect the new correction applied.

Detail Graph allows a number of data-selection options:

Table 3-17: Detail Graph Button Options













| Buttons | Name | Description |
|---|--------------------|--|
|  | Copy Graph | Copies Graph in the Detail Graph to the clipboard so it can be pasted in a different program such as Microsoft Word. |
|  | Copy Graph Data | Copies graph data in the Detail Graph to the clipboard so it can be pasted in a different program such as Microsoft Excel. |
|  | Zoom Mode | Click to enable zoom capability in the graph. Once the Zoom button is pressed, drag over a region in the graph to view just that area enlarged to the full size of the graph. Alternatively, drag the cursor along a part of the x or y axis to zoom in on just that axis. |
|  | Auto Scale | Manual prompt to automatically scale the Detail Graph and Multi-Well Graphs to include all data points of the desired traces. |
|  | Auto Scale Always | Automatically scales the Detail Graph and Multi-Well Graph to include all data points of the desired traces without manual prompting. |
|  | Undo Zoom | Rescales a graph to the original settings. |
|  | Manual Scale Graph | Opens the Manual Scale Graph dialog to set the maximum and minimum values for the X- and Y-axis of the Detail Graph and Multi-Well Graph. |
|  | Show/Hide Data | Toggle button that shows or hides the relative light units values for the selected traces at each read. |
|  | Show Point Labels | Show Point Labels writes the value of each data point (for example, relative light units value) beside the point on the graph. |
|  | Show/Hide Legends | A toggle button that displays or hides the Detail Graph legend. |

Table 3-17: Detail Graph Button Options (continued)

| Buttons | Name | Description |
|---|------------------------------|--|
|  | Average/Overlay Trace | <p>A toggle button that displays the average or overlay traces for selected groups.</p> <p>Average Group—Displays average time point values for all wells in the selected group. A separate trace is generated for each read mode.</p> <p>Note: When Well Selection Mode selected in the Multi-Well Graph, all traces displayed in the Detail Graph are averaged regardless of the group they are assigned.</p> <p>Overlay Group—Displays traces for all the wells in selected groups.</p> <p>Note: Traces are color-coded by group.</p> |
|  | Show/Hide Standard Deviation | <p>A toggle button that enables or hides the standard deviation for the average group trace.</p> <p>Note: Average Group must be enabled to access this function.</p> <p>When Well Selection Mode is selected in the Multi-Well Graph, all traces displayed in the Detail Graph are averaged regardless of the group they are assigned, and the standard deviation for this average is displayed.</p> |

Individual Well Selections

To view a single well in the **Detail Graph**, double-click or drag the mouse cursor inside the Multi-Well Graph over a well of interest.

To view the traces of multiple wells overlaid in the **Detail Graph**, drag the cursor inside the Multi-Well Graph over the multiple wells of interest. Data from any rectangular block of wells within the plate can be viewed in this way. Additional wells can be added to the overlay by double-clicking or dragging the mouse cursor over the wells of interest.

Group Selections

You can also select data to view in the **Detail Graph** using group membership (see [Grouping on page 86](#)). Use **Select Group** or **Group Selection Mode** functions found in the Multi-Well Graph.

Analyzing Data

The **Grouping** and **Corrections** dialogs are used to analyze data produced during an experiment. **Settings** made in these dialogs in a protocol file affect the way that data, in data files generated with the protocol, are displayed when first viewed in ScreenWorks Software. They also affect data in the automatic output option **Auto Export** that occur when the protocol is run. See [Exporting Data on page 157](#). Data files created by ScreenWorks Software retain all raw data readings, so the analysis options in the **Grouping** and **Correction** dialogs apply to acquired data, modifying the display of these, irrespective of the analysis settings configured in the original protocol. See [Grouping on page 86](#) and [Corrections on page 91](#).

Grouping

Groups are used to select data to be viewed in the detail graph, and the data itself can be corrected relative to responses to positive and/or negative control groups, in the **Correction** dialog.



Define groups in the **Analysis** > **Configure Groups** > **Grouping** dialog.

A group is a selection of wells with a common characteristic—usually in the type of compound added to the wells. Three groups—background fluorescence correction, positive controls and negative controls—are defined by default, but no wells are assigned to them. You specify the well assignment for these groups.

You can create additional groups to represent wells with a specific ligand, or ligand concentration, or any other characteristic you want to classify data by. Group definition includes specific input for compound concentration, as this is commonly used for group definition. Additionally, the **Series** option allows you to create a series of groups, where a number of different concentrations are automatically assigned to each group. This can be useful for creating groups for an IC_{50} or EC_{50} experiment.

The **Grouping** dialog contains a grid representing the 96-well, 384-well or 1536-well read plate, and above this a list box where groups are defined and selected. The basic mode of operation is to select a group from the list box, then drag the cursor over the wells in the grid that belong to that group. Additional groups can be defined to add to the list.

Table 3-18: Grouping Dialog Options

| Options | Description |
|----------------------|---|
| Groups | <p>Displays the list of defined groups, including group name (typically compound name), concentration and notes associated with the group. Wells assigned to a group in the grid below have the same color as the group in the list.</p> <p>Positive Controls—Use this group to assign positive control wells, typically the maximum response to a concentration of an agonist. Note: Positive Control wells need to be defined in order to use the Positive Control Scaling Correction.</p> <p>Negative Controls—Use this group to assign negative control wells, typically buffer addition controls. Note: Negative control wells need to be defined in order to use the Negative Control Correction feature.</p> <p>BF Controls—Use this group to assign wells in which to measure background fluorescence. Measurements for each time point from these wells are used for Background Fluorescence Correction for ratiometric data. See Ratiometric Options on page 93.</p> <p>_add new group—Use this to define a new individual group or a series of groups that are user-defined.</p> |
| Add/Edit Group | <p>Opens the Edit Group dialog to edit the group selected in the Groups list. This could be an existing group, or create a new group by selecting the Add New Group row. See Adding a New Group on page 88.</p> <p>Shortcut: Open the Edit Group dialog by double-clicking on a group in the Groups list.</p> |
| Delete Group | <p>Deletes the group selected in the Groups list. Any wells assigned to the group lose their coloring, indicating that they no longer belong to any group.</p> |
| Delete All Groups | <p>Deletes all user-defined groups. Note: Positive, negative and background fluorescence control groups cannot be deleted, however this command deletes all wells assigned to these groups.</p> |
| Clear All Selections | <p>Removes all group assignments from all wells. Group names in the top list are not affected.</p> |
| Undo Last Selection | <p>Removes group membership from the last wells that were assigned to a group.</p> |

Assigning Wells to Groups

To assign wells to a group:

1. Select the appropriate group from the **Groups** list, for example, **Positive Controls**.
2. Drag the cursor over the wells from the plate layout that you want to include in the selected group. Alternatively, the row or column title can be selected to include all the respective wells in the group.



Note: Wells in a group do not need to be contiguous. Multiple areas of the plate can be selected.

You can deselect individual or neighboring wells by right clicking in one corner of a well to be deselected, then dragging over the well or wells to deselect.

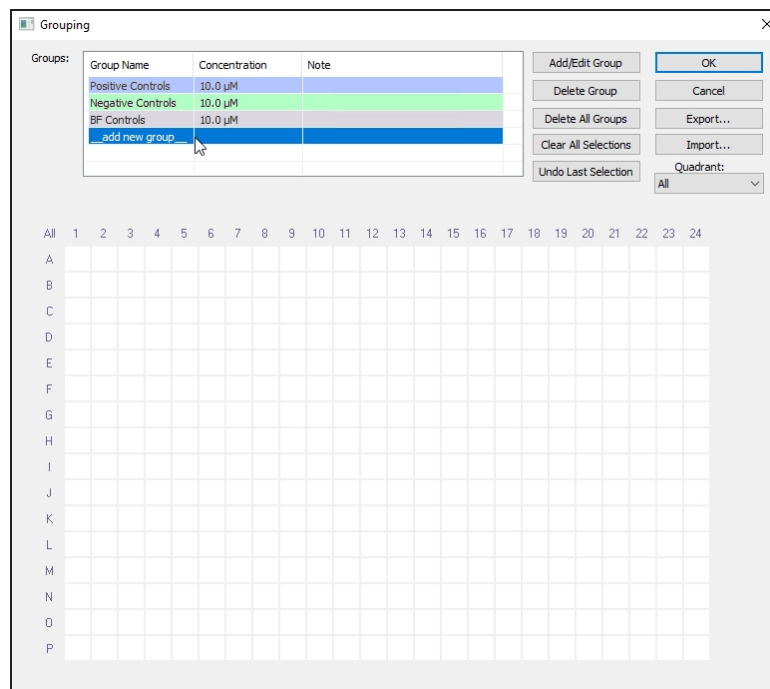
3. Repeat Steps 1–2 to assign other wells to different groups.

Adding a New Group

To add a new group:



1. Define groups in the **Analysis > Configure Groups > Grouping** dialog.
2. Select **_add new group** from the **Groups** list and click **Add/Edit Group**, or double-click on the group name.



- When the **Select Type of Group** dialog opens, choose to configure a single group with specific settings in the **Group** tab, or to configure a series of groups, with incrementing concentration values, in the **Series** tab.

Group Tab

To assign a single group concentration:

- Click **Group**.
- Type a name for the group (for example, compound name) in the **Group Name** field.
- In the **Concentration** field, select or type the concentration value.
- In the **Units** drop-down list, select **nM**, **µM**, **mM**, **M**, or **Log** concentration units.
- In the **Notes** field, type in any notes that you want associated with the specified group and data file.
- Click **OK** to finish save the defined group.

Series Tab

When a series is created, assign wells to it. See [Assigning Wells to Groups on page 88](#).

Successive rows or columns are assigned different concentrations and group color automatically, as configured in the series.



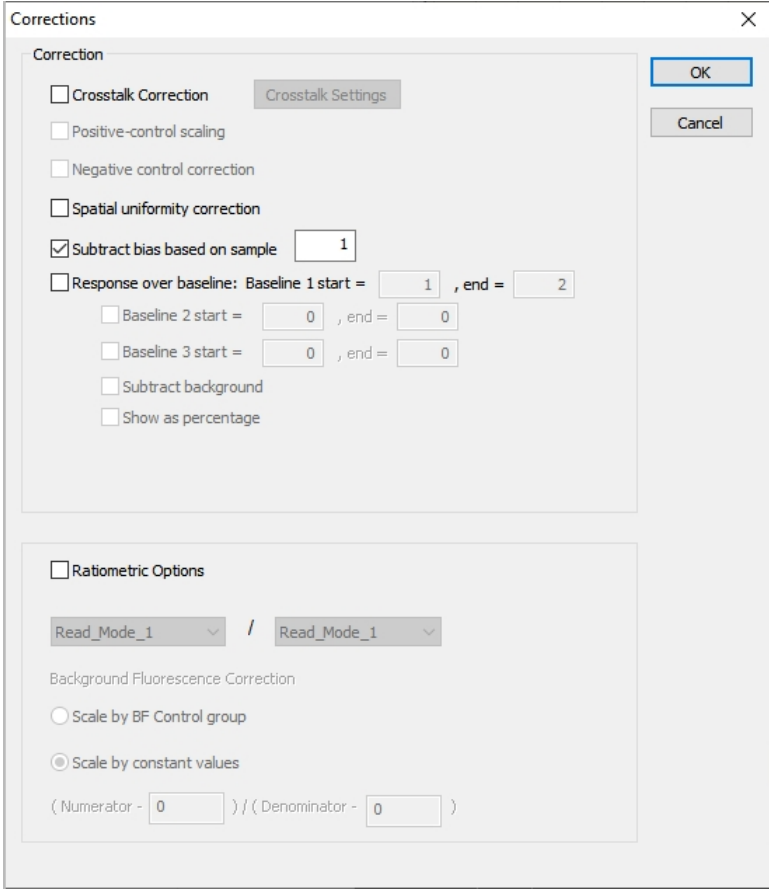
Note: The **Series** option is typically used for assigning a dose-response curve with regular increment steps.

To assign a series of concentrations:

1. Click **Series**.
2. In the **Group Name** field, type a name for the group, for example, compound name.
3. In the **Starting Value** field, select or type the value for the lowest or highest concentration of the series.
4. In the **Step Increment** field, select or type the value for the change in concentration between concentrations in the series.
5. In the **Units** drop-down list, select **nM**, **μM**, **mM** or **M** concentration units.
6. Specify the **Step Increment** for the **Starting Value** to create the series.
7. In the **Operation** drop-down list, select **Plus**, **Minus**, **Multiply** or **Divide**.
8. If you have **Replicates**, select whether they are aligned in **Row** or **Column** format on the plate.
 - If replicates align in both row and column formats, indicate the number of replicates, for example, replicates in rows indicates all wells in the selected rows have the same concentration.
 - If two rows are to include the same concentration, type 2.
 - If there are no replicates, select **None**.
9. Select the direction in which the series increments, either **Left**, **Right**, **Top** or **Bottom**.
10. In **Notes**, type details you want associated with the specified group and data file.
11. Click **OK** to finish save the defined group.

Corrections

Use the **Corrections** dialog to apply data-correction algorithms and to view ratiometric data.



Changes affect:

- Data displayed in the Multi-Well Graph and **Detail Graph**.
- Data displayed in group statistic table and graphs.
- Data in files are exported manually or automatically. See [Exporting Data on page 157](#).

Exported data cannot be changed, so settings made in this dialog in protocol files are important for these output options. For viewing data in ScreenWorks Software, however, since raw data are always kept, correction options selected in protocol files can be altered when the data are viewed in resulting data files.

Table 3-19: Corrections Dialog Options

| Options | Description |
|-------------------------------|---|
| Crosstalk Correction | Only available with Aequorin, ICCD camera. Check this option to correct all wells for crosstalk from neighboring wells. The crosstalk function is based on well-to-well distance and the measured percentage of crosstalk. |
| Positive Control Scaling | Check to average each maximum relative light units value from the positive control group and normalize all samples to this value (set at 100%). This function is useful when graphing a dose-response curve or when comparing data between experiments. |
| Negative Control Correction | Check this option to have an average negative control well value and ratio-to-well value calculated for each sample interval (time point). This ratio is calculated for each sample interval and applied to all wells. This function provides a good correction for signal drift and artifacts. |
| Spatial Uniformity Correction | <p>Check this option to use the initial signal to normalize signal in each well, removing fluctuations due to well-to-well variation in cell density and dye loading. The processing algorithm averages the initial signal from all wells together. It then scales each individual well relative to the average. This correction is particularly recommended when you want to normalize cell number, type and dye-loading conditions throughout a plate.</p> <p>Note: Spatial uniformity should only be used when all wells in a plate are treated the same prior to an experiment (for example, dye loading, cell numbers, etc.).</p> |
| Subtract Bias Based On Sample | Check this option to subtract the relative light units measured at a selected read number from all other time points in each well. This option enables you to set the Y-axis scale so that at the time point specified, the Y-axis values for all data graphs is zero. Select the read number to be subtracted in the field to the right of the check box. The default is read number 1. |

Table 3-19: Corrections Dialog Options (continued)

| Options | Description |
|------------------------|---|
| Response Over Baseline | <p>Check this option to display the trace as a ratio of the response to the average of a set of predefined reads.</p> <p>Baseline Start—First read to be included in the averaged baseline value.</p> <p>Baseline End—Last read to be included in the average baseline value.</p> <p>Show As Percentage—Displays response as a percent increase over the average baseline value.</p> |

For more details, see [Data Processing Algorithms on page 249](#).

Each of the correction options can be applied alone or in combination with the others. When a combination is selected they are applied in the following order:

1. Crosstalk
2. Spatial Uniformity Correction or Response Over Baseline
3. Negative Control Correction
4. Positive Control Scaling
5. Subtract Bias



Note: Spatial uniformity correction and percent baseline cannot be applied at the same time.

Ratiometric Options

Where two read modes are configured, you can view the data as a ratio of one read mode to the other, for each data point.

When this option is enabled, traces of the calculated ratio can be displayed in the Detail Graph pane.

Before calculating the ratio, ratiometric data is corrected for fluorescence background by subtracting estimated background values from both numerator and denominator for each time point. After the fluorescence background is subtracted, the ratio is then calculated. Additional selected corrections are applied to the ratio, beginning with negative control correction, positive control scaling, followed by subtract bias.



Note: Spatial uniformity correction and response over baseline are disabled when **Ratiometric Options** are active.

Table 3-20: Ratiometric Options

| Options | Description |
|---------------------|--|
| Ratiometric Options | <p>Enable ratiometric data viewing by selecting the check box. Specify how the ratio is defined by selecting the read modes to be used as numerator and denominator.</p> |

Table 3-20: Ratiometric Options (continued)

| Options | Description |
|------------------------------------|--|
| Background Fluorescence Correction | <p>Scale by BF Control Group—For each time point and read mode in the ratio, the average value for that time point in the background fluorescence control wells (BF Controls, in the Grouping dialog) is subtracted. The ratio is calculated after subtraction of the average value from both numerator and denominator.</p> <p>Scale by Constant Values—Constant values to subtract from the numerator and denominator for each time-point ratio are entered by the user.</p> <p>You must enter these constant values in the numerator and denominator fields when this option is selected.</p> |

Group Statistics

Below the Multi-Well and Detail Graphs on the **Analysis** process page, the **Group Statistics Table** is used to analyze data produced during an experiment. Only groups defined in the **Grouping** dialog can be analyzed in this section. In addition, settings made in the **Correction** dialog influence the group statistics that are reported. See [Corrections on page 91](#).

To analyze data in the **Group Statistics Table**, groups must first be defined in the **Grouping** dialog. See [Grouping on page 86](#). When defined, the **Configure Kinetic Reduction** button is selected in Multi-Well Graph to define the reads and reduction type to be analyzed.

Table 3-21: Group Statistics Table Button Options




| Buttons | Name | Description |
|---|---------------------|---|
|  | Copy Table Data | Copies data in the Group Statistic Table to the clipboard so it can be pasted in a different program such as Microsoft Excel. |
|  | Collapse All Groups | <p>Collapses the Group Statistics Table to only display the group name and relative statistics.</p> <p>This works as a toggle button with Expand All Groups which displays the individual wells and kinetic reduction values that comprise each group statistic.</p> |
|  | Expand All Groups | <p>Expands the Group Statistics Table to display all the individual wells and kinetic reduction values that comprise each group statistic.</p> <p>This works as a toggle button with Collapse All Groups, which displays only the group name and individual statistics.</p> |

Table 3-21: Group Statistics Table Button Options (continued)




| Buttons | Name | Description |
|---|----------------------|---|
|  | Auto Fit All Columns | This is used to collapse the column width so it is no wider than the title of the column or largest data point. |
|  | Select Statistics | <p>Opens the Choose Statistics dialog for defining groups statistics, for example, Average, Standard Deviation and Z-score, are displayed in the table.</p> <p>Group statistics available include the following:</p> <p>Concentration—Numerical concentration values assigned to each group in the Grouping dialog.</p> <p>Units—Concentration units (for example, uM) assigned to each group in the Grouping dialog.</p> <p>Notes—Comments assigned to each group in the Grouping dialog.</p> <p>Average—Numerical average of the kinetic reduction values for a given group.</p> <p>Maximum—Highest value (a single number) of all kinetic reductions within a group.</p> <p>Maximum-Minimum—Result of subtracting the minimum kinetic reduction value from the maximum count kinetic reduction value for a single group.</p> <p>Minimum—Lowest kinetic reduction value within a group.</p> <p>Sum—Numerical sum of all kinetic reduction values within a group.</p> <p>Standard Deviation—Defines the numerical value associated with one standard deviation from the average.</p> <p>Standard Deviation+1—Equals the average group value plus one standard deviation.</p> <p>Standard Deviation-1—Equals the average group value minus one standard deviation.</p> |

Table 3-21: Group Statistics Table Button Options (continued)

| Buttons | Name | Description |
|--|-------------------------------|---|
|  | Select Statistics (continued) | <p>Z Score—Used to evaluate quality or performance of the assay and is dependent on the concentration evaluated¹.</p> <p>Z-factor—Used to evaluate the quality or performance of the assay at a given concentration. This is typically used for all concentrations, not including the positive control. Calculation used:</p> $Z = 1 - [3 * (\text{std dev of GROUP} + \text{std dev of neg ctrls}) / \text{abs}(\text{mean of GROUP} - \text{mean of neg ctrls})]$ <p>Z'-factor—A characteristic parameter for the quality of the assay itself. This is typically performed using data only from the positive control concentration. Calculation used:</p> $Z' = 1 - [3 * (\text{std dev of pos ctrls} + \text{std dev of neg ctrls}) / \text{abs}(\text{mean of pos ctrls} - \text{mean of neg ctrls})].$ |
| <p>1. Zhang, J., Chung, T. D. Y. and Oldenburg, K. R. (1997). A Simple Statistical Parameter for Use in Evaluation and Validation of High Throughput Screening Assays. Journal of Biomolecular Screening 4(2):67:73.</p> | | |

The statistics defined in the **Group Statistic Table** can be plotted in the **Group Statistic Graph**, located behind the **Detail Graph**, to quickly view a dose-response curve. All groups are included in the graph, unless the user right-clicks on the desired group and selects **Exclude Groups From Statistic Chart**. Groups not used in the graph have a green line through the entire row of the group. To include the group in the graph, right-click the group and select **Include Groups In Statistic Chart**. Data displayed in the **Group Statistic Graph** is defined by selecting a specific group statistic from the **Group Statistic Table**. For Example, if the standard deviation is the needed data to be evaluated, select the **Standard Deviation** column header to highlight this data in the **Group Statistic Table** and it displays in the graph.

Table 3-22: Group Statistic Graph Options



| Buttons | Name | Description |
|---|-----------------|--|
|  | Copy Graph | Copies data in the Group Statistic Graph to the clipboard so it can be pasted in a different program such as Microsoft Word. |
|  | Copy Graph Data | Copies Graph Data in the Group Statistic Graph to the clipboard so it can be pasted in a different program such as Microsoft Excel. |

Table 3-22: Group Statistic Graph Options (continued)














| Buttons | Name | Description |
|---|--------------------------|---|
|  | Zoom Mode | Press to enable zoom capability in the graph. Once the Zoom button is pressed, drag over a region in the graph to view just that area enlarged to the full size of the graph. Alternatively, drag the cursor along a part of the X- or Y-axis to zoom in on just that axis. |
|  | Auto Scale | Manual prompt to automatically scale the Group Statistic Graph to include all data points of the desired traces. |
|  | Auto Scale Always | Automatically scales the Group Statistic Graph to include all data points of the desired traces without manual prompting. |
|  | Undo Zoom | Rescales graph to the original settings. |
|  | Manual Scale Graph | Opens the Manual Scale Graph dialog to set the maximum and minimum values for the X- and Y-axis graph. |
|  | Show/Hide Data | Toggle button that shows or hides the data for the selected traces at each read. |
|  | Show Point Labels | Show Point Labels writes the value of each data point, such as relative light units value, next to the point on the graph. |
|  | Show/Hide Legends | A toggle button that displays or hides the detail graph legend. |
|  | Show/Hide Smoothed Curve | A toggle button that displays or hides the 4-parameter curve fit. This feature can be displayed at the same time as Show/Hide Original Trace . |
|  | Show/Hide Original Trace | A toggle button that displays or hides the trace of the graph, which connects each data point. This feature can be displayed at the same time as Show/Hide Smooth Curve . |

Table 3-22: Group Statistic Graph Options (continued)


| Buttons | Name | Description |
|---|-----------------------------|---|
|  | Show/Hide Real Data Points | A toggle button that displays or hides the data points to create the graph. Note: When the data points are hidden, Molecular Devices recommends to activate either the Original Trace or Smoothed Curve function. |
|  | Display X-Axis in Log Scale | A toggle button that converts the X-axis from a concentration to a log concentration scale. |
|  | Show/Hide EC/ICXX Value | A toggle button that displays or hides the effective/inhibition concentration from being displayed. When activated, you define the percentage of activation/inhibition that the graph plots. |

Exporting Data

You can export data the following two ways:

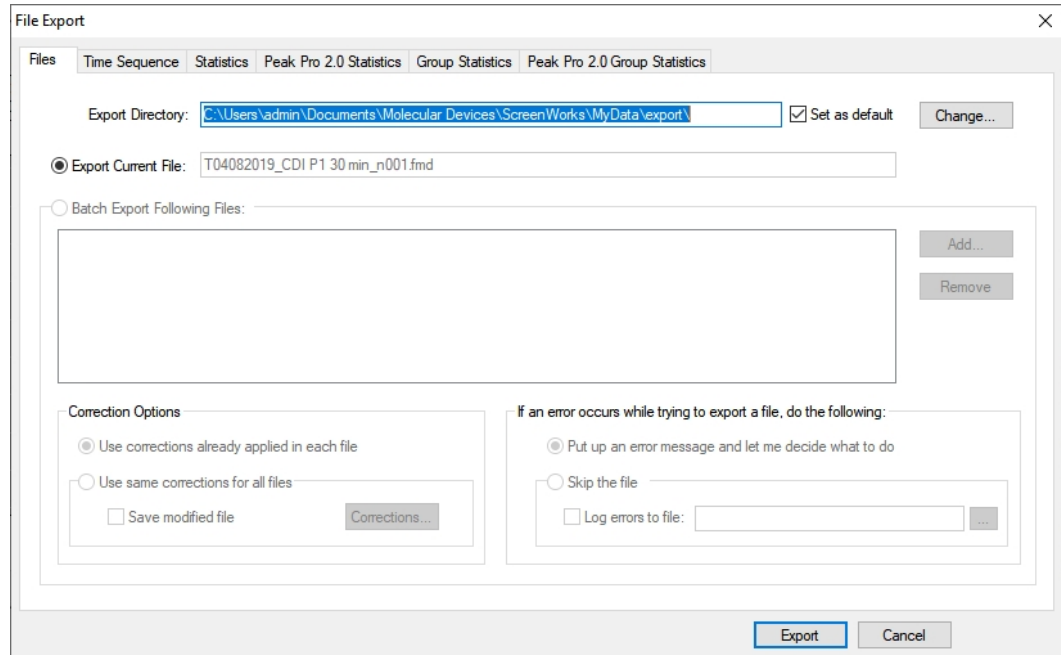
- Automatic export when a protocol is run.
- Export from already acquired data.

To configure the data export:

- In protocol files, in the **Analysis** screen, click  **Configure Auto Export**.
- For data files, use **File > Export, File** or **Batch Export**.



Note: **Auto-Export** and **Batch Export** do not support the ScreenWorks Peak Pro Software version 2.0.



Data is exported as ASCII text format files with a separate file exported for each measurement configuration specified. When you export data, you specify a folder in which to write the output files. The default export folder is:

C:\Documents and Settings\[your_user_name]\My Documents\Molecular Devices\ScreenWorks\MyData

The **File Export** dialog has the following tabs:

- **Files**—Export directory setting or batch files selections.
- **Time Sequence**—Exports time-point measurements for selected read modes. The measurement values that are exported have any corrections configured in the **Correction** dialog applied. If there are two read modes and **Ratiometric Options** is selected in the **Correction** dialog, you can also export the ratio for each time point. Files have a *.seqn extension, where n increments for each file generated in the export.
- **Statistics**—Exports averages, maximums, and other kinetic reduction values for selected numbers of reads for each well. Files have a *.statn extension, where n increments for each file generated in the export.
- **Peak Pro 2.0 Statistics**—You must run the Peak Pro 2.0 analysis before you can do the export. After running the analysis, this export works the same way as Statistics. See [Exporting Data With Peak Pro 2.0 Analysis on page 180](#).
- **Group Statistics**—Exports the group statistical values for selected numbers of reads for each group and are based on user-defined kinetic reduction settings. Files have a *.gstatn extension, where n increments for each file generated in the export.
- **Peak Pro 2.0 Group Statistics**—You must run the Peak Pro 2.0 analysis before you can do the export. After running the analysis, this export works the same way as Group Statistics. See [Exporting Data With Peak Pro 2.0 Analysis on page 180](#).



CAUTION! Data created in ScreenWorks Software version 2.0 cannot be viewed or exported by ScreenWorks Software version 5.0.

Exporting Time Sequence

Configure parameters for export of time-sequence data in the **Export Time Sequence** tab. The exported files contain a value for every read interval for each well, for the read mode or ratio selected. This contrasts with the Statistics file, which contains one kinetic reduction per well.

Table 3-23: Export Time Sequence Options

| Options | Description |
|----------------------|--|
| Enable Time Sequence | Check this option to have a time sequence export file created. |
| File Name Options | Choose a name for the exported time sequence file. Use name of data file —Export files with the data or protocol file name, with *.seq <i>n</i> extension, where <i>n</i> is an integer ≥ 1 . Use user-defined name —Enter your own name for the export files (maximum of 25 characters). Files are given *.seq <i>n</i> extension, where <i>n</i> is an integer ≥ 1 . |
| Data Format | Format the output from this group of options: Show Labels —Check to include information about the processing options selected (for example, name of *.fmd file and well labels). Any corrections prior to export (e.g., negative control correction) are reported at the top of the exported ASCII file. Order of Well Data —Order the data by column (for example, A1, B1, C1) or by row (for example, A1, A2, A3). Set this option in accordance to the way your spreadsheet or database handles well data. Order Time Data —Select vertical or horizontal ordering of the data. |
| Misc. Parameters | Select experimental parameters and group statistics to export with the data. |

Table 3-23: Export Time Sequence Options (continued)

| Options | Description |
|----------------------------|---|
| Individual File Parameters | <p>Use the bottom section of the dialog to choose the number and type of time sequence files to be created from the same data set.</p> <p>Number of Files—Click the Add Statistic button to add to the number of statistic files to be created from the same data set.</p> <p>Read Mode—Enter the read mode (or ratiometric data) from which you want to export data.</p> <p>From Quadrant—You can choose a specific quadrant from which to export, or export data from all quadrants. This enables you to export data from one plate into four different spreadsheet files, if desired.</p> <p>Note: This option is only available for data created in 384-well or 1536-well formats.</p> |

Exporting Statistics

Use the **Export Statistics** tab to configure kinetic reductions, for example, averages and maximum and minimum values, for selected numbers of reads within each well. Multiple kinetic reductions can be configured, each resulting in another export file. Within each file only one value is reported per well.

Table 3-24: Export Statistics Options

| Options | Description |
|-------------------|---|
| Enable Statistics | Check this box to have a statistic file created. |
| File Name Options | <p>Choose a name for the exported statistic file.</p> <p>Use name of data file—Export files with the data or protocol file name, with *.stat<i>n</i> extension, where <i>n</i> is an integer ≥ 1.</p> <p>Use user-defined name—Enter your own name for the export files (maximum of 25 characters). Files are given *.stat<i>n</i> extension, where <i>n</i> is an integer ≥ 1.</p> |
| Data Format | <p>Format the output from this group of options:</p> <p>Show Labels—When checked, the output file contains information about the processing options selected (for example, name of *.fmd file and well labels). Any corrections done prior to exporting the data (e.g., negative control correction) are indicated in text format at the top of the exported ASCII file.</p> <p>Order of Well Data—Order the data by column (for example, A1, B1, C1) or by row (for example, A1, A2, A3). Set this option in accordance to the way your spreadsheet or database handles well data.</p> |

Table 3-24: Export Statistics Options (continued)

| Options | Description |
|----------------------------|---|
| Misc. Parameters | Select experiment parameters and group statistics to export with the data. |
| Individual File Parameters | <p>Use the bottom section of the dialog to designate the number and type of statistic files to be created from the same data set.</p> <p>Number of Files—Click the Add Statistic button to add to the number of statistic files to be created from the same data set.</p> <p>Compute—Select a kinetic reduction type. See Kinetic Reduction Types on page 81</p> <p>Start Read—Enter the first read number to be included in the kinetic reduction.</p> <p>End Read—Enter the last read number to be included in kinetic reduction. This can equal the Start Read, if you want to extract values from a single read.</p> <p>Read Mode—Enter the read mode (or ratiometric data) from which the data processes.</p> <p>From Quadrant—You can choose a specific quadrant from which to export, or export data from all quadrants. This enables you to export data from one plate into four different spreadsheet files, if desired.</p> <p>Note: This option is only available for data created in 384-well or 1536-well formats.</p> |

Exporting Group Statistics

Use the **Export Group Statistics** tab to configure a group statistics export report, for example, average, standard deviation and z-scores, based on the kinetic reduction types defined for a select number of reads as set up on the Analysis screen.



Note: If there are no groups defined on the Analysis screen, an empty report exports.

Table 3-25: Export Group Statistics Options

| Options | Description |
|-------------------------|---|
| Enable Group Statistics | Check this box to have a group statistic file created. |
| File Name Options | <p>Choose a name for the exported group statistic file.</p> <p>Use name of data file—Export files with the data or protocol file name, with *.gstat<i>n</i> extension, where <i>n</i> is an integer ≥ 1.</p> <p>Use user-defined name—Enter your own name for the export files (maximum of 25 characters). Files are given *.gstat<i>n</i> extension, where <i>n</i> is an integer ≥ 1.</p> |

Table 3-25: Export Group Statistics Options (continued)

| Options | Description |
|-------------------|--|
| Kinetic Reduction | <p>Configures the parameters used to define the kinetic reduction.</p> <p>Reduction Type—Defines the reduction to be applied to the kinetic data traces exported. See Kinetic Reduction Types on page 81 for details.</p> <p>Start Read—Define the first read to be used to determine the kinetic reduction.</p> <p>End Read—Define the last read used to determine the kinetic reduction.</p> <p>Read Mode—Select the read mode to apply the kinetic reduction.</p> |
| Misc. Parameters | Select experiment parameters to export with the data. |

Batch Exporting



Note: Batch Export does not support the separately licensed analysis module ScreenWorks Peak Pro Software version 2.0.

Manual export of files is accessed through **File > Export, File or Batch Export**. **Manual Export** allows the same files as defined above to be exported.

Table 3-26: Batch Export Options

| Options | Description |
|------------------------------|---|
| Export Directory | <p>User defines where export files are to be sent. The default export folder is:</p> <p>C:\Documents and Settings\[your_user_name]\My Documents\Molecular Devices\ScreenWorks\MyData</p> |
| Batch Export Following Files | <p>Selecting Add opens the Open File dialog to choose data files to be exported.</p> <p>Note: Hold down the SHIFT or CTRL key to select data files.</p> |
| Correction Options | <p>During export, either the existing or new corrections can be applied.</p> <p>Use corrections already applied in each data file—Applies the existing corrections saved with each data file during export.</p> <p>Use same corrections for all files—Applies and saves new corrections to all data files. Useful when exporting all data files with same parameters or saving the same correction to multiple data files. Details regarding available corrections can be found under Corrections on page 91.</p> |

Table 3-26: Batch Export Options (continued)

| Options | Description |
|--|--|
| If an error occurs while trying to export a file, do the following | <p>Put up an error message and let me decide what to do— Enables the user to decide how to proceed when an error is encountered.</p> <p>Skip the file—Does not include the data in the export, but continues to export data. The option to write the error to a log file is available.</p> |

When batch exporting data, the export files produced vary depending on the options you select in the individual export statistic, time sequence and group statistic sections. If a single export file is needed for each data file exported, select **Use name of data file** in the respective section you want to export. Export files created using this method are labeled with the name of the data file from which it was created. However, if a single export file is needed that contains information for multiple data files, select **Use user-defined name** in the respective section. In this instance, all exported information combines into one file labeled with the desired user-defined name. When the file is open, the individual data file names which the information was exported from are used as the header for each data set within the export file.

Image Display

Use the **Image Display** dialog, opened from the **Analysis** process page, to view images saved in data files when **Save Images** was enabled during a **Read with TF** step. This option stores a total of 100 images per experiment; the number of images per dispense is user-defined. The default is set to one image before and nineteen images after the fluid addition is initiated. Saved *.tif files can be played back in sequence, frame-by-frame or as a video.

Image Display is useful for diagnosing problems. For example, if cells are blown off during a fluid dispense, you see dark holes in the middle of the cell layer after the fluid addition. If the entire well decreases in relative light units, it is likely that extracellular dye has been diluted.

Notes

In protocol files, open the **Protocol Notes** dialog where you can enter comments that you want kept with all data files generated with the protocol. The dialog contains a simple text editor for you to format your comments.

In data files the same dialog shows the comments that were added in the protocol file, now in read-only format, so they cannot be edited.

Transfer Fluid Process

This process enables you to program reagent or cell additions and to execute them in conjunction with other processes.

The default protocol has one **Transfer Fluid** process.

A blue **Transfer Fluid** icon indicates that the process occurs in series with other blue-colored processes in the protocol. A purple icon indicates that the execution of the process is synchronized to a **Read** or **Mix Fluid** process, which also has a purple icon.

The **Transfer Fluid** process has different options for pin tools and standard pipettors. In each case you must configure aspiration and dispense of compound separately.

For a standard pipettor you can configure more than one aspiration or dispense (but not both) within the process, with the appropriate 'multiple' **Fluid Transfer Type**.

With a standard pipettor selected, aspiration and dispense steps are configured by selecting the appropriate row in the aspiration or dispense table and pressing the **Edit** button to open a configuration dialog. With the Pin Tool, aspiration and dispense configuration is done directly in the **Transfer Fluid** page.

To enable additional dispense steps with a standard pipettor, select the multiple dispense option in the **Fluid Transfer Type**. Then configure the first dispense with a volume less than the volume aspirated. An additional dispense is automatically enabled for the remaining volume. Similarly, for multiple aspirates, configure the first aspirate with less volume than the dispense.



Note: The entire volume of fluid aspirated is always dispensed.

Aspirate Configuration (Standard Pipettor)

Pipettor type: 384 Settings Analysis Mix With TF Transfer Fluid Read With TF

Transfer Fluid

Fluid Transfer Type: Single aspirate - Single dispense Read

| Aspirate | Plate | Volume... | Height(µl) | Speed(...) | Tip Up ... | HoldVo... | Quadrant | Done | Fill Res... | Drain R... | Drain R... |
|--|----------------|-----------|------------|------------|------------|-----------|----------|------|-------------|------------|------------|
| <input checked="" type="checkbox"/> Aspirate | Source Plate 2 | 13.5 | 20.0 | 20.00 | 20 | 0.0 | NONE | NO | N/A | N/A | N/A |
| <input type="checkbox"/> Aspirate | Source Plate 1 | 0.0 | 4.6 | 20.00 | 20 | 1.0 | NONE | NO | N/A | N/A | N/A |
| <input type="checkbox"/> Aspirate | Source Plate 1 | 0.0 | 4.6 | 20.00 | 20 | 1.0 | NONE | NO | N/A | N/A | N/A |
| <input type="checkbox"/> Aspirate | Source Plate 1 | 0.0 | 4.6 | 20.00 | 20 | 1.0 | NONE | NO | N/A | N/A | N/A |

Put Tips In Well Before Read Mix Fluid before Aspirate Edit Aspirate

| Dispense | Plate | Volume... | Height(µl) | Speed(...) | ExpeVol(µl) | Pause(s) | Removal Spe... | Quadrant | Done | Hold Pi... |
|--|------------|-----------|------------|------------|-------------|----------|----------------|----------|------|------------|
| <input checked="" type="checkbox"/> Dispense | Read Plate | 13.5 | 20.0 | 30.0 | 0.0 | 0 | 20 | NONE | NO | NO |
| <input type="checkbox"/> Dispense | Read Plate | 0.0 | 46.0 | 20.0 | 1.0 | 0 | 20 | NONE | NO | NO |
| <input type="checkbox"/> Dispense | Read Plate | 0.0 | 46.0 | 20.0 | 1.0 | 0 | 20 | NONE | NO | NO |
| <input type="checkbox"/> Dispense | Read Plate | 0.0 | 46.0 | 20.0 | 1.0 | 0 | 20 | NONE | NO | NO |

Mix Fluid after Dispense Unload Tips After Fluid Transfer Edit Dispense

Table 3-27: Aspirate Configuration Options for Standard Pipettor Use

| Option | Description |
|---------------------|---|
| Fluid Transfer Type | <p>Select a type of fluid transfer:</p> <p>Single Aspirate–Single Dispense—Aspirates one liquid volume and dispenses that entire liquid volume in one action into a plate.</p> <p>Single Aspirate–Multiple Dispense—Aspirates one liquid volume and dispenses up to four distinct liquid volumes into four distinct quadrants or plates. Add additional dispenses by making the dispense volume less than the aspirate volume.</p> <p>Note: The last dispense expels all liquid remaining in the tip. Tips to improve Single Aspirate–Multiple Dispense (by quadrature) pipetting precision include:</p> <ul style="list-style-type: none"> * When pipetting do not use hold and expel volumes. In addition use the slowest removal speed (2 mm/s). * Aspirate 5–10% more than is needed from the source plate (for example, aspirate 110 μL when dispensing 25 μL into four quadrants). * Include a mix step with 3 strokes in the desired source plate to pre-wet tips. Mixing of fluid in the source plate primes the internal walls of the tips, reducing surface tension between the fluid and tip plastic. <p>Multiple Aspirate–Single Dispense—Aspirates up to four liquid volumes from up to four aspirate plates and dispenses entire tip contents into a single plate. To add additional aspirates, make the aspirate volume less than the dispense volume.</p> |
| Read | <p>Check this box to link a Read process to the Transfer Fluid process. This automatically adds a Read with TF step to the protocol, and opens the configuration page for that step.</p> <p>See Read Process on page 116 for more information about read process options.</p> <p>Note: This option is required if you intend to read a plate while transferring liquid during a kinetic cell-based assay (for example, calcium mobilization).</p> |
| Aspirate | <p>Check the check box at the start of a row to activate an aspirate sequence.</p> <p>If Multiple Aspirate has been selected as Fluid Transfer Type, enter an aspirate volume less than the dispense volume to add another aspirate step.</p> <p>Note: Only checked aspirate sequences are active in the experimental protocol.</p> |

Table 3-27: Aspirate Configuration Options for Standard Pipettor Use (continued)

| Option | Description |
|--|---|
| Edit Aspirate | <p>Configure an aspirate step in the table by selecting it and pressing the Edit Aspirate button, or double-click on the row. This opens the Edit Aspirate dialog.</p> <p>Use this dialog to define aspiration parameters reported in the table.</p> <p>Source Plate—Select the plate from which you want an aliquot of fluid removed. Typically, this is a plate containing an agonist, antagonist or if the Cell Suspension option is installed, cells in suspension.</p> <p>Quadrant—Select quadrant from which volume is aspirated.</p> <p>Note: Quadrant selection is only available when liquid is aspirated from a plate using a pipettor head of lower density, for example, 384-well plate using a 96 pipettor head.</p> <p>Volume—Enter volume (in μL) to aspirate.</p> <p>Speed—Enter speed at which to aspirate (in $\mu\text{L/s}$).</p> <p>Fill Reservoir Speed—Select the speed with which fill the Cell Reservoir from the spinner flask.</p> <p>Drain Reservoir Speed¹—Part of the Cell Suspension System. Select the speed with which to drain the cells from Cell Reservoir to drain destination.</p> <p>Drain Destination¹—Select a destination where cells should be drained from Cell Reservoir after pipettor aspiration.</p> <p>Height—Enter distance (measured in μL) from the bottom of the well the tips should start aspirating.</p> <p>Note: Tips move downward full volume of aspiration, for example, aspirate 10 μL with a start height of 35 μL moves the pipettor height to 25 μL after aspiration. Pipettor head movement helps prevent well overflow when tips are submerged in the well. If downward pipettor motion is not desired, set tip height to minimum value. Tip movement downward halts when the tips reach 1/30th well volume.</p> <p>Hold Volume—Enter size of air gap (in μL) following aspiration. Air gaps prevent liquid from leaking out of the tip before dispensing.</p> |
| Edit Aspirate (continued) | <p>Removal Speed—Enter rate (in mm/s) at which tips are pulled from well immediately after aspiration. This value is adjusted depending on fluid viscosity or volume of fluid in the well.</p> <p>Note: For low volume transfers, a slow up-speed may improve precision.</p> |
| Put Tips in Well Before Read | <p>Check this option to place tips in the read plate prior to beginning read intervals.</p> <p>This option is suggested for assays in which placing tips below the fluid surface can prevent excessive turbulence.</p> <p>Enabling this option helps ensure accurate timing of fluid dispense relative to reads.</p> |
| <p>1. These options are only available when Cell Reservoir is installed and selected in Source Plate 3 settings.</p> | |

Dispense Configuration (Standard Pipettor)

Pipettor type: 384 Settings Analysis Transfer Fluid Read With TF

Transfer Fluid

Fluid Transfer Type: Single aspirate - Single dispense Read

| Aspirate | Plate | Volume... | Height(µl) | Speed(...) | Tip Up ... | Hold/Vo... | Quadrant | Done | Fill Res... | Drain R... | Drain R... |
|--|----------------|-----------|------------|------------|------------|------------|----------|------|-------------|------------|------------|
| <input checked="" type="checkbox"/> Aspirate | Source Plate 2 | 12.5 | 5.0 | 20.00 | 20 | 1.0 | NONE | NO | N/A | N/A | N/A |
| <input type="checkbox"/> Aspirate | Source Plate 1 | 0.0 | 4.7 | 20.00 | 20 | 1.0 | NONE | NO | N/A | N/A | N/A |
| <input type="checkbox"/> Aspirate | Source Plate 1 | 0.0 | 4.7 | 20.00 | 20 | 1.0 | NONE | NO | N/A | N/A | N/A |
| <input type="checkbox"/> Aspirate | Source Plate 1 | 0.0 | 4.7 | 20.00 | 20 | 1.0 | NONE | NO | N/A | N/A | N/A |

Put Tips In Well Before Read Mix Fluid before Aspirate Edit Aspirate

| Dispense | Plate | Volume... | Height(µl) | Speed(...) | ExpelVol(µl) | Pause(s) | Removal Spe... | Quadrant | Done | Hold Pt... |
|--|------------|-----------|------------|------------|--------------|----------|----------------|----------|------|------------|
| <input checked="" type="checkbox"/> Dispense | Read Plate | 13.5 | 40.0 | 50.0 | 0.0 | 0 | 20 | NONE | NO | NO |
| <input type="checkbox"/> Dispense | Read Plate | 0.0 | 46.7 | 20.0 | 1.0 | 0 | 20 | NONE | NO | NO |
| <input type="checkbox"/> Dispense | Read Plate | 0.0 | 46.7 | 20.0 | 1.0 | 0 | 20 | NONE | NO | NO |
| <input type="checkbox"/> Dispense | Read Plate | 0.0 | 46.7 | 20.0 | 1.0 | 0 | 20 | NONE | NO | NO |

Mix Fluid after Dispense Unload Tips After Fluid Transfer Edit Dispense

Table 3-28: Dispense Configuration Options for Standard Pipettor Use

| Options | Description |
|---------------|---|
| Dispense | <p>Check the check box at the start of a row to activate a dispense sequence.</p> <p>If Multiple Dispense has been selected as Fluid Transfer Type, enter a dispense volume less than the aspirate volume to add another dispense step.</p> <p>Note: Only checked dispense sequences are active in the experimental protocol.</p> |
| Edit Dispense | <p>Configure a dispense step in the table by selecting it and clicking Edit Dispense, or double-click on the row to open the Edit Dispense dialog.</p> <p>Use this dialog to define dispense parameters reported in the table.</p> <p>Target Plate—From the list, select the plate to receive the fluid dispense. Typically, this is the Read Plate.</p> <p>Quadrant—Select quadrant from which volume is dispensed.</p> <p>Note: Quadrant selection is only available when liquid is aspirated from a plate using a pipettor head of lower density for example, 384-well plate using a 96-pipettor head.</p> <p>Volume—Enter volume (in µL) to dispense.</p> <p>Note: In the last dispense, the entire contents in the tip is dispensed.</p> |

Table 3-28: Dispense Configuration Options for Standard Pipettor Use (continued)

| Options | Description |
|---------------------------|---|
| Edit Dispense (continued) | <p>Height—Enter the distance (measured in μL) from the bottom of the well the tips should be inserted prior to dispensing.</p> <p>Note: Tips move upward the full volume of dispense during dispense, for example, a 10 μL dispense with a start height of 35 μL moves the pipettor to a height of 45 μL after dispensing. This movement helps prevent well overflow when tips are submerged in the well. Tips do not move higher than the full volume of the well.</p> <p>Height value needs to be adjusted when transferring from previous FLIPR Tetra Systems.</p> <p>Speed—Enter speed at which to dispense (in $\mu\text{L}/\text{s}$). Ideal dispense speed takes into account the volume added, how fast the signal increases in response to compounds and strength of the attached cells at the bottom of the plate.</p> <p>Expel Volume—Enter an additional volume (in μL) the pipettor dispenses over the dispense volume. Typically, this value is equal to the aspirate Hold Volume value.</p> <p>In a multiple dispense, only the last dispense can have an Expel Volume.</p> <p>If Mix Fluid After Dispense is selected, Expel Volume automatically changes to zero for the dispense associated with the fluid transfer, to ensure that tips remain in the well prior to the mix step. However, the expel volume designated in the dispense is transferred to the Mix with TF step, to help ensure liquid is properly expelled from the tips after mixing.</p> <p>Removal Speed—Enter rate (in mm/s) at which tips are pulled from well immediately after the dispense. This value is adjusted for fluid viscosity, fluid volume in the well or if the transfer is performed to a dry well.</p> <p>Pause in Well—Enter time (in seconds) to pause with the tips in the wells after dispensing and mixing (if specified) before the next pipettor move.</p> <p>Hold pipettor during dispense—Select this check box to hold the pipettor at your specified dispense height.</p> <p>Read intervals might occur simultaneously with this step. Typically, the pipettor is paused during a read to prevent disturbances in the imaging process.</p> |

Table 3-28: Dispense Configuration Options for Standard Pipettor Use (continued)

| Options | Description |
|----------------------------------|--|
| Mix Fluid After Dispense | <p>Check this box to link a Mix Fluid process to the Transfer Fluid process.</p> <p>This automatically adds a Mix with TF step to the protocol, and opens the configuration page for that step.</p> <p>See Mix Fluid Process on page 113 for more information about mixing options.</p> <p>Note: This option is only available when Fluid Transfer Type has a single dispense.</p> |
| Unload Tips after Fluid Transfer | <p>Instructs the instrument to unload tips after the Fluid Transfer process. Typically, this is used when changing the tips during an experiment or when reducing down-time between experiments.</p> <p>Note: This option is only available when Load Tips Position is selected on the Settings process page. Using this feature also requires that no source plate is placed in plate position 3 during the experiment.</p> |

Aspirate Configuration (Pin Tool)

Pipettor type: 384pin tool Settings Analysis Transfer Fluid Read With TF

Aspirate

Source Plate: Source Plate 1 default384 v

Height (µL): -12.00 = -1.00 mm Float pins

Quadrant: None v

Down Speed (mm/s): 10.0

Removal Speed (mm/s): 10.0

Pause In Well (s): 0

Fill Reservoir Speed: v Mix Fluid before Aspirate

Drain Destination: v

Drain Reservoir Speed: v

Dispense

Target Plate: Read Plate default384 v

Height (µL): 1.00 = 0.08 mm Float pins

Quadrant: None v

Down Speed (mm/s): 10.0

Removal Speed (mm/s): 10.0

Pause In Well (s): 0 Read

After Dispense

Mix Fluid Unload Pin Tool

Table 3-29: Aspirate Configuration Options for Pin Tool Use

| Options | Description |
|---------------|--|
| Source Plate | Select the plate from which you want an aliquot of fluid removed. Typically, this is a plate containing an agonist or antagonist. |
| Height | Enter distance (measured in µL) from the bottom of the well to place the pins. The distance equivalent in millimeters is displayed as well. Float pins —Moves the pipettor down to a position at which the pins should be 1 mm below the bottom of the wells. As pins come to rest on the well bottoms additional downward motion of the tip block is absorbed by the tips moving into the block. This ensures that all pins are submerged in the same volume of fluid across the plate. The Height is designated as -1.00 mm. |
| Quadrant | Select the quadrant from which volume is aspirated. Note: Quadrant selection is only available when liquid is aspirated from a plate using a pipettor head of lower density, for example, 1536-well plate using a 384 pin tool head. |
| Down Speed | Enter the speed (in mm/s) at which the pin tool is inserted into the source plate. |
| Removal Speed | Enter the speed (in mm/s) at which the pin tool is removed from the source plate. The speed of removal determines the volume of fluid remaining on the pins. Note: Increasing the removal speed increases the volume of compound on the pin after aspiration. To decrease the fluid volume, slow down the removal speed. It is recommended that assay development be done to determine the volume transferred based on removal speed. |

Table 3-29: Aspirate Configuration Options for Pin Tool Use (continued)

| Options | Description |
|---------------|---|
| Pause in Well | Enter a duration (in seconds) to hold the pin tool in the source plate after full insertion and before removal. |

Dispense Configuration (Pin Tool)



Table 3-30: Dispense Configuration Options for Pin Tool Use

| Options | Description |
|---------------|---|
| Target Plate | Select the plate to receive the fluid dispense; typically, the Read Plate. |
| Height | Enter the distance (measured in µL) from the bottom of the well the tips should be inserted. The distance equivalent in millimeters is displayed to the right of the µL value. |
| Quadrant | Select the quadrant of the target plate into which compound is to be dispensed. Note: Quadrant selection is only available when liquid is aspirated from a plate using a pipettor head of lower density, for example, 1536-well plate using a 384 pin tool head. |
| Down Speed | Enter the speed (in mm/s) at which the pin tool is inserted into the target plate. |
| Removal Speed | Enter the speed (in mm/s) at which the pin tool is removed from the target plate. The speed of removal determines the volume of fluid remaining on the pins. |
| Pause in Well | Enter a duration (in seconds) to hold the pin tool in the dispense plate after full insertion and before removal. |

Table 3-30: Dispense Configuration Options for Pin Tool Use (continued)

| Options | Description |
|-----------------|--|
| Read | <p>Check this box to link a Read process to the Transfer Fluid process. This automatically adds a Read with TF step to the protocol, and opens the configuration page for that step.</p> <p>See Read Process on page 116 for more information about read process options.</p> <p>Note: This option is required if you intend to read a plate while transferring liquid during a kinetic cell-based assay (for example, calcium mobilization).</p> |
| Mix Fluid | <p>Check this box to link a Mix Fluid process to the Transfer Fluid process.</p> <p>This automatically adds a Mix with TF step to the protocol, and opens the configuration page for that step.</p> <p>See Mix Fluid Process on page 113 for more information about mixing options.</p> |
| Unload Pin Tool | Automatically unload the pin tool after the dispense. |

Mix Fluid Process

Fluid mixing steps can be added to protocols as independent protocol steps, **Mix Fluid** or linked to a **Transfer Fluid** step, **Mix with TF**. Independent Mix Fluid steps can be configured to mix fluid in source or read plates, while Mix with TF steps immediately follow the Transfer Fluid step they are linked to, mixing fluid in the read plate where compound was dispensed. Different mix options are offered for standard pipettors and the pin tool, described in the following sections.

Mix Fluid (Standard Pipettor)

Fluid mixing with a standard pipettor consists of a series of aspirations and dispenses in and out of the tips.

The following table describes the options available on the **Mix Fluid** process page:

| Options | Description |
|-----------|--|
| Mix Plate | Select the plate that has the fluid to be mixed. Typically, it is a source plate from which you transfer fluid. |
| Quadrant | <p>If appropriate, select a quadrant from which fluid is mixed.</p> <p>Note: Quadrant selection is only available when liquid is dispensed to a plate using a pipettor head of lower density, for example, 384-well plate using a 96 pipettor head.</p> |
| Volume | Enter volume (in μL) to mix. |
| Speed | Enter speed (in $\mu\text{L/s}$) to mix the fluid. |

| Options | Description |
|-----------------------|--|
| Height | Enter distance (in μL) from the bottom of the well that the tips will be inserted prior to mixing. Note: Tips move up and down with aspirate and dispense commands. These movements help to prevent well overflow when the tips are submerged. |
| Expel Volume | Enter an additional volume (in μL) the pipettor dispenses, beyond the volume entered in the Volume field above. This helps to ensure that all fluid is expelled from the tips after mixing. |
| Strokes | Enter number of times to aspirate and dispense mix volume. One stroke is equal to one aspiration and one dispense. |
| Removal Speed | Enter rate (in mm/s) at which tips are pulled from well immediately after the mix. This value is adjusted for fluid viscosity or fluid volume in the well. |
| Pause In Well | Enter time (in seconds) to pause with the tips in the well after mixing and before the next pipettor head move. |
| Unload Tips After Mix | Automatically unloads tips after the mix step. |

Mix Fluid (Pin Tool)

Fluid mixing with a pin tool consists of a sequence of up and down motions of the pin block.

| Options | Description |
|---------------|---|
| Mix Plate | Select the plate that has the fluid to be mixed. Typically, it is a source plate from which you transfer fluid. |
| Top Height | Enter the height (in μL) for the top of the stroke during mixing. |
| Quadrant | If appropriate, select a quadrant in which fluid is mixed. Note: Quadrant selection is only available when liquid is dispensed to a plate using a pipettor head of lower density, for example, 1536-well plate using a 384 pipettor head. |
| Bottom Height | Enter the height (in μL) for the bottom of the stroke during mixing. |
| Strokes | Enter the number of times the pins move within the mix step. One stroke is equal to one up-stroke and one down-stroke. |
| Down Speed | Enter the speed (in mm/s) of the down-stroke. |
| Removal Speed | Enter rate (in mm/s) at which pins are pulled from the well immediately after the mix. This value is adjusted for fluid viscosity or fluid volume in the well. The speed of removal determines the volume of fluid remaining on the pins. |
| Up Speed | Enter the speed (in mm/s) of the up-stroke. This is not the rate which pins remove from the well. |

| Options | Description |
|-----------------|--|
| Pause in Well | Enter a duration to hold the pin tool in the well plate once mixing is completed, before withdrawal. |
| Unload Pin Tool | Check to have the pin tool unloaded after mixing. |

Mix With TF

Fluid can be mixed immediately after dispense to ensure proper mixing of agonists and antagonists with the cell monolayer. Typically, this is necessary when transferring fluids one-tenth or less of the total fluid volume of a well.

To add a mix after dispense, check the **Mix After Fluid Dispense (Mix Fluid for the Pin Tool)** option in the **Transfer Fluid** process dialog. This inserts a mix step and the **Mix Fluid** process icon turns purple, indicating that mixing occurs immediately after dispense and concurrently with a Read process, if selected.



Note: Any **Expel Volume** configured in the dispense step prior to a mix is locked if **Mix Fluid After Dispense** is selected. Tips are raised during Expel Volume, but this could lead to inadequate mixing, so mixing occurs with the tips at the level they finish the basic dispense at, and **Expel Volume** occurs after mixing is completed.

Mixing parameters for a **Mix with TF** are similar to a **Mix Fluid** process, however **Mix Plate, Quadrant, Hold Volume** and **Tip Up Speed**, are not configurable. These options can only be accessed in the **Transfer Fluid** process to which the mix is linked.

To remove a **Mix with TF** process, deselect it in the **Transfer Fluid** process dialog.

Read Process

Read steps can be added to protocols as independent protocol steps, **Read** or linked to a **Transfer Fluid** step, **Read with TF**. Configuration options for the two types of step is similar, except **Read with TF** allows you to time the read precisely with the time that compound is dispensed, and it also gives you the option to save the images used to measure the response.



Note: When using the HS EMCCD camera, depending on your **Protocol > Settings > Edit read mode > Camera Mode** setting, when using a fast Read time interval value, the plate matrix panel updating might be delayed until the end of the data acquisition rather than updating during the data acquisition.

Specifically, the following settings result in the plate matrix panel updating at the end of the data acquisition:

For **Camera Mode > HighSpeed**—a **Read time interval(s)** setting of 0.018 seconds or shorter

For **Camera Mode > Normal or Sensitivity**—a **Read time interval(s)** setting of 0.029 or shorter

For **Camera Mode** setting details, see [Edit Read Mode on page 73](#).

Read

To add a Read process to a protocol that is independent of a liquid transfer, drag the **Read** process icon from the **Process Explorer** to the Experiment window. The Read process page opens when the icon is released. When added to the protocol, the icon is orange, indicating it is an optics function and occurs in parallel with fluid processes during the experiment.

A Read consists of a sequence of images taken at a predefined rate, from which relative light values for each well are measured. A single Read process can be divided into two intervals with different read rates. Typically this is used to take images at a high rate while compound is being added, and shortly thereafter, to best capture kinetic effects, then images can be taken at a less frequent rate as cells settle and less change occurs.



Note: The read rate you configure for the Read process is the same for all the read modes you have enabled. For example, if you set the Read Interval at one second, and have two read modes, two images are taken every second—one for each read mode.

Pipettor type: 384

Settings Analysis Transfer Fluid Read

First Interval

Read time interval (s):

Number of reads:

Second Interval

Read time interval (s):

Number of reads:

The following table describes the setting options in the dialog:

Table 3-31: Read Process Configuration Options Dialog

| Options | Description |
|-----------------|--|
| First Interval | <p>Configure the first series of reads taken during the Read process.</p> <p>Read Interval—Enter time (in seconds) between reads. The same interval is used for each read mode (for example, the instrument takes images twice as fast if a second read mode is enabled).</p> <p>Note: Time should be no less than exposure time plus read-out time (see Setup Read Mode on page 73 for more information). If it is less, then the experiment speed is determined by the capability of the computer.</p> <p>Number of Reads—Enter the total number of reads to be taken in the first interval.</p> |
| Second Interval | <p>Configure the second series of reads taken during the Read process. To skip a second series with a different read interval, make Number of Reads zero.</p> <p>Read Interval—Enter time (in seconds) between reads. Must be a multiple of the first read time (for example, if read interval 1 is 1 second, then read interval 2 should be 2 seconds or 3 seconds).</p> <p>Note: Time should be no less than exposure time plus read-out time (see Setup Read Mode on page 73 for more information). If it is less, then the experiment speed is determined by the capability of the computer.</p> <p>Number of Reads—Enter the total number of reads to be taken in the second interval.</p> |

Read With TF

Plate reading concurrent with fluid transfer is necessary to collect data for kinetic cell-based assays as the response time is very rapid. To link a Read process to a **Transfer Fluid** process, select **Read** in the **Transfer Fluid** page. Inserting a read this way turns the **Read** process icon purple, indicating that the read occurs concurrently with liquid addition.

Read with TF configuration is similar to that of an independent Read process, but with the addition of more options; **Number of Reads Before Dispense** and **Save Images**.



Note: To remove a **Read With TF** process, deselect the **Read** option in the **Transfer Fluid** process dialog.

Pipettor type: 384

Settings Analysis Transfer Fluid **Read With TF**

First Interval

Read time interval (s):

Number of reads before dispense:

Number of reads after dispense:

Total number of reads:

Save images

Number of images before dispense:

Number of images after dispense:

Second Interval

Read time interval (s):

Number of reads:

The following table describes the setting options in the dialog:

| Options | Description |
|---------------------------------|---|
| Number of Reads Before Dispense | Enter the number of reads to be taken before fluid is added to the plate. This provides a cellular response baseline before the addition of agonist or antagonist. Number of Reads Before Dispense is included in the total number of first interval reads entered in the Number of Reads field, for example, if the Number of Reads is 60 and the number of Reads Before Dispense is five, there are 55 reads after the dispense is initiated. |
| Number of Reads After Dispense | Enter the number of reads to be taken after fluid is added to the plate. This captures the cellular response after the addition of agonist or antagonist. Number of Reads After Dispense is included in the total number of first interval reads entered in the Number of Reads field |
| Save Images | <p>When checked, up to 100 images are saved in an image file (*.tif) associated with an experiment. These images can be used for more than one dispense as long as the total images saved do not exceed 100 for the experiment.</p> <p>Number of Images Before Dispense—Enter the total number of images to be saved prior to initiating the dispense.</p> <p>Number of Images After Dispense—Enter the total number of images to be saved after initiating the dispense.</p> <p>By default one image prior to and nineteen after dispense are saved for review in Image Display on page 104.</p> <p>Note: Images may be useful to troubleshoot problems with your cell plate such as cells detaching during fluid dispense.</p> |

Wash Tips or Pins Process

This process is normally used to eliminate unwanted agonist or antagonists prior to another addition within an experiment or between experiments. In many cases, it reduces consumable costs and downtime associated with replacing tips.

For standard pipettors, tips are inserted into the wash location and wash fluid aspirated and dispensed. This can be done a number of times (strokes) in the same wash fluid, then the whole process repeated up to five times (wash cycle) with fresh wash fluid.

For the pin tool, up and down motion is used to force fluid in and out of the capillary slot of the pins. A typical wash protocol uses an initial solvent to remove compound from the pins, followed by a second solvent to remove the first solvent. It is recommended that pins are blotted at a blotting station to remove excess liquid from the pins between washes.

Wash Tips (Standard Pipettor)

Pipettor type: 384

Settings
Analysis
Transfer Fluid
Read With TF
Wash Tips

Wash Tips

Fluid type: A ▼
 Pump Speed: Fast ▼

Wash Cycles: 2 ▼
 Strokes: 5 ▼

Volume/Stroke (µL): 28.0
 Hold Time (s): 0.0

Aspirate Speed (µL/s): 20.00 ▼
 Dispense Speed (µL/s): 20.0

Expel Volume (µL): 2.0 ▼

Unload Tips After Washing

Table 3-32: Wash Tips Configuration Options for Standard Pipettor Use

| Options | Description |
|----------------|--|
| Fluid Type | Select the fluid in the wash reservoir during tip washing, from wash fluid container A or B. |
| Pump Speed | Select the speed (slow or fast) you want the pump to fill and empty the wash reservoir. Note: Molecular Devices recommends you use Fast except with volatile solvents. |
| Wash Cycles | From the list, select the number of times that fluid cycles through the wash reservoir. For each cycle, the pipettor executes the number in the Strokes field. |
| Strokes | From the list, select the number of times to aspirate and dispense wash fluid. |
| Volume/Stroke | Enter the volume (in µL) of fluid to be aspirated and then expelled for each stroke. |
| Hold Time | Enter the length of time (in seconds) to pause between aspirating and expelling fluid. |
| Aspirate Speed | From the list, select the speed (in µL/s) to aspirate fluid. |

Table 3-32: Wash Tips Configuration Options for Standard Pipettor Use (continued)

| Options | Description |
|---------------------------|--|
| Dispense Speed | From the list, select the speed (in $\mu\text{L/s}$) to dispense fluid. |
| Unload Tips After Washing | Check this option to have tips removed from the pipettor head after washing is complete. |

Wash Pins (Pin Tool)

Pipettor type: 384pintool

Settings Analysis Transfer Fluid Read With TF Wash Pins

Wash Pins

Fluid type: A ▼ Pump Speed: Fast ▼

Wash Cycles: 2 ▼ Strokes: 5 ▼

Down Speed (mm/s): 10.0 Stroke Distance (mm): 5.0

Up Speed (mm/s): 10.0 Removal Speed (mm/s): 10.0

Pause In Well (s): 0.0 Unload Pin Tool After Wash

Table 3-33: Wash Pins Configuration Option for Pin Tool User

| Options | Description |
|-------------|--|
| Fluid Type | Select the fluid in the wash reservoir during tip washing, from wash fluid container A or B. |
| Pump Speed | Select the speed (slow or fast) you want the pump to fill and empty the wash reservoir. Note: Molecular Devices recommends you use Fast except with volatile solvents. |
| Wash Cycles | From the list, select the number of times that fluid cycles through the wash reservoir (up to 5). For each cycle, the pipettor executes the number in the Strokes field. |
| Strokes | From the list, select the number of times to raise and lower the pins. One stroke is equal to one upstroke and one down-stroke. |
| Down Speed | Set the speed of the pin tool during the downstroke. |

Table 3-33: Wash Pins Configuration Option for Pin Tool User (continued)

| Options | Description |
|----------------------------|--|
| Stroke Distance | Set the height of one stroke that the pin tool travels. |
| Up Speed | Set the speed of the pin tool during the up-stroke. |
| Removal Speed | Set the speed that the pin tool is withdrawn from the wash location. The faster the withdrawal speed the more fluid remains on the pins. |
| Pause in Well | Set a time to hold the pin tool at the bottom of the wells in the wash solution after the wash is completed. |
| Unload Pin Tool After Wash | Check to have the pin tool removed after washing is complete. |

Blot Pins Process

The **Blot Pins** process is only available for the pin tool. This function is used to dry the pins between transfers or washes, to prevent carry-over.

Pipettor type: 384pintool

Settings
Analysis
Transfer Fluid
Read With TF
Wash Pins
Blot Pins

Wash Pins

Fluid type: A ▼ Pump Speed: Fast ▼

Wash Cycles: 2 ▼ Strokes: 5 ▼

Down Speed (mm/s): 10.0 Stroke Distance (mm): 5.0

Up Speed (mm/s): 10.0 Removal Speed (mm/s): 10.0

Pause In Well (s): 0.0 Unload Pin Tool After Wash

Table 3-34: Blot Pins Configuration Options

| Options | Description |
|----------------------------|---|
| Target Plate | Select the plate. |
| Height | <p>Enter distance (measured in μL) from the bottom of the well to place the pins. The distance equivalent in millimeters is displayed as well.</p> <p>Float pins—Moves the pipettor down to a position at which the pins would be 1 mm below the bottom of the wells. As pins come to rest on the blotting surface additional downward motion of the tip block is absorbed by the tips moving into the block. This ensures that all pins are in contact with the blotting surface. The Height is designated as -1.00 mm.</p> |
| Pause in Well | Enter a duration during which the pin tool is held motionless at the set height. |
| Unload Pin Tool After Blot | Automatically unload the pin tool once blotting is complete. |

Pause Pipettor Process

Pause Pipettor steps are included in a protocol to prevent future pipettor motion until the pause is complete. **Read** processes can continue to execute, unless stopped by the linked action “read with transfer”. This can be useful to stop the pipettor from picking up compound if it sits for an extended period in the air. Also, it can be used to provide time for solvents (such as EtOH) to evaporate off the pins after a wash.

This process has only one configurable setting, the duration of the pause, entered in seconds.

Wash Cell Reservoir Process

This process is only available when the Cell Reservoir is installed and is recommended for use before and after using cells in suspension.

This process can be used as a way of priming the reservoir with cells prior to an assay or for cleaning purposes.

No pipetting is associated with this step.

The following table describes the options available in the **Wash Cell Reservoir** process page.

| Options | Description |
|--|---|
| Pre-coat Tubes | Select to pump the fluid up to the sensor before filling the reservoir. |
| Fluid Source | Select which fluid to use for the wash process. All Fluid 1–4 bottles are available as well as the Cell Flask. |
| Drain Reservoir Destination ¹ | Select a drain destination for the fluid left in the reservoir after the wash. Waste bottles, Cell Flask, or any of the Fluid bottles are available. |
| Fill Reservoir Speed ¹ | Select a pump rate for filling the Cell Reservoir; the range is 1–10. |
| Drain Reservoir Speed ¹ | Select a pump rate for draining the Cell Reservoir; the range is 1–10. |
| Wash Cycles | Set how many wash cycles with the above settings should be done. |
| Hold Time | This option soaks the reservoir in the pumped fluid by leaving the fluid in the reservoir for a specified time; the range is 0–300 seconds (0–5 minutes). |
| 1. This option is only available when the Cell Reservoir is installed and selected as the source plate for aspiration. | |

Finish With Source Process

In **Remote Mode**, the **Finish with Source** process instructs the instrument to notify the robot controller to remove the desired source plate. This option is required when using multiple plates in a single source location.



Note: This process is not active in **Manual Mode** or for the read plate position.

Creating New Protocols

In the ScreenWorks Software, creating a protocol involves combining processes represented in the protocol builder as colored block icons at the top of the Experiment window. The available processes depend on your instrument configuration. See [Instrument Configuration Tab on page 69](#). The color of the process icon depend on the function in the protocol. See [Protocol Process Icon Colors on page 71](#).

A red box around the process in the protocol list indicates the process selection and indicates which configuration settings dialog is open.

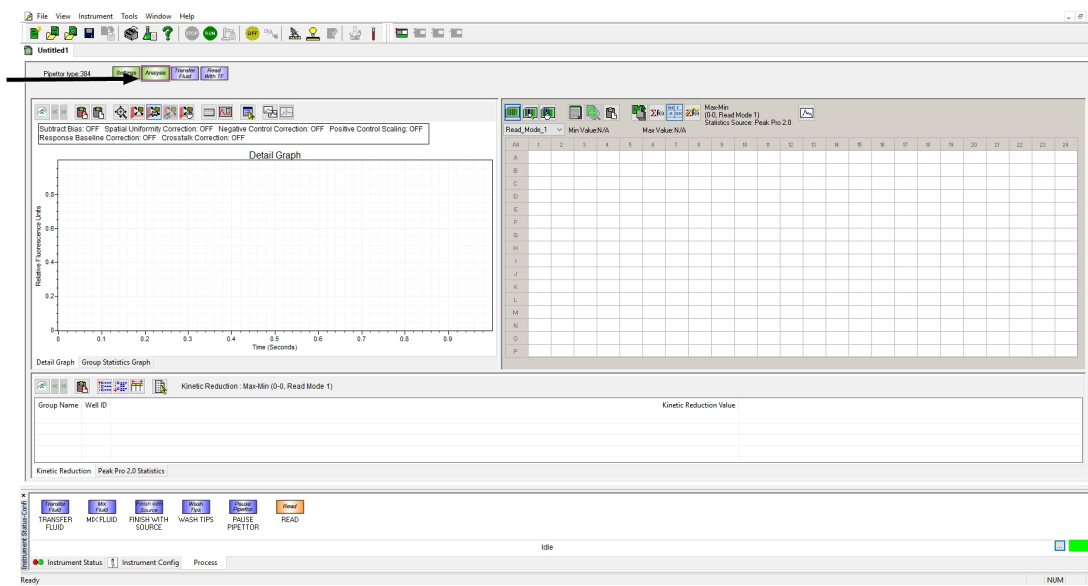


Figure 3-9: Example New Protocol Build

Protocols contain combinations of the following processes. Available options depend on the components installed in your instrument.

- **Settings**—required for every protocol, automatically included at the beginning of every new protocol file, and cannot be removed.
- **Analysis**—required for every protocol, automatically included at the beginning of every new protocol file, and cannot be removed.
- **Transfer Fluid**—automatically included in new protocols and have the following selectable linked processes:
 - **Read with TF**—automatically included in new protocols after Transfer Fluid, and timed precisely to coincide with compound addition to the read plate. This setting can be deselected from within the Transfer Fluid configuration settings dialog.
 - **Mix with TF**—optional selection within the Transfer Fluid configuration settings dialog.
- **Mix Fluid**
- **Wash Reservoir**—only available when the **Cell Suspension** option is installed.
- **Wash Tips**—available only with pipettor configuration.
- **Wash Pins**—available only with pintool configuration.
- **Blot Pins**—available only with pintool configuration.
- **Pause Pipettor**
- **Finish with Source**
- **Read**

To create a new protocol:

1. Open the **Process** tab panel. See [Process Tab on page 70](#).
2. In the **Process** tab panel, add the assays steps. Do one of the following:
 - To place the new process at the end of your developing protocol list, drag and drop the process anywhere in the Experiment window.
 - To place the new process somewhere specific in your developing protocol list, drag and drop the process where you want it. You are notified when the position is invalid.



3. When the configuration screen opens for the added process, adjust settings as needed. See [Understanding Processes on page 72](#).
4. Repeat steps 2 and 3 until you are finished building your protocol.
5. Define the **Analysis** parameters.
Click the **Analysis** process, and then click **Grouping**, **Correction**, and **Export** to specify the analysis parameters for the experiment data. See [Analysis Process on page 78](#).
6. Click on any other process icon to open its configuration screen and make changes as needed.



Note: When using the HS EMCCD camera, depending on your **Protocol > Settings > Edit read mode > Camera Mode** setting, when using a fast Read time interval value, the plate matrix panel updating might be delayed until the end of the data acquisition rather than updating during the data acquisition.

Specifically, the following settings result in the plate matrix panel updating at the end of the data acquisition:

For **Camera Mode > HighSpeed**—a **Read time interval(s)** setting of 0.018 seconds or shorter

For **Camera Mode > Normal or Sensitivity**—a **Read time interval(s)** setting of 0.029 or shorter

7. Select **File > Save**.
8. In the **Save As** dialog, in the **File name** field, type a name for your protocol, and then click **Save**.



Tip: Protocol files are saved as an ***.fmp** file. Unless you change the **Save in** location for your protocols, the default save location is **C:\Documents\Molecular Devices\ScreenWorks\MyProtocols**.

9. Run your experiment.

Refer to the *FLIPR Penta High-Throughput Cellular Screening System Protocol Guide* for specific protocol settings.

Deleting Processes from the Protocol

Protocol processes are either linked to the Transfer Fluid process or not linked.

To delete processes that are not linked:

- In the protocol list, select a non-purple process icon, and then on your keyboard, press **DELETE**.

To delete a Transfer Fluid linked process:

1. In the protocol list, select the purple **Transfer Fluid** icon.
2. In the configuration settings dialog, deselect the linked process you want deleted, either **Read** or **Mix Fluid**.

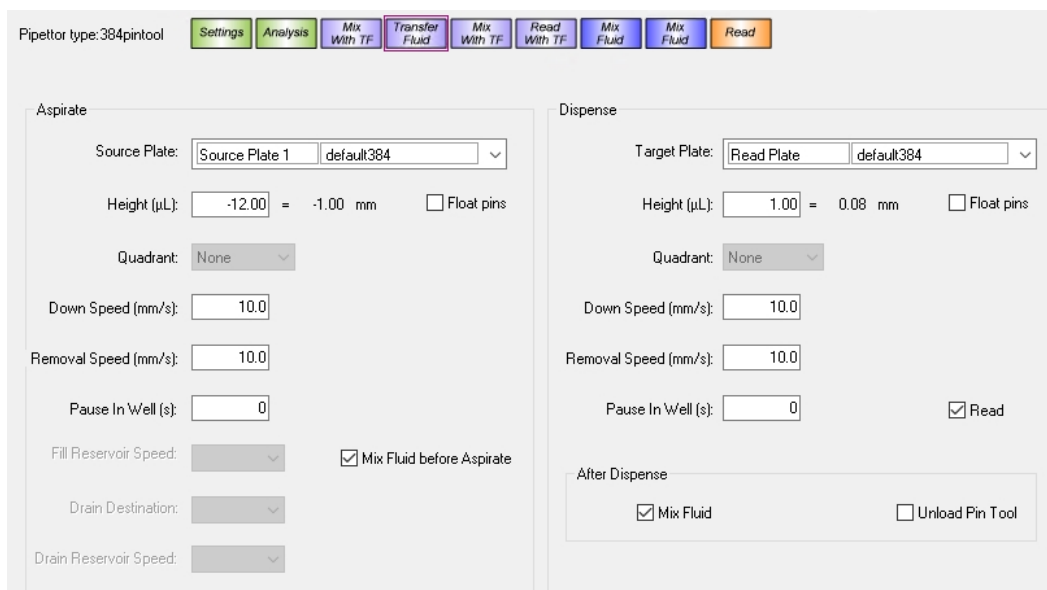


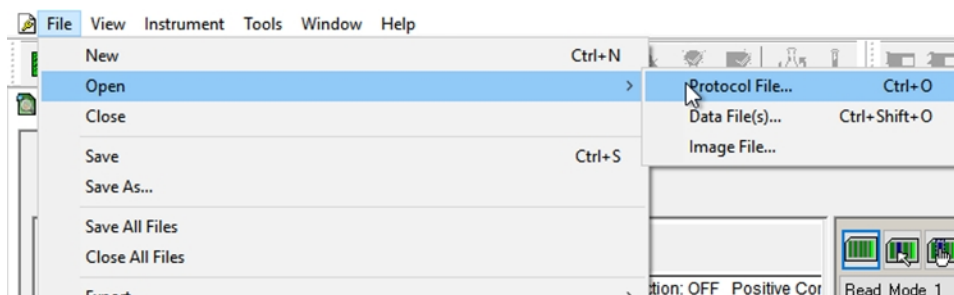
Figure 3-10: Transfer Fluid Configuration Settings Dialog With Linked Processes

Opening Protocols

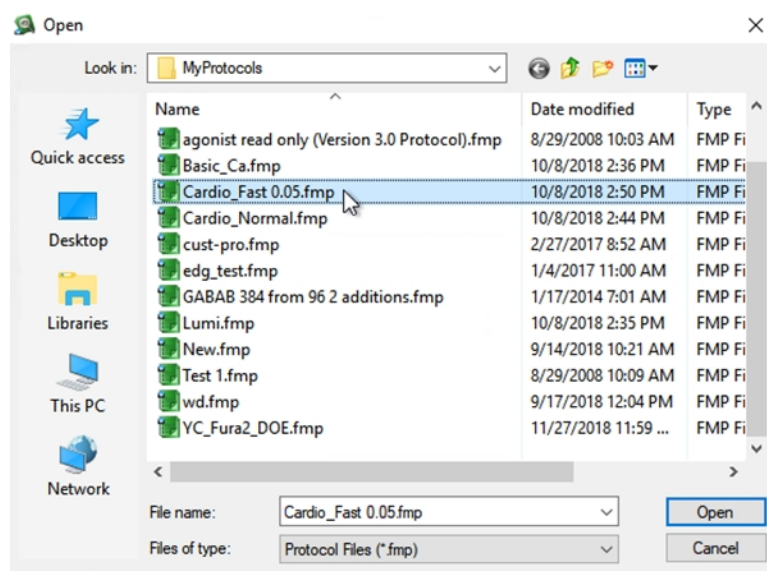
If you have saved protocol files to use, you can open an existing protocol file, rather than create a new one in the ScreenWorks Software.

To open a protocol file:

1. Select **File > Open > Protocol File**.



- In the **Open** dialog, select your protocol file name, then click **Open**.



Deleting Protocols

You cannot delete a protocol from within ScreenWorks Software.

To delete a protocol:

- In your Windows File Explorer navigate to where the protocols are saved.

*** Tip:** During the creation process, unless you changed the **Save in** location for your protocols, the default location for all of your protocol files is **C:\Documents\Molecular Devices\ScreenWorks\MyProtocols**.

- Select the ***.fmp** file you want removed, and then on your keyboard, press **DELETE**.

Saving Data Files as Protocol Files

Protocol information stored in data files (*.fmd) cannot be edited, nor used to run a new experiment; however, you can extract this information to a new protocol file. Experiments run using this new protocol file have exactly the same protocol steps as used to create the data file.

*** Tip:** When using a protocol file (*.fmp), you can add, remove, or change processes in the file, then save the amended protocol and run it.

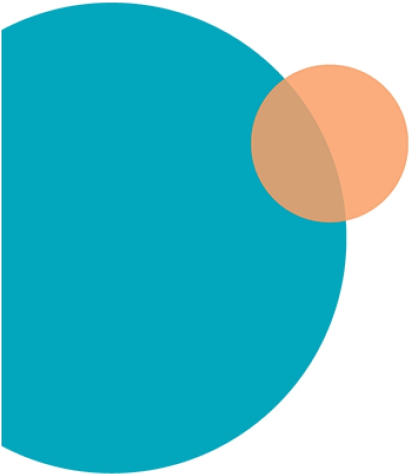
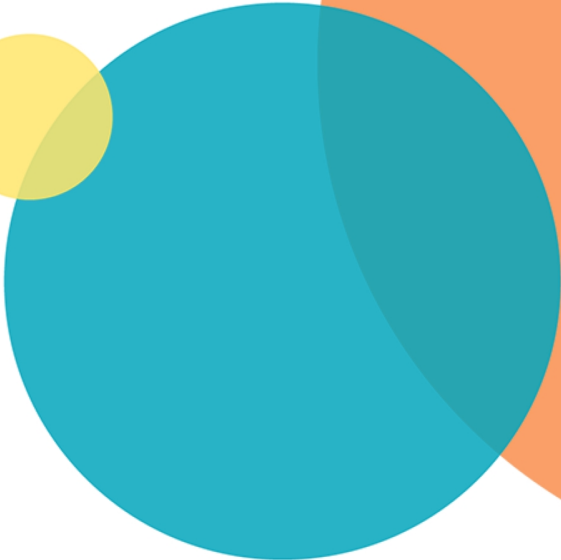
To store a data file as a protocol file:

- Select **File > Save As**.
- In the **File name** field, type a protocol name.
- In the **Save as type** field, select **Protocol Files (*.fmp)** option and click **Save**.

The stored file is stripped of all data and only associated protocol information is stored in the protocol file.



Note: Saved changes only affect the protocol. The data file from which the protocol was derived remains intact.



Chapter 4: Calibrating and Signal Testing

The following procedures and information are required to calibrate the optics and run a signal test of the FLIPR Penta System:

- [Calibrating Optics](#)
- [Running Signal Tests on page 135](#)

Calibrating Optics

Calibrating the optics of your FLIPR Penta System requires taking an image of the 96-well, 384-well, or 1536-well plate and using it as a mask to locate the wells. Typically, calibrations are run with the following procedures:

- After exchanging LED banks. See [Exchanging LED Modules on page 198](#).
- After exchanging the pipettor or pin tool head. See [Installing the Pipettor or Pin Tool Head on page 190](#)
- After a new plate is added to the plate library. See [Adding a Read Plate to Plate Library](#).
- After a new filter is added to the system. See [Exchanging Filters on page 201](#).

Adding a Read Plate to Plate Library

The FLIPR Penta System has default plate masks for 96-well, 384-well, and 1536-well read plate types. If you want to use a plate type that varies significantly from the default formats, you must create new dimensions and mask for that plate type. Use the description of the Plate Library menu item, see [Plate Process Definition on page 59](#), as a reference guide for options and parameters available when adding a plate to the software.

To add a plate and mask to the software:

1. Prepare a 1×10^{-8} M fluorescein solution.
2. Ensure the microplate bottom is clean and free of scratches and dust.
3. Dispense enough fluorescein solution into each well so as to cover the bottom.
Dispensing half of the total well volume is typically a safe volume with which to work. Use the FLIPR Penta System pipettor.
4. Place the microplate containing fluorescein in the read position and select **Plate Library** from the **Tools** menu.
5. From the **Plate Process Definition** list, select a plate with a similar well format.
6. Click **Copy Plate**.
7. Enter the **Plate Specifications**:
 - Enter a **Plate Name**.
 - Select **Read Plate** as the **Plate Type**.
 - Type in the plate dimensions.
8. Click **Finish** to create a new read plate with mask.

The plate mask definition is complete when the green **Assay Finished (Unlocked)** light illuminates on the instrument door.

Calibrating the Optics

FLIPR Penta System identifies the serial number, wavelength range and position of each LED bank present in the instrument. This information, in addition to the wavelength range for the emission filters, is stored with the system calibration files. Any time you change LEDs or emission filters in the FLIPR Penta System, you must recalibrate the Flat-Field Calibration files for the new LED/Filter pairs. The following procedure steps you through the process of calibrating the system optics.

*** Tip:** Due to the improved light tightness of the new FLIPR Penta System, the background correction calibration available in the previous versions of this software used with black bottom plates is no longer required.

If you have any questions or concerns regarding the calibration procedure, contact Molecular Devices Technical Support. See [Obtaining Support on page 226](#).

Excitation LED

The **Excitation LED Calibration** is needed only if the **LED Calibration Status** shows **NotDone** when you select to use the **335–345 nm/475–535 nm** (340 LED Module).

To run an Excitation LED Calibration:

1. Select **Instrument > Calibration**, or click on the **Calibration** icon.
The **Calibration** dialog displays a list of the **Excitation/Emission Wavelength** (LED/Filter) combinations for the installed modules, and corresponding status, **LED Calibration** either **Done** or **NotDone**, and **Flat-Field Calibration** either **Done** or **NotDone**.
2. Select **335–345 nm/475–535 nm**.
3. Click **Excitation LED Calibration**.
4. When the progress bar dialog closes, continue with **Flat-Field Calibration**. See [Flat-Field Calibration on page 132](#).

Flat-Field Calibration

Before adjusting your exposure settings, you first need to have an accurate flat-field calibration. This calibration is used to correct for light falloff at the corners of the camera lens, and to correct for any gradients or falloff in the coverage of the light module. When an accurate flat-field correction is applied, the well values for all wells give the same count value for the same stimulus.

*** Tip:** When testing for HS EMCCD saturation in the Signal Test dialog, keep this flat-field correction in mind, because the central values hit saturation first, while the outer values continue to get larger as the exposure increases to beyond the maximum camera sensor limit, because they are being multiplied by the correction factor.

During a fluorescence assay, microplate illumination results in imperfect, non-uniform excitation distribution on the microplate surface. In the FLIPR Penta System, light distribution is more intense in the middle of the microplate than the edges, therefore a flat-fielding algorithm is applied to the microplate image to compensate for variations in light intensity across the microplate.

During luminescence assays, LEDs are not used. Flat-field corrections are preset during installation and do not require any adjustment.

To run Flat-Field Calibration:

1. Select the **Instrument > Calibration**, or click on the **Calibration** icon.
The **Calibration** dialog displays a list of the **Excitation/Emission Wavelength** (LED/Filter) combinations for the installed modules, and corresponding calibration status for **Flat-Field Calibration**, either **Done** or **NotDone**, and **LED Calibration** of the selected **Plate Format**.



Note: **None** indicates either no LEDs are illuminated or the instrument used a blank emission filter position to create the calibration file.

2. Select a **Plate Format** (96, 384, or 1536).
3. Select the **Excitation/Emission Wavelength** pair with the **NotDone** status from the **Calibration Status** list.
4. Place the **Flat-Field Calibration Plate** into the **Read Position**.

The provided Yellow Plate can be used to Flat-Field calibrate the following **Excitation/Emission Wavelength** pairs:

- 335–345 nm/475–535 nm
- 380–390 nm/475–535 nm
- 390–420 nm/565–625 nm
- 420–455 nm/475–505 nm
- 420–455 nm/515–575 nm
- 470–495 nm/515–575 nm
- 470–495 nm/565–625 nm
- 495–505 nm/526–586 nm
- 510–545 nm/565–625 nm
- 610–626 nm/646–706 nm

The following dye should be used to calibrate the respective **Excitation/Emission Wavelength** (LED/Filter) pair:

- 360–380 nm/400–460 nm with MQAE. See [MQAE on page 134](#).
- 390–420 nm/440–480 nm with Coumarin. See [Coumarin on page 134](#).

5. Click **Flat-Field Calibration**.
6. After the **Assay Finished (Unlock)** light illuminates on the instrument, and the **Excitation/Emission Wavelength** status shows **Done** in the **Flat-Field Calibration** column, click **Close** to exit.
7. Place your yellow signal test plate of the appropriate well format into your instrument.

8. Run a signal test by selecting **Instrument > Manual Operation > Yellow Plate Signal Test**, or using the corresponding  button. The yellow plate standard deviation should be 5% or less. See [Running the Yellow Plate Signal Test on page 136](#).

Coumarin

Required material:

- Black clear-bottom plate of needed format and brand for assay (96, 384, or 1536)
- Coumarin (7-diethylaminocoumarin-3-carboxylic acid)
Catalog # 36799-100MG-F, Sigma-Aldrich

To run a Coumarin calibration:

1. Make a 0.02 M stock solution of Coumarin in Dimethyl Sulfoxide (DMSO). Mix by vortexing tube.
2. Store small aliquots at -20°C (-4°F).
3. Within one hour of calibration, thaw an aliquot and make 20 mL 10⁻⁵ M solution in the same buffer to be used for dye loading. Adjust pH to 7.4. Mix by vortexing.
4. Pipette a uniform quantity of the Coumarin solution into each well of the microplate. The volume must be adequate to cover the bottom of the well uniformly. The following quantities are recommended:
 - 96-well plate, 100 µL/well
 - 384-well plate, 40 µL/well
 - 1536-well plate, 6 µL/well
5. Check the plate visually to make sure there are no bubbles or unfilled wells. Shake or tap the microplate to dislodge any bubbles. Keep the Coumarin plate covered and in the dark until used.
6. Use the Coumarin plate as the **Flat Field Calibration Plate** in [Flat-Field Calibration on page 132](#).

MQAE

Required material:

- Black clear-bottom plate of desired format and brand for assay (96, 384, or 1536)
- MQAE (N-(ethoxycarbonylmethyl)-6-methoxyquinolinium bromide)
Catalog # 46123-100MG-F, Sigma Aldrich

To run a MQAE calibration:

1. Make a 0.02 M stock solution of MQAE in distilled water. Mix by vortexing tube.
2. Store small aliquots at -20°C (-4°F).
3. Within one hour of calibration, thaw an aliquot and make 20 mL 10⁻⁵ M solution in water or in a chloride-free buffer to be used for dye loading. Adjust pH to 7.4. Mix by vortexing.

4. Pipette a uniform quantity of the MQAE solution into each well of the microplate. The volume must be adequate to cover the bottom of the well uniformly. The following quantities are recommended:
 - 96-well plate, 100 μL /well
 - 384-well plate, 40 μL /well
 - 1536-well plate, 6 μL /well



Note: Check the plate visually to make sure there are no bubbles or unfilled wells. Shake or tap the microplate to dislodge any bubbles. Keep the MQAE plate covered and in the dark until used.

5. Use the MQAE plate as the **Flat Field Calibration Plate** in [Flat-Field Calibration on page 132](#).

Running Signal Tests

Run the following signal tests, as specified:

- Once a day, to check the system optics calibration, run the **Yellow Plate Test**. See [Running the Yellow Plate Signal Test](#).
- Before running an experiment, to verify your protocol setting performance, run the **Protocol Signal Test**. See [Running a Protocol Signal Test on page 138](#).




Tip: For the HS EMCCD camera using the HighSpeed or the Sensitivity camera mode, run a Protocol Signal Test to determine your saturation level for your biology and adjust your exposure settings accordingly. This is important because, after a 4 times overexposure increase, there are no oversaturation warning messages from this camera.

Running the Yellow Plate Signal Test

Use the Yellow Plate Signal Test once a day to verify that the system optics are calibrated. Additionally use the Yellow Plate Signal Test after you recalibrate or change the optics.

To run a Yellow Plate Signal Test:

1. Place the yellow signal test plate on the stage in the read position.
2. In the ScreenWorks Software, select **Instrument > Manual Operation > Yellow Plate Signal Test** or click  **Yellow Plate Signal Test**.

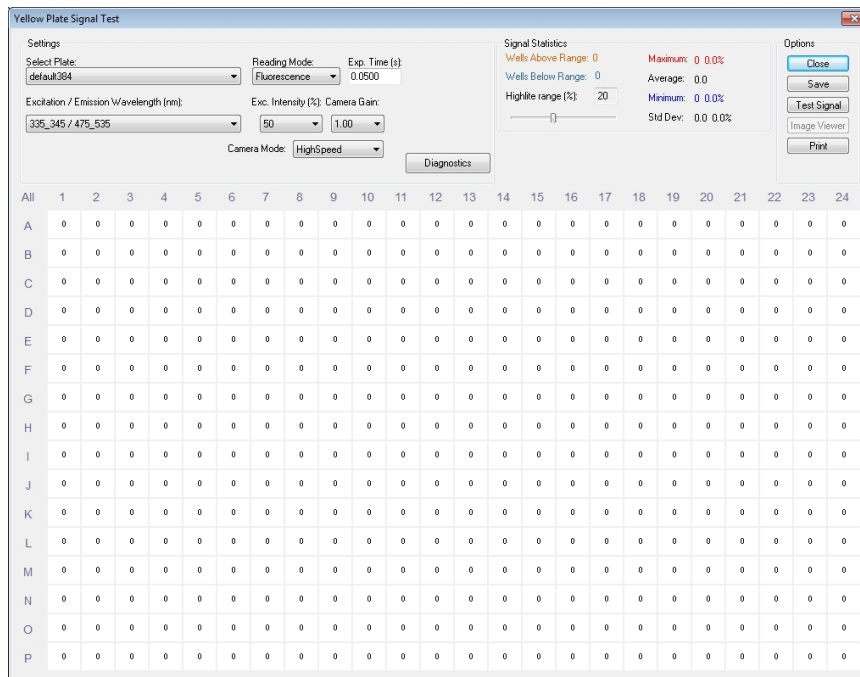


Figure 4-1: Yellow Plate Signal Test Dialog

3. Set the following parameters:
 - a. Select the correct plate from the **Select Plate** list.
 - b. Select **Reading Mode > Fluorescence** only.



CAUTION! Although both Fluorescence and Luminescence are available options in the test dialog, currently there is no easy way of testing signal strength in Reading Mode > Luminescence. Do not run any Reading Mode > Luminescence signal tests.

- c. Select the appropriate **Excitation/Emission Wavelengths**.
 - d. Set the **Excitation Intensity** to 50.
 - e. Set the **Exposure Time** to 0.1 seconds.
 - f. Adjust the following variable camera settings:
 - For the EMCCD camera, set the **Camera Gain** to 80.
 - For the ICCD camera, set the **Gate Factor** to 6%.
 - For the HS EMCCD camera:
 - In **Camera Mode > Normal**, set the **Camera Gain** to 1.5 and the **Exposure Time** to 0.005 seconds.
 - In **Camera Mode > HighSpeed**, set the **Camera Gain** to 20 and the **Exposure Time** to 0.001 seconds.
4. Take a picture by clicking **Test Signal**.

When the instrument is calibrated with the yellow test plate for the appropriate plate format and optics, normal test plate results are a relative standard deviation less than 5%.

If the test results are abnormal values, look at the **Image Display** for any anomalies on the plate. For additional information see, [Troubleshooting on page 209](#).

5. Print the results and keep them in a “Maintenance” folder by the instrument to track the standard deviation of the yellow signal test plate over time.



Note: The **Yellow Plate Signal Test** and **Image Display** results are not saved within a data file.

Alternatively, you can save the files on the hard drive in the signal test directory (**C:\Documents and Settings\[your_user_name]\My Documents\Molecular Devices\ScreenWorks\MySignalTests**) as a *.sig file, which can be opened in third-party spreadsheet software.



Note: The relative standard deviation should be less than 5% if the flat-field calibration was performed using the Flat Field Calibration Plate of the respective plate format (96-wells, 384-wells, or 1536-wells). See [Flat-Field Calibration on page 132](#).

Running a Protocol Signal Test

Use the Protocol Signal Test to optimize the camera settings used in your fluorescence assay protocols. Adjustments made to the settings in this test can be saved within your protocol.

Refer as needed to the fluorescence signal recommendations in [Optimizing Optics Hardware on page 151](#).



CAUTION! Although both Fluorescence and Luminescence are available options in the test dialog, currently there is no easy way of testing signal strength in Reading Mode > Luminescence.

Do not run any Reading Mode > Luminescence signal tests.

To run a fluorescence protocol signal test:

1. Place your cell plate on the stage in the read position.
2. In the ScreenWorks Software, open a fluorescence assay protocol you want to test.



3. Click **Protocol Signal Test**.
4. Click **Test Signal**.
5. Review the results and change settings as needed in the **Protocol Signal Test** dialog.

Depending on the installed camera and the read mode, different parameters are available for adjusting the signal strength. See [Signal Test Dialog Options on page 55](#) for more details.

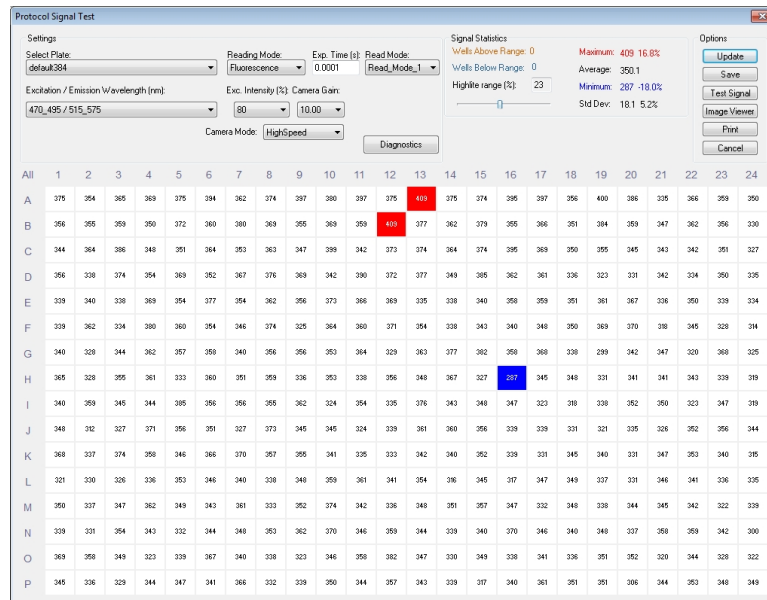


Figure 4-2: Example Protocol Signal Test Dialog

- Depending on the range of values acquired in your results, decide either to run the experiment or change the **Excitation Intensity**, **Exposure Time**, **Camera Gain**, or **Camera Gate** settings.

*** Tip:** For a calcium mobilization assay, Molecular Devices recommends starting between 200 to 1,500 counts.



CAUTION! When using the HS EMCCD camera, after a 4 times overexposure increase, there are no warning messages.

- Click **Update** to apply the new settings, and click **Test Signal** to recheck them.
- To keep the new settings as part of your open protocol, click **Save**.

Determining Protocol Saturation Levels for the HS EMCCD Camera

When using the HS EMCCD camera HighSpeed camera mode and Sensitivity camera mode, after about a 4x overexposure level increase, no saturation warning appears. This is a problem for detecting when the image measurements are saturated, and the system cannot give a reliable saturation warning for grossly overexposed well signals. For more details, see [Understanding HS EMCCD Camera Protocol Saturation on page 38](#).

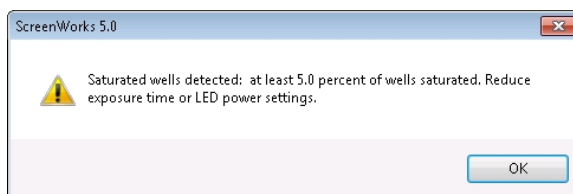
While you create a protocol that uses the HS EMCCD camera HighSpeed camera mode or Sensitivity camera mode, run a Protocol Signal Test using a cells test plate to determine your optimal exposure levels.

To determine the optimal exposures for your biology using HighSpeed and Sensitivity camera mode:

- Prepare a test plate to check for saturation of your maximum signals with the **Protocol Signal Test**. See [Running a Protocol Signal Test on page 138](#).
- Start with an initial estimated exposure setting, based on experience or as recommended in the *FLIPR Penta High-Throughput Cellular Screening System Protocol Guide*, then adjust your exposure to keep your maximum signals below saturation.

*** Tip:** The gain factors are approximately linear, so reducing the gain by half is similar to reducing exposure time by half.

- Click **Test Signal** to see the exposure setting results.
- If the test gives you a saturation warning, reduce your exposure by a factor of 4x and test again.



5. If the test does not give a saturation warning, click **Image Display**, and do the following:
 - See if the image is not showing bright wells, which confirms that you are below saturation.

*** Tip:** If you are not sure, reduce your exposure by 2x and test again to verify.

- See if the image is blooming. If yes, reduce your exposure by a factor of 8x and test again.

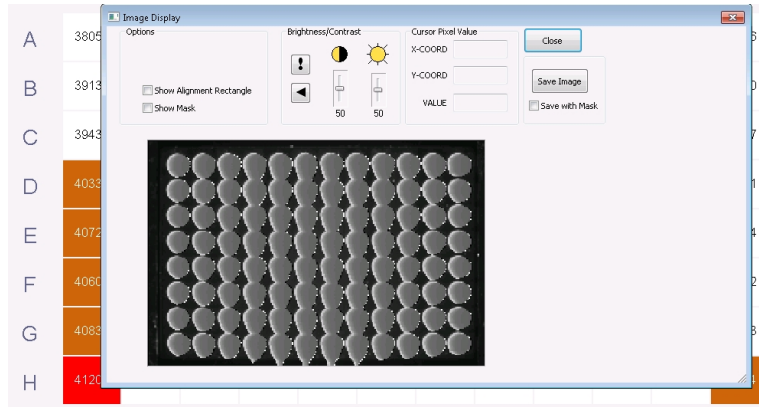


Figure 4-3: HighSpeed Camera Mode Blooming Image Display

- See if the image is reversing, showing wells with dark centers and bright outlines. If yes, reduce your exposure by a factor of 4x and test again.

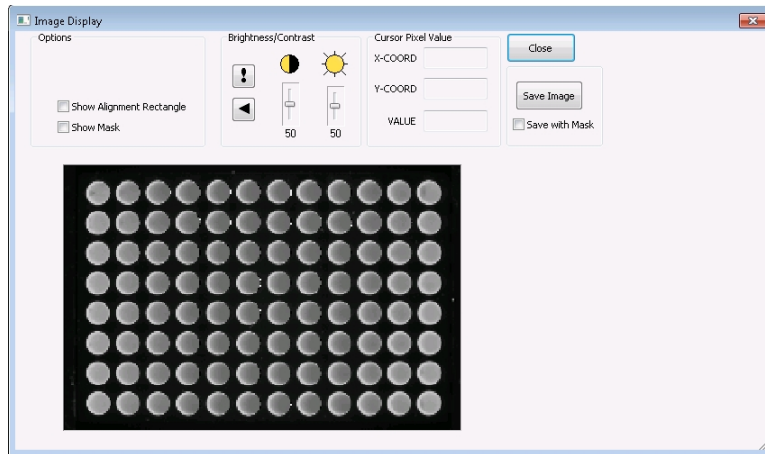



Figure 4-4: HighSpeed Camera Mode Reversing Image Display

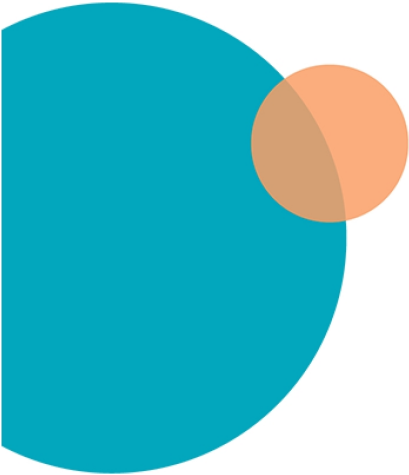
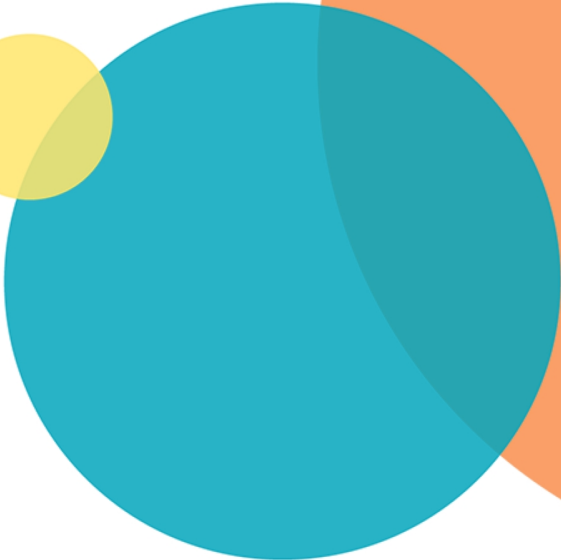
6. When you have a test which is below the saturation level, adjust your exposure linearly to give an average value that matches your target exposure. Keep the maximum signal below 43,000 counts.

*** Tip:** If your experiment could have brightness that exceeds your signal test sample plate, then reduce the exposure accordingly so that your brightest signal does not exceed 43,000 counts.

7. Leave enough room above the starting levels so that you can measure your experiment without camera saturation.

 **Tip:** If you do not expect your signal to drop during the experiment, start as low as you can to give enough room for the signal to rise without saturation.

8. To keep the new settings as part of an open protocol, click **Save**.



The following provides a starting point for setting up and running a kinetic cell-based assay on the FLIPR Penta System. Refer to the *FLIPR Penta High Throughput Cellular Screening System Protocol Guide* for specific assay protocols.

To run your experiment, use the following workflow processes:

1. [Preparing Cells](#)
2. [Starting the System on page 145](#)
3. [Checking the System on page 146](#)
4. [Dye-Loading the Cells for Fluorescence Assays on page 146](#)
5. [Preparing a Source and Compound Plate on page 147](#)
6. [Setting Up an Assay Protocol on page 149](#)
7. [Optimizing the Optics and Fluid Dispensing on page 150](#)
8. [Starting the Assay Run on page 155](#)

Preparing Cells

Prepare your cells according to the following assay you need to run:

- [Preparing Cells for Adherent Assays](#)
- [Preparing Cells for Suspension Assays on page 144](#)

Preparing Cells for Adherent Assays

Location of Cells in the Plate

The excellent signal-to-background noise ratio of the FLIPR Penta System is made possible by the instrument's bottom illumination and detection system. The read plate is illuminated by the LED modules at an angle approximately 45°. It is recommended to have cells located at the bottom of the wells during an experiment. Adherent cells are typically grown overnight in the cell plate, whereas non-adherent cells are seeded either the night before, on a coated plate, or the day of an experiment and then centrifuged.

Cell Densities

Cell densities used in fluorescence assays vary because each cell type has different requirements. Cell densities range from 1,500 to 5,000 cells per well for 1536-well plates, 5,000 to 20,000 for 384-well plates and 20,000 to 80,000 cells per well for 96-well plates. Non-adherent assays require higher cell densities.

It is necessary to optimize the cell seeding density so that a uniform, 80–90% confluent monolayer is formed on the day of the experiment. Over- or under-confluent cell monolayer may result in reduced cellular response to the test compounds.

Cells that are normally maintained in culture at subconfluent levels should be seeded at relatively low densities. Depending on the individual cell line, attachment matrix-coated plates (such as the poly D-lysine coated plate) might be required to improve adherence and minimize cellular “blow-off” during compound addition. Assay development needs to be performed to determine the optimal seeding density and plate environment for each cell line.

Cell Seeding

Cells can be seeded in plates using a multi-channel pipettor or a liquid dispensing system. We recommend seeding a 1536-well or 384-well cell plate using an automatic instrument rather than seeding manually. Thin needles used in automatic liquid-dispensing instruments prevent air bubble formation in well bottoms—a problem commonly encountered when cells are seeded with a manual pipettor. 96-well plates can be seeded by either manual or automatic methods. Refer to your specific instrument user guide for instructions on how to dispense cell suspension into wells.

Cells are seeded in clear, flat-bottom, black-wall 1536-well, 384-well, or 96-well tissue culture treated plates. A flat plate bottom ensures that cellular fluorescence is localized to a single horizontal plane. Adherent cells typically are seeded one day before an experiment. Non-adherent cells are either plated one day before on a coated plate or the day of an experiment. All steps are carried out in the same black-wall 96-well, 384-well, or 1536-well plate.



Note: Flat-bottom, clear-wall tissue culture plates may also be used with the FLIPR[®] Calcium Assay Kit and Membrane Potential Assay Kit. Depending on signal intensity, either black-wall or white-wall plates can be used when running aequorin assays.

Preparing Cells for Suspension Assays

Location of Cells in the Plate

Suspension cells can be used for luminescence experiments. In this case, cells are kept in uniform suspension by the Cell Suspension option, which consists of a cell suspension module and reservoir system. Suspended cells are first pumped to the reservoir and then delivered to the read plate by the FLIPR Penta System pipettor head.

These assay methods are also described in the literature, for example, in Boie et al., *Eur. J. Pharmacol.*, 340(2-3):227-241 (1997), and in United States Patent 6,872,536 and European Patent 1,145,002. Users interested in the patented methods may wish to consult legal counsel in evaluating these patents.

Cell Densities

Cell densities in luminescence suspension assays vary. Depending on the bioluminescence signal from the cells, the density can be adjusted to reach appropriate signal window. Bright cells can be diluted, while dim cells might require a higher concentration. The suggested cell densities are 1,000–5,000 cells per well for 1536-well plates, 2,500–10,000 cells per well for 384-well plates, and 5,000–40,000 cells per well for 96-well plates.

Cell Seeding

The size of the experiment dictates the number of cells required. A typical CHO-based 100-plate suspension cell assay requires approximately $250\text{--}300 \times 10^6$ cells in the suspension cell flask. This can be accomplished by culturing cells in multiple layer flasks. Cells should be split 2–3 days prior to the assay and not be more than 80–90% confluent prior to antibiotic removal. All flasks should be incubated at 37°C (98.6°F) and 5% CO₂. One day prior to testing, spent media from each flask should be replaced with growth media that does not contain selection antibiotics.

Starting the System

To ensure proper system function, power the instrument well before you begin an experiment. The camera requires about five minutes to reach operating temperature, and if you are heating the plate stage, **Instrument > Manual Operation > Temperature Control**, allow 15 minutes for complete temperature equilibration. View the **Instrument Status** panel for the stage temperature.

To start the system:

1. Verify that the appropriate pipettor head, wash reservoir, LED banks and filters are in place before powering on the instrument.
2. Follow the procedure in [Powering On the System on page 18](#).
3. If incorrect components are installed, see the appropriate section in [Exchanging Hardware Components on page 186](#) for installation instructions.



CAUTION! Do not attempt to change pipettor heads, wash reservoirs, LED banks and filters while the instrument is running. All features must be exchanged in the appropriate instrument mode. When the instrument is started for the first time after a re-configuration of LEDs or emission filters, optics calibration is required.

Checking the System

Once a day, before running your first assay plate, do the following system checks:

1. Make sure that the system is turned on and the camera is cooled to proper operating temperature.
 - EMCCD camera: -70°C (-94°F) $\pm 2^{\circ}\text{C}$
 - HS EMCCD camera: -70°C (-94°F) $\pm 2^{\circ}\text{C}$
 - ICCD camera: -20°C (-4°F) $\pm 5^{\circ}\text{C}$
2. Run a Yellow Plate Signal Test to ensure the system is operating according to specifications. See [Running the Yellow Plate Signal Test on page 136](#).

A yellow signal test plate is provided with each corresponding pipettor head supplied with the system.

A signal test plate only needs to be run once a day unless system components such as pipettor heads or LEDs are changed.



CAUTION! To avoid damaging the yellow signal test plates:

Avoid scratching the plate bottom because scratches can affect the standard deviation.

Store the plates in a safe place away from bright light on an even surface.

Dye-Loading the Cells for Fluorescence Assays

Many cell-based responses require fluorescent dye loaded in cells to bind or change conformation in the presence of a ligand. These changes are monitored through a shift in emission wavelengths and captured by a CCD camera. Calcium sensitive dyes, such as the ones used in Molecular Devices FLIPR Calcium Assay kits, monitor intracellular calcium flux assays, or environmentally sensitive for membrane potential assays. Depending on the application evaluated, dye-loading typically lasts from 30 minutes to 2 hours, at either room temperature or 37°C (98.6°F).

Dye-Loading Duration and Temperature

Optimal dye-loading time depends on the cell type and presence of an anion exchange inhibitor. For most cell lines, especially for calcium mobilization assay development, start with a 60 minute to 120 minute loading time at 37°C (98.6°F).



CAUTION! Avoid exceeding the optimal loading time. If anion exchange inhibition does not enhance dye-loading conditions, it should not be used.



Tip: If loading for 30 minutes yields an acceptable fluorescence signal, as has been observed for some cell lines, use a shorter loading time.

In some cases, incubation at room temperature can work as well or better than 37°C (98.6°F).

Preparing a Source and Compound Plate

Preparation Time for the Source Plate

Depending on assay complexity, source plate preparation time may vary. To avoid conflicts, plan your experiment carefully to ensure source plates are ready before dye loading is complete.

Recommended Source Plates

Compound conservation can be achieved by reducing dead volumes and minimizing source plate adherence. Reducing source plate adherence help to ensure proper compound concentrations are delivered to a read plate. Polypropylene plates are most commonly used for this purpose because they are solvent-resistant, can withstand repeated freeze-thaws, and have a low retention. In addition, proteins are less likely to adhere to polypropylene plates rather than polystyrene surface.

Manufacturers offer a wide selection of plate bottom shapes (U-, V- or flat bottom) to lessen dead volumes. Dead volumes decrease as you move from a flat to U- to V-bottom plates. Contact the plate manufacturer for the dead volume in the plates you are using.

Concentration of Compounds in the Source Plate

Compound concentrations are prepared based on ratio of addition to initial well volume of the read plate. Common concentrations are 3X, 4X or 5X because ratio of addition to read plate volumes are 1:3, 1:4 or 1:5, respectively. Volume of addition is dependent on compound mixing efficiency, cell adherence, and kinetics of cellular response. Assay development should always be performed to determine optimal addition volume and concentrations prior to screening.

Addition and Mixing of Compounds to the Cell Plate

Prompt mixing is required to initiate robust cellular kinetics associated with signal transduction assays. Proper addition parameters (for example, dispense speed, height, and volume) initiate a rapid response upon addition. Slow signal increases or variation in signal between wells may be signs that mixing is too slow or not uniform. Optimal addition parameters typically consist of a fast dispense speed, low pipettor height, and large addition volume. Assay optimization is always recommended to determine the proper assay conditions prior to running a screen.

If addition parameters are not optimized, artifacts occur affecting data quality. One common problem is adding compounds too rapidly. Weakly adherent cells may dislodge from the monolayer, causing a decrease in fluorescence upon addition. Cell displacement can be reduced by slowing dispense speed while decreasing addition volume and increasing height. Larger dispense volumes mix rapidly because they encounter less surface tension as compared to smaller volumes. However, small sample volumes are less disruptive to a cell monolayer and reduce compound consumption. To avoid dislodging weakly adherent cells, compounds should be added in small volumes and at relatively low dispense speeds.

*** Tip:** Mixing is not typically used during addition in calcium or membrane potential assays. Signal artifacts, such as additional noise in the fluorescent trace, can be attributed to cell disruption induced by rapid pipettor movement. However, it is not uncommon to use mixing before aspirating, to resuspend reagents that have fallen out of solution in the source plate.

For temperature-controlled assays, cell and source plates should be brought to temperature outside of the instrument. It takes approximately 45 to 60 minutes in a 95% humid, CO₂ incubator (15 minutes in a heating block) to warm plates to equilibrium at 37°C (98.6°F).

Compound Plates for Suspension Assays

Unlike adherent assays, suspension cell assays require cell addition directly to the compound in the read plates. Compounds are suggested to be prepared at 2X concentration in the same diluent as used for cells, since equal volumes of cells are added to the read plate containing the compound.

In 384-well assay 25 µL cells are added to 25 µL 2X compound; in a 1536-well assay 2 µL cells are added to 2 µL compound. Plates should be stored covered to prevent evaporation.

Setting Up an Assay Protocol

*** Tip:** Refer to the *FLIPR Penta High Throughput Cellular Screening System Protocol Guide* for specific assay protocols.

An effective way to reduce time and ensure consistent screening parameters is to pre-configure the protocol (*.fmp) files. You can define and save protocol files for each cell line, project, or user. Saved protocol files are opened from **File > Open > Protocol File**. In addition, the last 10 files (6 data and 4 protocol) opened in ScreenWorks Software are listed in the bottom portion of the **File** menu.

A protocol file defines the following experiment parameters:

- Instrument conditions, for example, read mode, plate positions, file names and file storage
- Assay steps, for example, fluid transfer and reads
- Analysis options, for example, groups, corrections and export parameters

To create a new protocol:

1. Open the **Process** tab panel. See [Process Tab on page 70](#).
2. In the **Process** tab panel, add the assays steps. Do one of the following:
 - To place the new process at the end of your developing protocol list, drag and drop the process anywhere in the Experiment window.
 - To place the new process somewhere specific in your developing protocol list, drag and drop the process where you want it. You are notified when the position is invalid.



3. When the configuration screen opens for the added process, adjust settings as needed. See [Understanding Processes on page 72](#).
4. Repeat steps 2 and 3 until you are finished building your protocol.
5. Define the **Analysis** parameters.

Click the **Analysis** process, and then click **Grouping**, **Correction**, and **Export** to specify the analysis parameters for the experiment data. See [Analysis Process on page 78](#).

- Click on any other process icon to open its configuration screen and make changes as needed.



Note: When using the HS EMCCD camera, depending on your **Protocol > Settings > Edit read mode > Camera Mode** setting, when using a fast Read time interval value, the plate matrix panel updating might be delayed until the end of the data acquisition rather than updating during the data acquisition.

Specifically, the following settings result in the plate matrix panel updating at the end of the data acquisition:

For **Camera Mode > HighSpeed**—a **Read time interval(s)** setting of 0.018 seconds or shorter

For **Camera Mode > Normal or Sensitivity**—a **Read time interval(s)** setting of 0.029 or shorter

- Select **File > Save**.
- In the **Save As** dialog, in the **File name** field, type a name for your protocol, and then click **Save**.



Tip: Protocol files are saved as an ***.fmp** file. Unless you change the **Save in** location for your protocols, the default save location is **C:\Documents\Molecular Devices\ScreenWorks\MyProtocols**.

- Run your experiment.

Refer to the *FLIPR Penta High-Throughput Cellular Screening System Protocol Guide* for specific protocol settings.

Optimizing the Optics and Fluid Dispensing

Optimizing the Hardware Settings

Prior to screening, the following hardware components should be optimized for signal output and detection:

- Excitation intensity
- Exposure time
- Number of reads
- Camera gain for EMCCD cameras
- Gate Open for ICCD camera
- Pipettor height
- Dispense speed
- Dispense volume
- Cell parameters for suspension assays


Optimizing Optics Hardware

Optimization is available to help amplify weak signals or reduce saturation. Modifying read settings might not alter the quality of your data because it only regulates the amount of light emitted or detected by the system. For instance, doubling the excitation intensity doubles the RLU's detected by the system, but your signal-to-noise ratio and Z-factor can remain the same.

To optimize the fluorescence signal:

1. Place the cell plates on the read positions.
2. With the **Camera Mode > Normal** setting, start with the excitation intensity, camera gain and exposure time settings shown below. These settings are the most frequently used to measure basal fluorescence signal.

| Camera Type | Calcium Assay | | | Membrane Potential Assay | | |
|----------------------------|---------------|----------|---------|--------------------------|----------|---------|
| | EMCCD | HS EMCCD | ICCD | EMCCD | HS EMCCD | ICCD |
| Excitation wavelength (nm) | 470–495 | 470–495 | 470–495 | 510–545 | 510–545 | 510–545 |
| Emission Wavelength (nm) | 515–575 | 515–575 | 515–575 | 565–625 | 565–625 | 565–625 |
| Excitation Intensity (%) | 50 | 40 | 50 | 50 | 50 | 50 |
| Exposure Time (s) | 0.4 | 0.05 | 0.533 | 0.4 | 0.005 | 0.533 |
| Camera Gain | 130 | 6.5 | 2000 | 50 | 1.5 | 2000 |
| Gate Open | — | — | 6% | — | — | 6% |

3. Run a  **Protocol Signal Test** to evaluate the basal fluorescence signal.

The fluorescence signal for intracellular calcium assays vary depending on the installed camera and settings. Look for the following fluorescence counts:

- EMCCD camera: 200–1,200
 - ICCD camera: 500–5,000
 - HS EMCCD camera: 4,000–6,0000
4. If the basal fluorescence signal is substantially out of these ranges, adjust the excitation intensity, exposure time, gain, or gate. Use the following suggestions to adjust the basal fluorescence:

Basal fluorescence too low:

- Increase the **Excitation Intensity**. During the assay, the **Excitation Intensity** ranges between 20% and 100%.
- Increase the **Exposure Time**. When you increase the **Exposure Time**, the read interval must be increased to a minimum of exposure time + 0.1 s. It takes 0.1 s to integrate the data.
- Increase the **Camera Gain**.
 - EMCCD camera: The gain varies between 1, extremely bright fluorescence, to 240, luminescence.
 - HS EMCCD camera: The gain varies between 1 and 128.
- Increase the **Gate Open**.
 - ICCD camera: The gate varies between 0.06%, extremely bright fluorescence, to 35%, very dim fluorescence.

Basal fluorescence too high:

- Decrease the **Excitation Intensity**.
- Decrease the **Exposure Time**.
- Decrease the **Camera Gain** (EMCCD and HS EMCCD camera).
- Decrease the **Gate Open** (ICCD camera).

Adjusting the Pipettor Height

To ensure prompt compound delivery, the pipettor height can be adjusted. It is based on pipettor mechanics and the physics of liquid. For Fluorescence an air gap, called the **Hold Volume**, is drawn into the tips after fluid aspiration from the source plate. This air gap ensures fluid doesn't leak out during pipettor movement. The hold volume is the first item to leave the pipette tip during dispense and enters the target well as a bubble. These bubbles can cause random light reflections and spurious signals. To avoid bubbles, dispense fluid with tips just above, but touching the initial well fluid level. This leaves the tips touching the meniscus after the addition has been completed to ensure complete sample dispensing.



Note: For Luminescence suspension assays, no Hold Volume is recommended along with low and fast addition.

If pipette tips are in the air at the end of fluid delivery, a drop can form on the tip end due to surface tension resulting in reproducibility problems. Therefore, the pipettor height should be set just above, but touching the starting fluid volume in the wells.

The fluid surface in a 384-well or 1536-well plate is a deeply curved meniscus. Set the pipettor height somewhere below the starting fluid volume for the 384-pipettor. There is no air gap used with the 1536-pipettor.

Adjusting the Fluid Dispensing Speed

Default dispense speed is 50 $\mu\text{L/s}$, 20 $\mu\text{L/s}$ and 2 $\mu\text{L/s}$ when dispensing to 96-well, 384-well, and 1536-well plates respectively. The table below provides recommended fluid dispense speeds for all three plate formats.

| Cell Conditions | 96-Well Plate ($\mu\text{L/s}$) | 384-Well Plate ($\mu\text{L/s}$) | 1536-Well Plate ($\mu\text{L/s}$) |
|--|-----------------------------------|------------------------------------|-------------------------------------|
| Slow dispensing speed for weakly adherent cells or non-adherent cells. | 10–50 | 5–20 | 1–6 |
| Fast dispensing speed for strongly adherent cells. | 100–200 | 25–50 | 5–10 |

These values must be experimentally determined for each cell type, but it is generally preferable to dispense as fast as possible to enhance mixing of the compounds in the wells. The trade-off is that the pipetting speed must not be so forceful as to dislodge cells from the well.

Optimizing Fluid Volume

The fluid volume parameters have the following range:

- 96-well plate = 5–200 μL
- 384-well plate = 1–25 μL
- 1536-well plate = 0.5–3 μL

Large volumes mix more rapidly into the wells than smaller ones. However, smaller sample volumes cause less disruption to the cell layer and allow the same source plate to be used for multiple cell plates. To avoid dislodging weakly adherent cells, smaller compound volumes should be added to the cell plate, and the pipettor should dispense the compounds at a relatively low speed.

Optimizing Pin Tool Delivery

To ensure proper transfer of compound using a pin tool to your target plate, evaluate the following factors during assay development. Controlling and standardizing the following factors for a given application helps you deliver reproducible volumes:

- Pin diameter
- Surface tension of the liquid being transferred
- Surface tension of the pin
- Speed of removal from source liquid
- Speed of pin striking recipient plate
- Depth to which the pin is submerged in the source plate
- Depth to which the pin is submerged in the target plate
- Volume of slot in pin
- Surface tensions of plate and dwell time



Note: The FLIPR Penta instrument and the FLIPR Tetra instrument use the same pin tools.

For more information regarding optimization of pin tools, refer to http://vp-scientific.com/molecular_devices.htm or contact:

V&P SCIENTIFIC, INC.

9823 Pacific Heights Boulevard, Suite T

San Diego, CA 92121

Toll Free: +1-800.455.0644

Phone: +1-858.455.0643

Fax: +1-858.455.0703

Email: sales@vp-scientific.com


Optimizing Cell Delivery

During suspension assays, cells are pumped from the external Cell Flask to the Cell Reservoir and pipetted to the read plate. There are several user adjustable settings for suspension cell optimization. These parameters need to be optimized for each cell line as each cell line is different. The settings include:

- Stir speed in the Cell Flask
- Fill pump speed
- Drain pump speed
- FLIPR Penta System aspirating/dispensing parameters

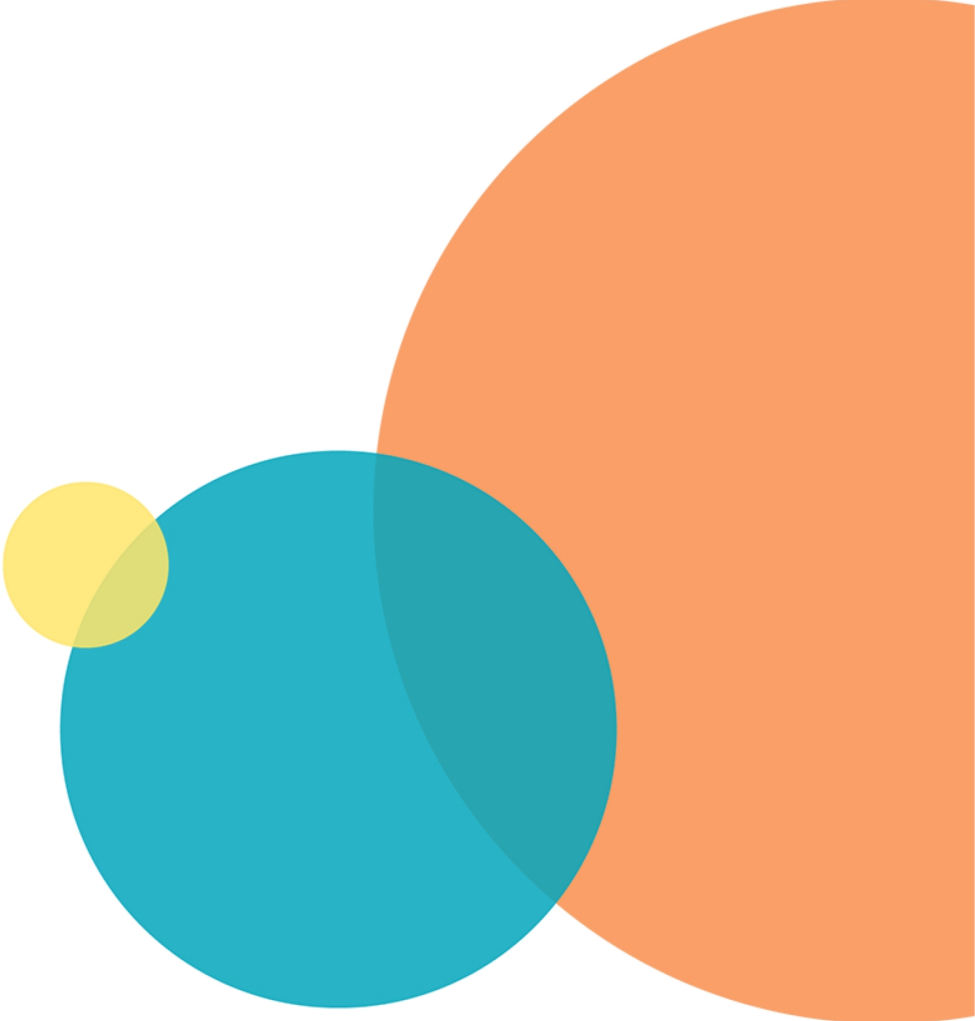
Starting the Assay Run

To start the assay:

1. Check the Experiment screen to verify that the **Experiment Summary** is accurate and that the appropriate Protocol *.fmp file is being used.
2. Verify the plates and tips are loaded in the instrument.
3. Save the protocol.
4. To start, click  **Run** on the toolbar.



CAUTION! The upper and lower doors automatically lock before any part in the FLIPR Penta System moves. These doors must remain closed until the end of the run, including through the duration of the pipettor tip wash.




Chapter 6: Exporting Data

You can export data the following two ways:

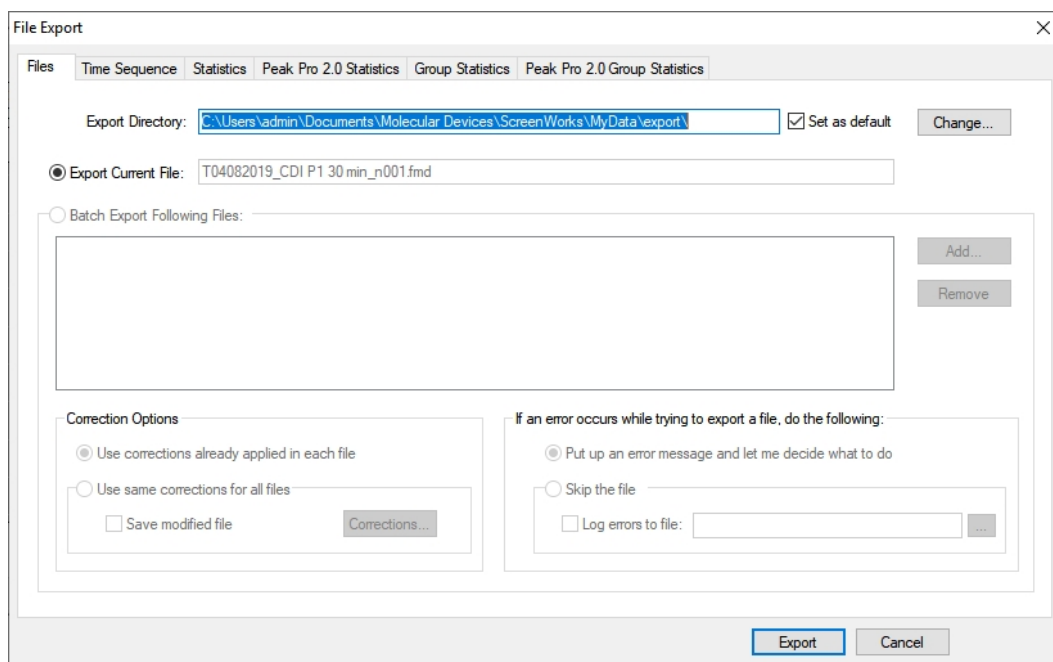
- Automatic export when a protocol is run.
- Export from already acquired data.

To configure the data export:

- In protocol files, in the **Analysis** screen, click  **Configure Auto Export**.
- For data files, use **File > Export, File** or **Batch Export**.



Note: **Auto-Export** and **Batch Export** do not support the ScreenWorks Peak Pro Software version 2.0.



Data is exported as ASCII text format files with a separate file exported for each measurement configuration specified. When you export data, you specify a folder in which to write the output files. The default export folder is:

C:\Documents and Settings\[your_user_name]\My Documents\Molecular Devices\ScreenWorks\MyData

The **File Export** dialog has the following tabs:

- **Files**—Export directory setting or batch files selections.
- **Time Sequence**—Exports time-point measurements for selected read modes. The measurement values that are exported have any corrections configured in the **Correction** dialog applied. If there are two read modes and **Ratiometric Options** is selected in the **Correction** dialog, you can also export the ratio for each time point. Files have a *.seqn extension, where n increments for each file generated in the export.
- **Statistics**—Exports averages, maximums, and other kinetic reduction values for selected numbers of reads for each well. Files have a *.statn extension, where n increments for each file generated in the export.
- **Peak Pro 2.0 Statistics**—You must run the Peak Pro 2.0 analysis before you can do the export. After running the analysis, this export works the same way as Statistics. See [Exporting Data With Peak Pro 2.0 Analysis on page 180](#).
- **Group Statistics**—Exports the group statistical values for selected numbers of reads for each group and are based on user-defined kinetic reduction settings. Files have a *.gstatn extension, where n increments for each file generated in the export.
- **Peak Pro 2.0 Group Statistics**—You must run the Peak Pro 2.0 analysis before you can do the export. After running the analysis, this export works the same way as Group Statistics. See [Exporting Data With Peak Pro 2.0 Analysis on page 180](#).



CAUTION! Data created in ScreenWorks Software version 2.0 cannot be viewed or exported by ScreenWorks Software version 5.0.

Exporting Time Sequence

Configure parameters for export of time-sequence data in the **Export Time Sequence** tab. The exported files contain a value for every read interval for each well, for the read mode or ratio selected. This contrasts with the Statistics file, which contains one kinetic reduction per well.

Table 6-1: Export Time Sequence Options

| Options | Description |
|----------------------------|---|
| Enable Time Sequence | Check this option to have a time sequence export file created. |
| File Name Options | Choose a name for the exported time sequence file. Use name of data file —Export files with the data or protocol file name, with *.seq <i>n</i> extension, where <i>n</i> is an integer ≥ 1 . Use user-defined name —Enter your own name for the export files (maximum of 25 characters). Files are given *.seq <i>n</i> extension, where <i>n</i> is an integer ≥ 1 . |
| Data Format | Format the output from this group of options: Show Labels —Check to include information about the processing options selected (for example, name of *.fmd file and well labels). Any corrections prior to export (e.g., negative control correction) are reported at the top of the exported ASCII file. Order of Well Data —Order the data by column (for example, A1, B1, C1) or by row (for example, A1, A2, A3). Set this option in accordance to the way your spreadsheet or database handles well data. Order Time Data —Select vertical or horizontal ordering of the data. |
| Misc. Parameters | Select experimental parameters and group statistics to export with the data. |
| Individual File Parameters | Use the bottom section of the dialog to choose the number and type of time sequence files to be created from the same data set. Number of Files —Click the Add Statistic button to add to the number of statistic files to be created from the same data set. Read Mode —Enter the read mode (or ratiometric data) from which you want to export data. From Quadrant —You can choose a specific quadrant from which to export, or export data from all quadrants. This enables you to export data from one plate into four different spreadsheet files, if desired. Note: This option is only available for data created in 384-well or 1536-well formats. |

Exporting Statistics

Use the **Export Statistics** tab to configure kinetic reductions, for example, averages and maximum and minimum values, for selected numbers of reads within each well. Multiple kinetic reductions can be configured, each resulting in another export file. Within each file only one value is reported per well.

Table 6-2: Export Statistics Options

| Options | Description |
|-------------------|---|
| Enable Statistics | Check this box to have a statistic file created. |
| File Name Options | <p>Choose a name for the exported statistic file.</p> <p>Use name of data file—Export files with the data or protocol file name, with *.stat<i>n</i> extension, where <i>n</i> is an integer ≥ 1.</p> <p>Use user-defined name—Enter your own name for the export files (maximum of 25 characters). Files are given *.stat<i>n</i> extension, where <i>n</i> is an integer ≥ 1.</p> |
| Data Format | <p>Format the output from this group of options:</p> <p>Show Labels—When checked, the output file contains information about the processing options selected (for example, name of *.fmd file and well labels). Any corrections done prior to exporting the data (e.g., negative control correction) are indicated in text format at the top of the exported ASCII file.</p> <p>Order of Well Data—Order the data by column (for example, A1, B1, C1) or by row (for example, A1, A2, A3). Set this option in accordance to the way your spreadsheet or database handles well data.</p> |
| Misc. Parameters | Select experiment parameters and group statistics to export with the data. |

Table 6-2: Export Statistics Options (continued)

| Options | Description |
|----------------------------|---|
| Individual File Parameters | <p>Use the bottom section of the dialog to designate the number and type of statistic files to be created from the same data set.</p> <p>Number of Files—Click the Add Statistic button to add to the number of statistic files to be created from the same data set.</p> <p>Compute—Select a kinetic reduction type. See Kinetic Reduction Types on page 81</p> <p>Start Read—Enter the first read number to be included in the kinetic reduction.</p> <p>End Read—Enter the last read number to be included in kinetic reduction. This can equal the Start Read, if you want to extract values from a single read.</p> <p>Read Mode—Enter the read mode (or ratiometric data) from which the data processes.</p> <p>From Quadrant—You can choose a specific quadrant from which to export, or export data from all quadrants. This enables you to export data from one plate into four different spreadsheet files, if desired.</p> <p>Note: This option is only available for data created in 384-well or 1536-well formats.</p> |

Exporting Group Statistics

Use the **Export Group Statistics** tab to configure a group statistics export report, for example, average, standard deviation and z-scores, based on the kinetic reduction types defined for a select number of reads as set up on the Analysis screen.



Note: If there are no groups defined on the Analysis screen, an empty report exports.

Table 6-3: Export Group Statistics Options

| Options | Description |
|-------------------------|---|
| Enable Group Statistics | Check this box to have a group statistic file created. |
| File Name Options | <p>Choose a name for the exported group statistic file.</p> <p>Use name of data file—Export files with the data or protocol file name, with *.gstatn extension, where <i>n</i> is an integer ≥ 1.</p> <p>Use user-defined name—Enter your own name for the export files (maximum of 25 characters). Files are given *.gstatn extension, where <i>n</i> is an integer ≥ 1.</p> |

Table 6-3: Export Group Statistics Options (continued)

| Options | Description |
|-------------------|--|
| Kinetic Reduction | <p>Configures the parameters used to define the kinetic reduction.</p> <p>Reduction Type—Defines the reduction to be applied to the kinetic data traces exported. See Kinetic Reduction Types on page 81 for details.</p> <p>Start Read—Define the first read to be used to determine the kinetic reduction.</p> <p>End Read—Define the last read used to determine the kinetic reduction.</p> <p>Read Mode—Select the read mode to apply the kinetic reduction.</p> |
| Misc. Parameters | Select experiment parameters to export with the data. |

Batch Exporting



Note: Batch Export does not support the separately licensed analysis module ScreenWorks Peak Pro Software version 2.0.

Manual export of files is accessed through **File > Export, File or Batch Export. Manual Export** allows the same files as defined above to be exported.

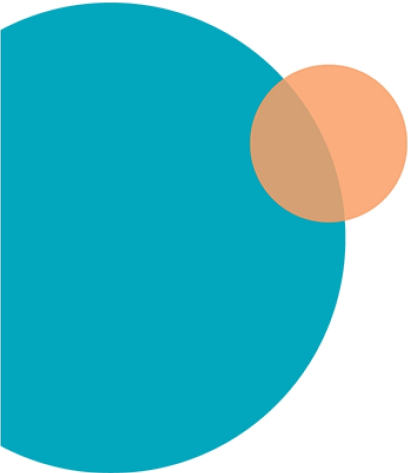
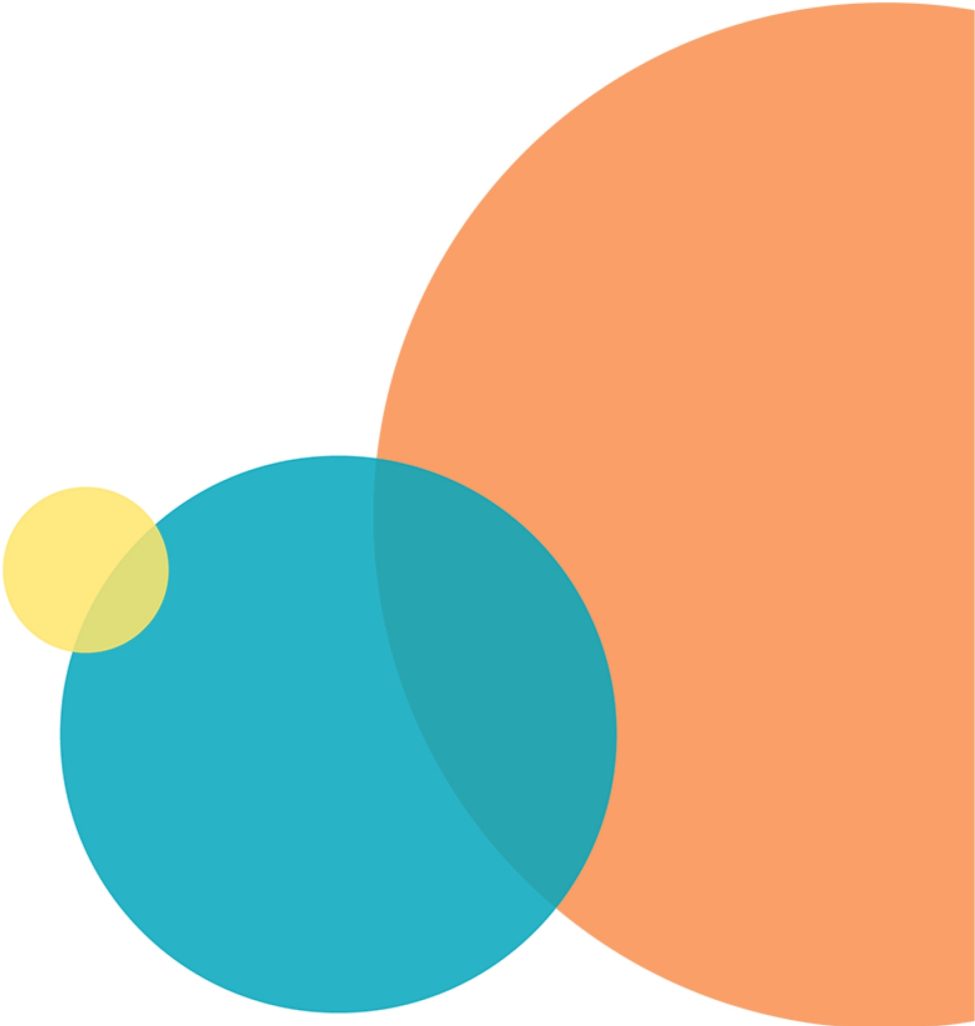
Table 6-4: Batch Export Options

| Options | Description |
|------------------------------|---|
| Export Directory | <p>User defines where export files are to be sent. The default export folder is:</p> <p>C:\Documents and Settings\[your_user_name]\My Documents\Molecular Devices\ScreenWorks\MyData</p> |
| Batch Export Following Files | <p>Selecting Add opens the Open File dialog to choose data files to be exported.</p> <p>Note: Hold down the SHIFT or CTRL key to select data files.</p> |
| Correction Options | <p>During export, either the existing or new corrections can be applied.</p> <p>Use corrections already applied in each data file—Applies the existing corrections saved with each data file during export.</p> <p>Use same corrections for all files—Applies and saves new corrections to all data files. Useful when exporting all data files with same parameters or saving the same correction to multiple data files. Details regarding available corrections can be found under Corrections on page 91.</p> |

Table 6-4: Batch Export Options (continued)

| Options | Description |
|--|---|
| If an error occurs while trying to export a file, do the following | <p>Put up an error message and let me decide what to do— Enables the user to decide how to proceed when an error is encountered.</p> <p>Skip the file— Does not include the data in the export, but continues to export data. The option to write the error to a log file is available.</p> |

When batch exporting data, the export files produced vary depending on the options you select in the individual export statistic, time sequence and group statistic sections. If a single export file is needed for each data file exported, select **Use name of data file** in the respective section you want to export. Export files created using this method are labeled with the name of the data file from which it was created. However, if a single export file is needed that contains information for multiple data files, select **Use user-defined name** in the respective section. In this instance, all exported information combines into one file labeled with the desired user-defined name. When the file is open, the individual data file names which the information was exported from are used as the header for each data set within the export file.



The ScreenWorks® Peak Pro™ Software Version 2.0 requires the purchase of a separate module license after a 14-day trial, which adds advanced peak detection and event characterization measurements and analysis functionality. This data analysis capability is especially useful with cardiomyocyte and neuronal protocols. Contact your Molecular Devices Sales Representative for purchasing. Detect and measure the activity of biological events that exhibit several components of interest, for example, multiple excitations, regular or irregular anomalous events such as early afterdepolarization-like events (EAD), irregular amplitudes or frequencies, and migrating threshold levels.



Note: For relatively fast events that have been acquired with long sampling intervals, use ScreenWorks Peak Pro Software version 1.0. See [Kinetic Reduction Types on page 81](#).

The Peak Pro 2.0 analysis algorithm reduces false detections without the need to condition the raw data. Using search vectors, you optimize the lengths of these vectors for specific data sets. They are generated with the midpoint of the vector at each point in the data set and are fitted by linear regression such that the slope at each point is known and the ups and downs of the adjacent noise events are dampened, which reduces the contribution of random noise to the overall movement of the vector.

For example, as shown in [Figure 7-1](#) and [Figure 7-2](#), adjusting the search vector length and specifying the minimum noise and amplitude values, very few false events appear.

In the following example, using a moderately noisy data set, using a search vector length setting of 3 results in a lot of false peaks:

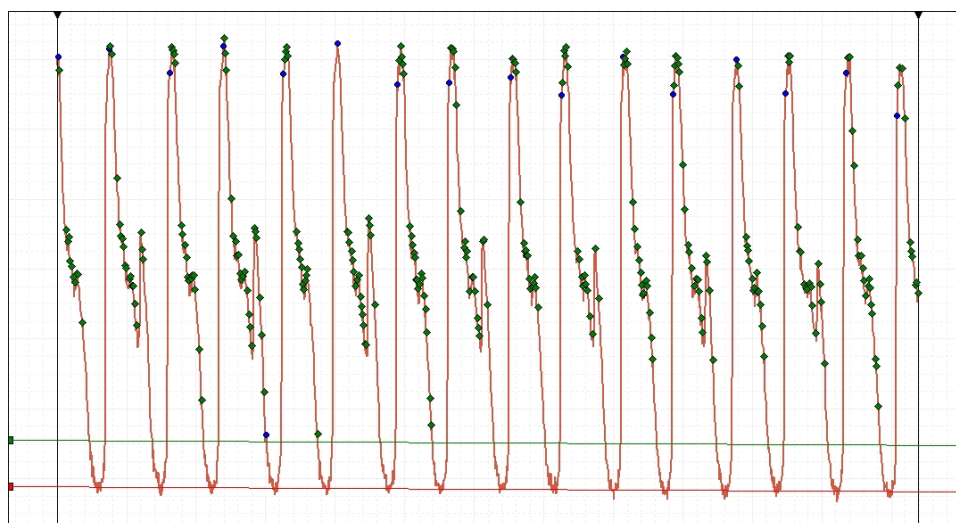


Figure 7-1: Search Vector Length of 3 is Noisy

Each marker represents a peak. The first peak of each event is a blue dot and the subsequent peaks are green diamonds.

Changing the search vector length setting to 11, and running the analysis again results in more precise data:

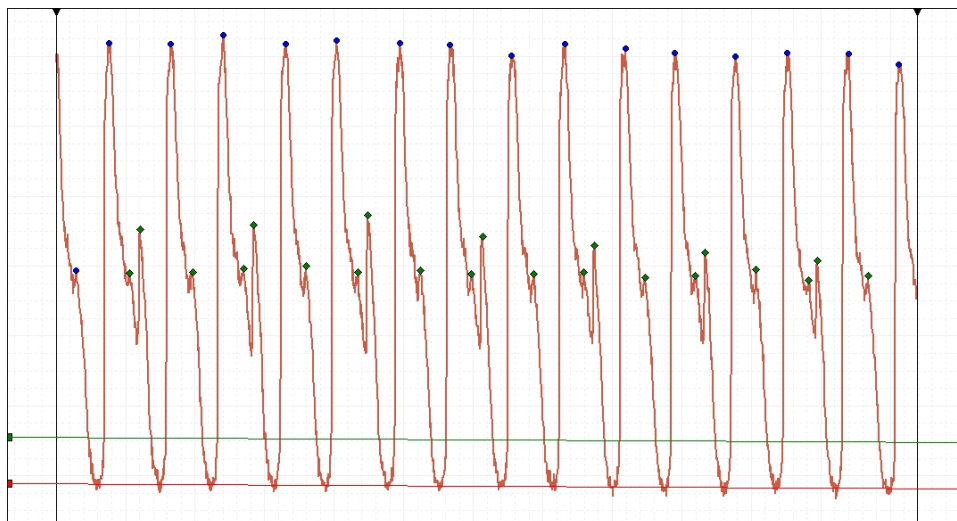


Figure 7-2: Search Vector Length of 11 Reduces Noise

The vector length did not need to be increased by much to see a substantial noise reduction improvement. The major peaks and minor peaks now appear to be real biological events.

The following sections explain how to use the Peak Pro 2.0 analysis software module in more detail. Topics include:

- [Interface Overview on page 167](#)
- [Running Peak Pro 2.0 Analysis on page 174](#)
- [Exporting Data With Peak Pro 2.0 Analysis on page 180](#)


Interface Overview

When you start the ScreenWorks Peak Pro Software version 2.0, a dialog appears with the open data file name in the title bar.

The Peak Pro 2.0 Analysis dialog has three tabs:

- [Options Tab Settings](#)
- [Measurements Tab Settings on page 171](#)
- [Configuration Tab Settings on page 174](#)

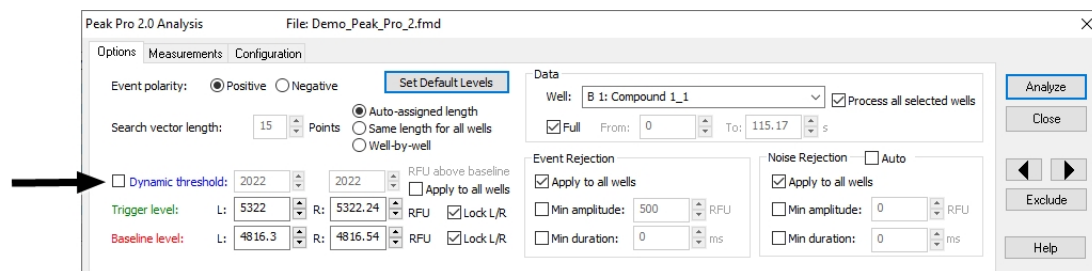
To open the Peak Pro 2.0 Analysis dialog:

1. Select **File > Open > Data File(s)**.
2. In the **Open** dialog, select your data file name, then click **Open**.
3. In the Multi-Well Graph, select individual wells, rows, or columns to analyze.
4. Click  to open the **Peak Pro 2.0 Analysis** dialog.

For more analysis procedures, see [Running Peak Pro 2.0 Analysis on page 174](#).

Options Tab Settings

The Options tab provides the peak detection search settings and data range settings.



The Options settings are organized into the following sections:

- [Data on page 169](#)
- [Event Rejection on page 170](#)
- [Noise Rejection on page 170](#)

- **Event polarity**—specifies the direction of the signal for detection, either Positive or Negative. Set by default for Positive.

*** Tip:** Negative-going polarity is rarely needed.

- **Set Default Levels**—resets the threshold, trigger, and baseline values to their default positions.
- **Search vector length**—determines the sensitivity of the peak detection. Longer vector lengths automatically filter out noise and short events, while shorter vector lengths detect lower amplitude transitions.

The search vector is generated from a linear regression across forward and backward points, centered on the mid-point of the vector, and computed for each data point. This vector slides along the data, checking the slope at each point, and determines peaks and valleys by noting the direction of the slope.

*** Tip:** This is the most important variable for good peak detection. Some experimentation with this setting might be necessary. You can specify a different vector length for each well.

- **Auto-assigned**—uses the sampling rate and an estimate of the noise amplitude to compute a vector length for each data point. By default, this check box is selected because it automatically determines the vector length suitable for most data sets.

*** Tip:** If needed for fine adjustment, deselect the Auto check box and manually enter Search vector length Points.

- **Same for all wells**—specifies one manually set vector length for the analysis of all the selected wells.
- **Well-by-well**—allows you to specify multiple manually set vector lengths for the analysis of wells. Each well can use a different vector length setting.
- **Dynamic threshold**—sets a lower limit for peak detection, rejecting all events below this threshold. Amplitude is relative to the baseline.
 - **Apply to all wells**—applies the specified threshold to all wells. When deselected, the value for an individual well is used.

- **Trigger level**—specifies when detection starts, with the first crossing of the level, and when the event ends, with the level re-crossing. By default, the level is positioned at 10% above the baseline, relative to the data range. Events that peak below the trigger level are ignored. Adjust as needed. This level can be set individually for each well. The signal values are relative fluorescence units (RFU).
 - **L**—specifies the left end of the trigger level line.
 - **R**—specifies the right end of the trigger level line.
 - **Lock L/R**—locks the left and right ends such that the level line moves as a unit when either the left or right values change.



Note: An event is the signal crossing the trigger level in both directions. The amplitude of the event, is measured relative to the baseline. The detection of the left and right antipeaks of the event is not limited by the trigger level and these measurements can extend as far as the baseline. The event duration is measured as the distance between the two antipeak positions.

- **Baseline level**—specifies the absolute end of the detection range. All amplitudes are measured relative to the baseline level and not the antipeak value. Adjust as needed. The signal values are relative fluorescence units (RFU). The algorithm automatically adapts to sloping baselines.
 - **L**—specifies the left end of the baseline level line.
 - **R**—specifies the right end of the baseline level line.
 - **Lock L/R**—locks the left and right ends such that the level line moves as a unit when either the left or right values change.

Data

- **Well**—individually lists all the selected wells from the plate matrix used to open the analysis dialog. This field also selects the graph that displays in the Detail Graph pane.
- **Process all selected wells**—includes in the process all of the selected wells in the Well list. When this check box is deselected, only the well visible in the Well field is processed.



Tip: You can select specific wells from the Well list and click Exclude for exceptions before analyzing all of the selected wells.

- **Full**—automatically selects the whole available data range.
- **From**—specifies the beginning point of your data range to be analyzed. Adjust as needed.
- **To**—specifies the end point of your data range to be analyzed. Adjust as needed.

Event Rejection

When you use Event Rejection, the amplitudes relative to the baseline that are lower than specified are rejected. Events with trough to trough times shorter than the specified duration are rejected.

- **Apply to all wells**—applies the specified amplitude and duration values to the data from all wells. When deselected, the value for each individual well is used.
- **Min amplitude**—specifies the lowest amplitude RFU value, measured from the baseline, for the acceptance of an event.
- **Min duration**—specifies the shortest duration ms value, duration, as measured between the left and right antipeak values, that is accepted.

Noise Rejection

When you use Noise Rejection, the amplitude is relative to the right trough on the rising phase and the left trough on the decay phase. Events with amplitudes lower than this are rejected. Events with right to left trough times that are less than the noise duration are rejected.



- **Auto** —Peak Pro 2.0 analysis measures the approximate noise level of the data averaging the difference between sequential amplitude values of all amplitudes within a distance of 10% of the baseline. Outliers are rejected using interquartile ranges. The noise estimate is then used to determine an approximate default noise level.
- **Apply to all wells**—applies the specified amplitude and duration values to the data from all wells. When deselected, the value for each individual well is used.
- **Min amplitude**—specifies the lowest amplitude RFU value before rejecting an event as noise. The measurement is taken between the lowest antipeak amplitude and the peak amplitude.
- **Min duration**—specifies the lowest duration time value before rejecting an event as noise. The measurement is taken between the antipeak time at the lowest antipeak and the time at the equivalent amplitude opposite the peak.

Measurements Tab Settings

The Measurements tab provides the properties of an event that can be selected for display in the Peak Pro 2.0 Statistics pane.

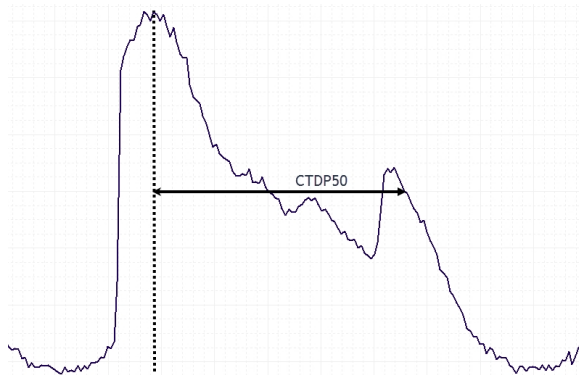
| Group Name | Well ID | Number of Peaks | Number of Peaks S.D. | Mean EAD-like peak Rate (PpM) | EAD-like peak Rate S.D. (PpM) | CTD @ 10% (s) | CTD @ 10% S.D. |
|-----------------------------|---------|-----------------|----------------------|-------------------------------|-------------------------------|---------------|----------------|
| Compound 1_1: Mean of wells | | 4.000 | 0.000 | | | 0.702 | 0.060 |
| B 1: Mean of events | B 1 | 4 | | | | 0.702 | 0.060 |
| Event 1 | B 1 | | | 170.795 | 118.585 | 0.660 | |
| Event 2 | B 1 | | | 301.752 | 330.202 | 0.744 | |
| Event 3 | B 1 | | | 244.533 | 249.135 | | |
| Event 4 | B 1 | | | 175.054 | 148.435 | | |
| Compound 1_8: Mean of wells | | 29.000 | 0.000 | | | 0.544 | 0.029 |
| O 2: Mean of events | O 2 | 29 | | | | 0.544 | 0.029 |
| Event 1 | O 2 | | | | | 0.576 | |
| Event 2 | O 2 | | | | | 0.563 | |
| Event 3 | O 2 | | | | | 0.537 | |
| Event 4 | O 2 | | | | | 0.584 | |
| Event 5 | O 2 | | | | | 0.579 | |
| Event 6 | O 2 | | | | | 0.563 | |
| Event 7 | O 2 | | | | | 0.568 | |

Figure 7-3: Peak Pro 2.0 Statistics Pane After Analysis With Expanded Group Data

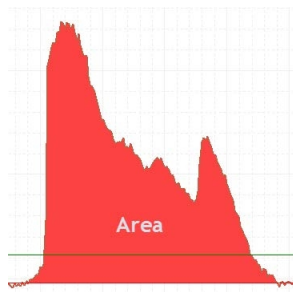
When you run an analysis, all of the following measurements are calculated:

- **Mean peak amplitude**—average amplitude of the main peaks of the events detected in the well data, measured from the peak to the baseline and expressed in relative fluorescence units (RFU).
- **Number of peaks**—total number of peaks.
- **Mean peak rate**—the numbers of peaks per minute (PpM). For cardiac data, this is equivalent to beats per minute (BPM).
- **EAD-like count**—reports the early afterdepolarization-like event (EAD-like) in an event. Includes Number of EAD-like events, Mean Number of EAD-like events, and Number of EAD-like events S.D..
- **Mean EAD-like rate**—average rate of early afterdepolarization-like event of the events detected in the well data, expressed in peaks per minute (PpM).
- **10-90% CTD**—Calcium Transient Duration. The width of a peak in seconds at various peak heights between 10% at the top and 90% at the bottom.

- **10-90% CTD**—Calcium Transient Duration from peak position measured as a straight line from the peak to a percentage of the distance from the peak.



- **Area**—below the event measured from the baseline to the start and end of the event as defined by the left and right antipeaks, expressed in relative fluorescence unit seconds.



- **Peak Spacing**—approximates the regularity of the peak spacing, expressed in seconds. The measurement reports either uniform spacing of peaks (OK), or irregular spacing (IRREG).
- **Rise slope**—rate of rise over the specified range. Expressed in relative fluorescence units per second.
- **Decay slope**—rate of decay over the specified range. Expressed in relative fluorescence units per second. Computed on if there is no early afterdepolarization-like event.
- **Rise time**—time to rising phase over the specified range.
- **Decay time**—time to decay over the specified range.
- **Linear run decay slope**—slope of a straight line from the peak to a specified percentage of the peak. Expressed in relative fluorescence units per second.
- **Linear run slope time**—from the peak. Expressed in seconds.

The following illustrates the measurement options and terminology:

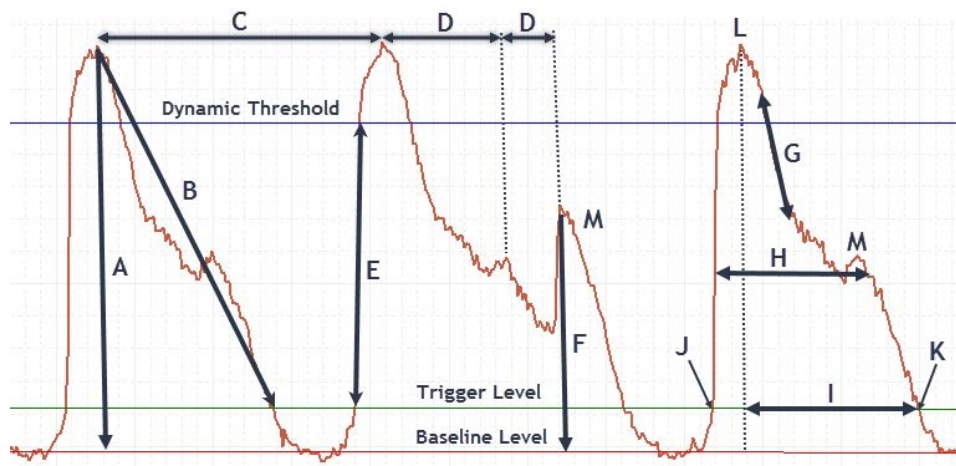
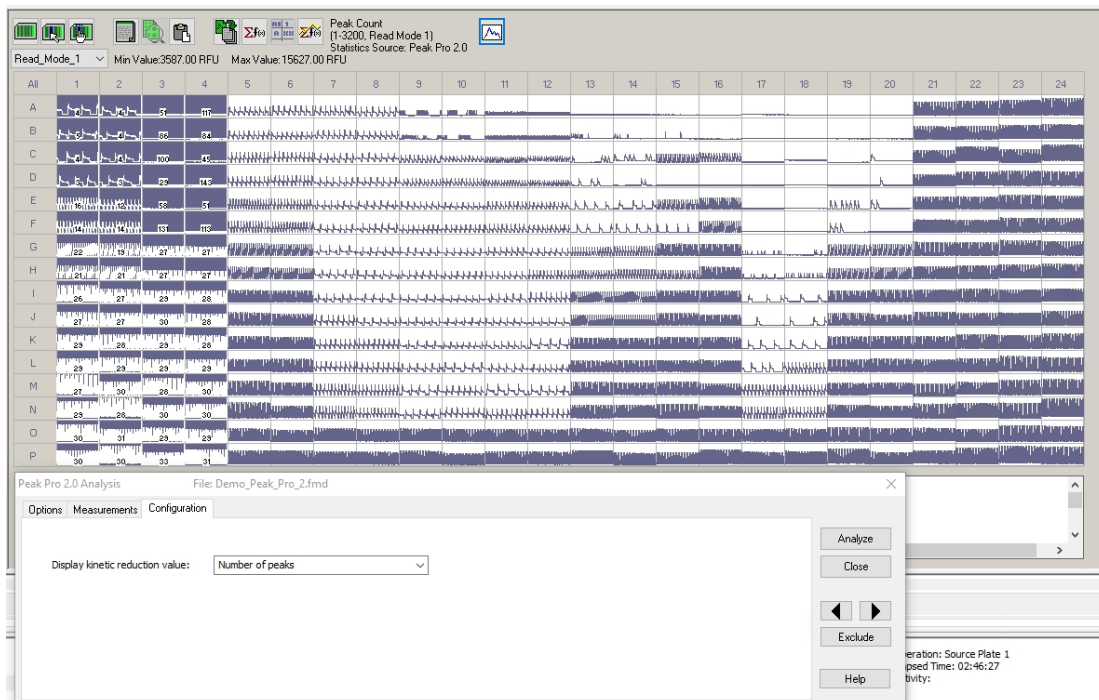


Table 7-1: Measurement Terminology Illustrated

| Item | Description |
|------|--|
| A | Main peak amplitude |
| B | Linear decay slope |
| C | Main peak interval used to calculate the peak rate, expressed in peaks per minute (PpM) |
| D | Early afterdepolarization-like event (EAD-like) intervals used to calculate the EAD-like event rate, expressed in peaks per minute (PpM) |
| E | Rise slope |
| F | EAD-like event amplitude |
| G | Decay slope |
| H | Calcium transient duration (CTD) |
| I | Calcium transient duration from peak position (CTDP90) |
| J | Start of an event |
| K | End of an event |
| L | Main peak |
| M | Early afterdepolarization-like event (EAD-like) |

Configuration Tab Settings

The Configuration tab provides the ability to display one kinetic reduction value in the Multi-Well Graph for your selected wells.



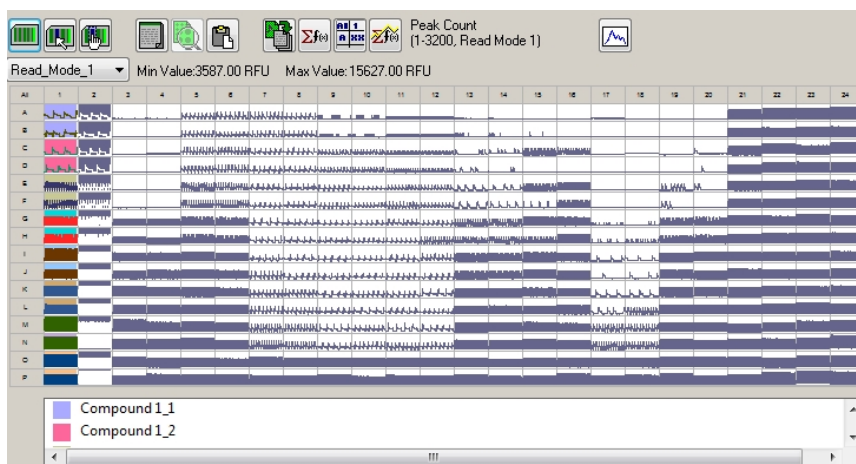
Running Peak Pro 2.0 Analysis


After you have saved experiment data files, you can run an analysis with the ScreenWorks Peak Pro Software Version 2.0 license. The Peak Pro 2.0 analysis functionality is best suited for cardiomyocyte, neuronal, and similarly fluxing assay data.


To run a Peak Pro 2.0 analysis:

1. Select **File > Open > Data File(s)**.
2. In the **Open** dialog, select your data file name, then click **Open**.


3. In the Multi-Well Graph, select the individual wells, rows, or columns to analyze.



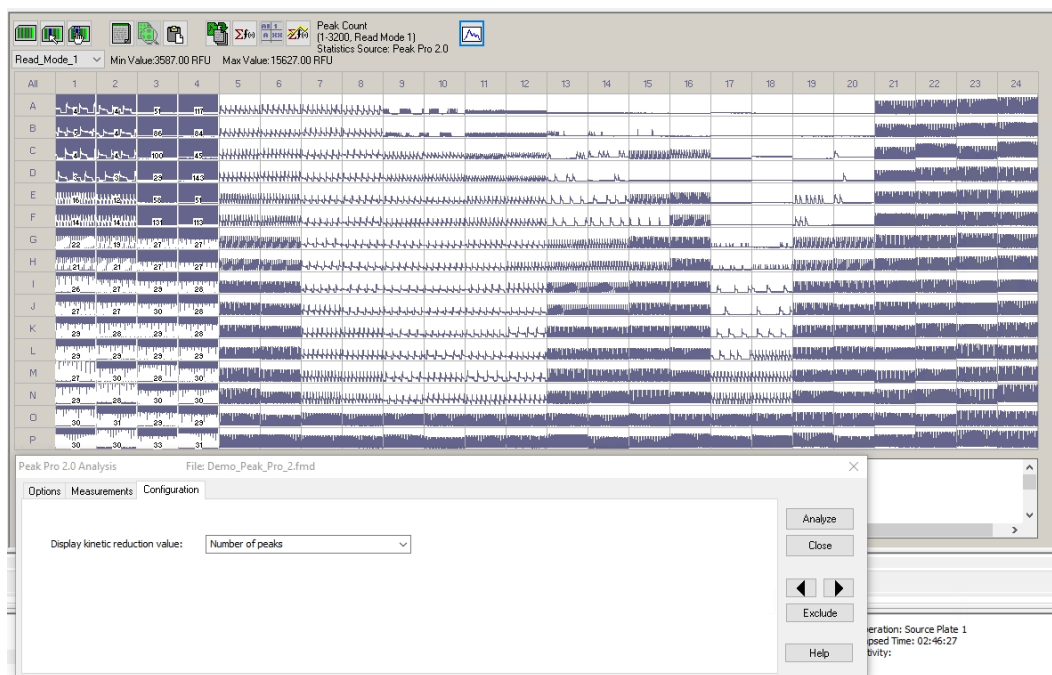
4. Click  to open the **Peak Pro 2.0 Analysis** dialog and collate the selected data, compute automatic baselines, thresholds, optimal search vector length, and format the results for subsequent analysis, sorting and recording.

 **Tip:** If needed, you can change your well selections.

5. In the **Options** tab, either run the default settings or manually adjust. See [Options Tab Settings on page 167](#).
6. In the **Measurements** tab, if needed, adjust ranges. See [Measurements Tab Settings on page 171](#).

 **Note:** All the listed measurements generate during the analysis.

- (Optional) In the **Configurations** tab, select one statistic type to **Display kinetic reduction value** in the Multi-Well Graph for your selected wells, for example, display **Number of peaks**.

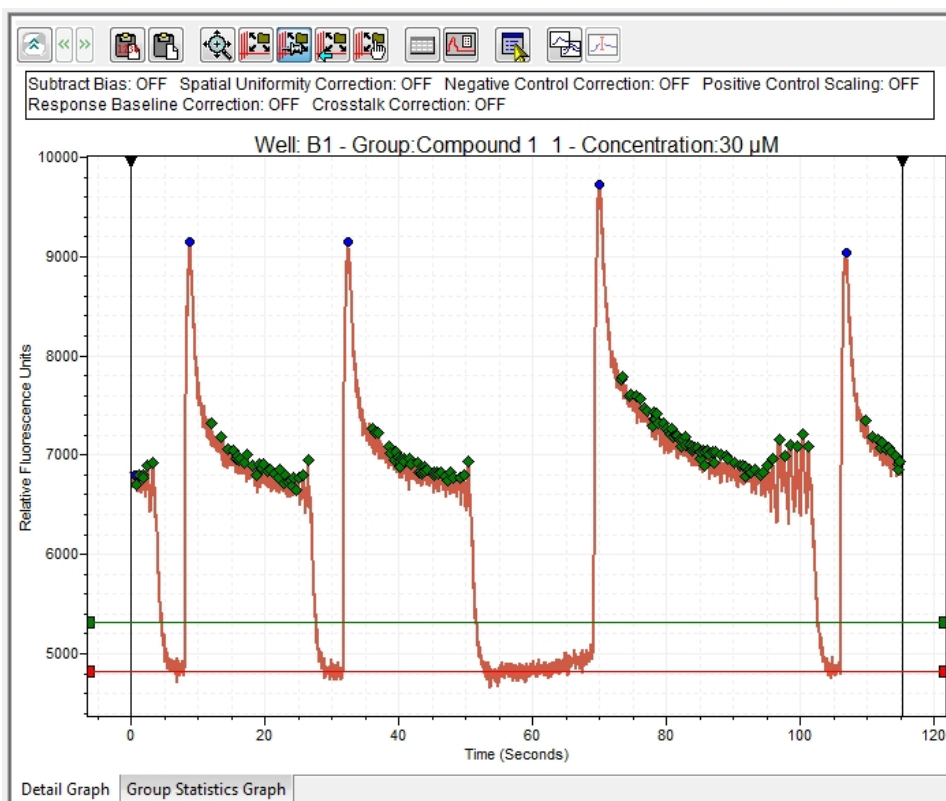


- To start the analysis, click **Analyze**.

9. Review the results in the **Detail Graph** pane.

Each major peak is marked with a blue dot. The green diamond markers are secondary peaks.

- * **Tip:** When there are a lot displayed, the secondary markers are questionably significant. Adjust the settings so that there is only one major peak and few secondary peaks per event.
- The major peak is the first peak in an event above the trigger level.
- An event is that portion of a signal that crosses the trigger level in both directions.

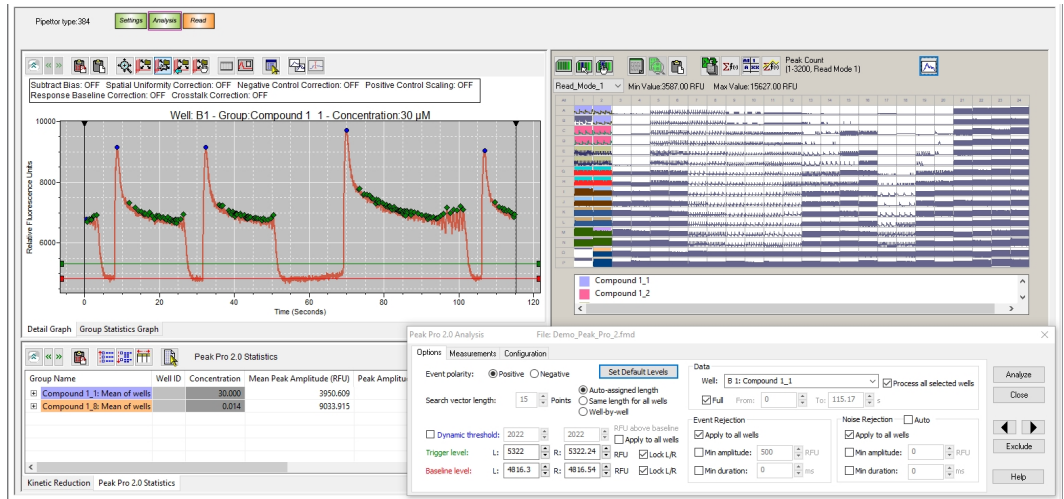


Using the **Peak Pro 2.0 Analysis** dialog arrow keys, you can scroll through each well graph to examine the result individually. Alternatively, you can select wells directly in the **Options** tab, in the **Data > Well** list.

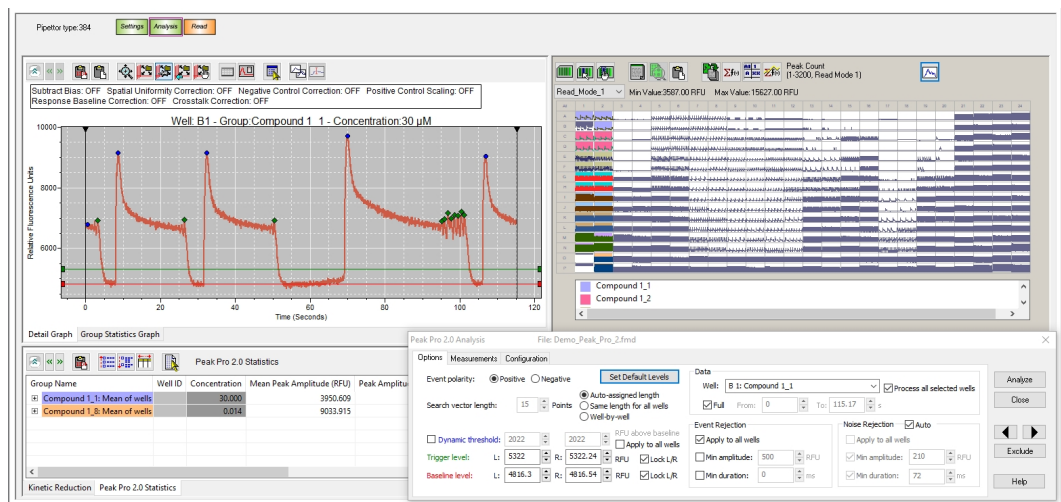
- If the first event marked is a partial signal that you want to skip in your analysis, adjust the analysis data time range.

To adjust the analysis data time range:

- Do one of the following:
 - In the graph, you can manually move the vertical start time range line.
 - In the **Options** tab, in the **Data** section, deselect **Full**, and type a **From** starting value.
- In the **Options** tab, click **Analyze**.




- To reduce the number of secondary peaks, which are likely noise and not biologically significant, adjust settings in the **Options** tab, click **Analyze**, and review the results again in the **Detail Graph** pane.

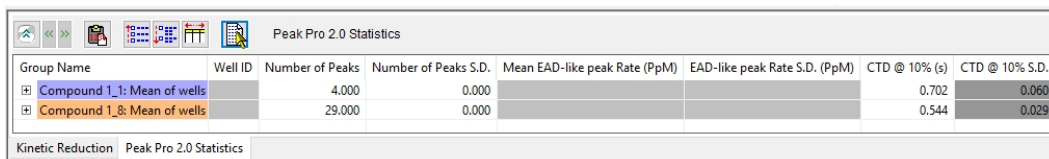


12. After you reduce all of the possible noise and have clearly identified the real biological events, you are finished with the analysis and can close the floating dialog.



Tip: To reopen the **Peak Pro 2.0 Analysis** dialog, click . Your last settings are retained in the dialog.

13. View the results in the **Peak Pro 2.0 Statistics** pane.



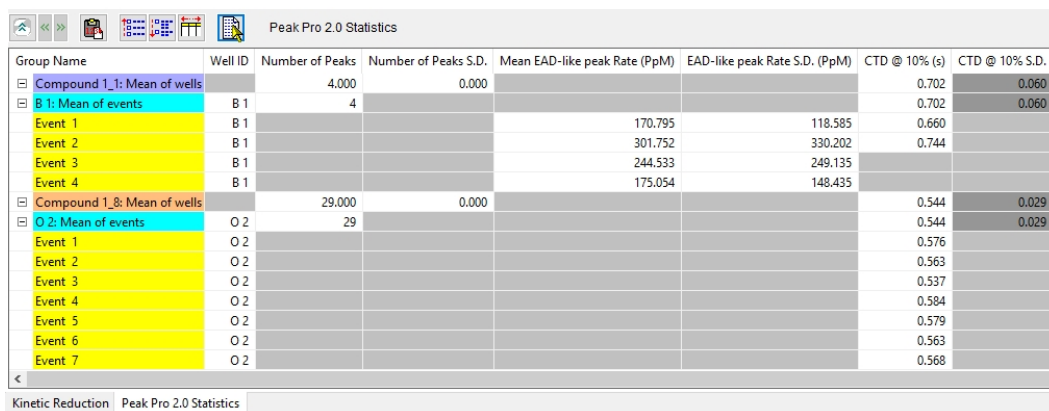
| Group Name | Well ID | Number of Peaks | Number of Peaks S.D. | Mean EAD-like peak Rate (PpM) | EAD-like peak Rate S.D. (PpM) | CTD @ 10% (s) | CTD @ 10% S.D. |
|-----------------------------|---------|-----------------|----------------------|-------------------------------|-------------------------------|---------------|----------------|
| Compound 1_1: Mean of wells | | 4.000 | 0.000 | | | 0.702 | 0.060 |
| Compound 1_8: Mean of wells | | 29.000 | 0.000 | | | 0.544 | 0.029 |

In the statistics pane, the results are sorted by **Group Name**. For help creating groups, see [Grouping on page 86](#).




Note: **Group Statistics Graph** functionality is not supported in ScreenWorks Software version 5.0.

14. Expand the **Group** rows and the **Mean** rows to see the **Event** rows of data.

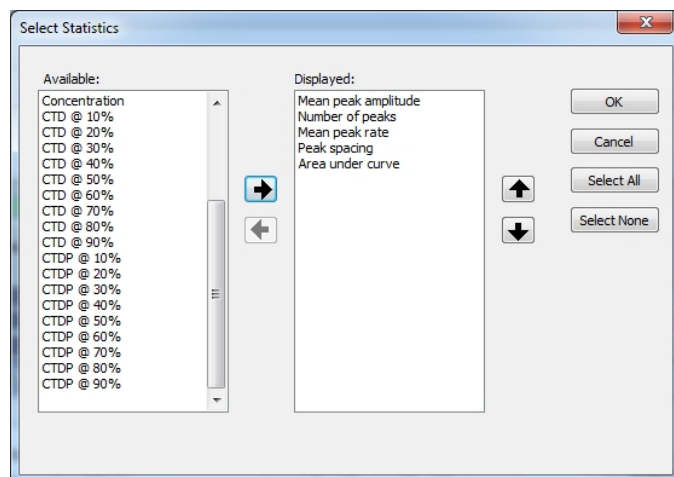


| Group Name | Well ID | Number of Peaks | Number of Peaks S.D. | Mean EAD-like peak Rate (PpM) | EAD-like peak Rate S.D. (PpM) | CTD @ 10% (s) | CTD @ 10% S.D. |
|-----------------------------|---------|-----------------|----------------------|-------------------------------|-------------------------------|---------------|----------------|
| Compound 1_1: Mean of wells | | 4.000 | 0.000 | | | 0.702 | 0.060 |
| B 1: Mean of events | B 1 | 4 | | | | 0.702 | 0.060 |
| Event 1 | B 1 | | | 170.795 | 118.585 | 0.660 | |
| Event 2 | B 1 | | | 301.752 | 330.202 | 0.744 | |
| Event 3 | B 1 | | | 244.533 | 249.135 | | |
| Event 4 | B 1 | | | 175.054 | 148.435 | | |
| Compound 1_8: Mean of wells | | 29.000 | 0.000 | | | 0.544 | 0.029 |
| O 2: Mean of events | O 2 | 29 | | | | 0.544 | 0.029 |
| Event 1 | O 2 | | | | | 0.576 | |
| Event 2 | O 2 | | | | | 0.563 | |
| Event 3 | O 2 | | | | | 0.537 | |
| Event 4 | O 2 | | | | | 0.584 | |
| Event 5 | O 2 | | | | | 0.579 | |
| Event 6 | O 2 | | | | | 0.563 | |
| Event 7 | O 2 | | | | | 0.568 | |

Each main row contains the average of measurements of each event as selected, and each event row contains the individual metrics of the individual peaks and related measurements.

15. To specify which of the data results display in the statistics pane, click  **Select Statistics**.

16. In the **Select Statistics** dialog, choose which statistics display and in which order.





17. When finished making changes, click **OK**.



Note: The **Peak Pro 2.0 Statistics** pane data never saves with the open data file used for analysis.

18. To save your analysis data, do either of the following:

- In the Multi-Well Graph pane, click  **Export**. See [Exporting Data With Peak Pro 2.0 Analysis](#).
- In the Statistics pane, click  **Copy Table Data to Clipboard** and then paste the delimited data into third-party software of your choice.

Externally, the saved data can be converted into graphs and reports.

Exporting Data With Peak Pro 2.0 Analysis

Exporting data for Peak Pro 2.0 is similar to the standard ScreenWorks Software exporting process, except that you must manually run a Peak Pro 2.0 analysis before you can export any Peak Pro 2.0 analysis data. For the standard exporting process, see [Exporting Data on page 157](#).




Note: An error message appears if you click **Export** before running your Peak Pro 2.0 analysis.



CAUTION! Because Peak Pro 2.0 analysis is a manually run process, **Auto-Export** and **Batch Export** functionality are not supported.

To export data with Peak Pro 2.0 analysis:

1. Run your Peak Pro 2.0 analysis. See [Running Peak Pro 2.0 Analysis on page 174](#).

2. Click  to open the **File Export** dialog.

- Click the tab for type of Peak Pro 2.0 analysis data export you want to do, **Peak Pro 2.0 Statistics** or **Peak Pro 2.0 Group Statistics**.
- Select **Enable**.



Note: After the 14-day trial expires, if you have not installed a ScreenWorks Peak Pro Software version 2.0 license, the **Peak Pro 2.0 Statistics** and **Peak Pro 2.0 Group Statistics** tab contents are disabled.

- Configure the data export in the selected tab. See [Exporting Peak Pro 2.0 Statistics on page 181](#) or [Exporting Peak Pro 2.0 Group Statistics on page 183](#).
- If you want to export data from the other tabs, repeat steps 3 through 5.
- Click **Export** to generate the data files from all of the enabled tabs.

Exporting Peak Pro 2.0 Statistics

After you run your Peak Pro 2.0 analysis, in the **File Export** dialog, use the **Peak Pro 2.0 Statistics** tab to configure kinetic reductions, for example, averages and maximum and minimum values, for selected numbers of reads within each well. Multiple kinetic reductions can be configured, each resulting in another export file, unless you select **Export all statistics to one file**. Within each file only one value is reported per well.

Table 7-2: Peak Pro 2.0 Statistics File Export Options

| Options | Description |
|-------------------|---|
| Enable Statistics | Select this box to have a statistic file created. |

Table 7-2: Peak Pro 2.0 Statistics File Export Options (continued)

| Options | Description |
|----------------------------|---|
| File Name Options | <p>Choose a name for the exported statistic file.</p> <p>Use name of data file—Export files with the data or protocol file name, with *.stat<i>n</i> extension, where <i>n</i> is an integer ≥ 1.</p> <p>Use user-defined name—Enter your own name for the export files (maximum of 25 characters). Files are given *.stat<i>n</i> extension, where <i>n</i> is an integer ≥ 1.</p> |
| Data Format | <p>Format the output from this group of options:</p> <p>Show Labels—When checked, the output file contains information about the processing options selected (for example, name of *.fmd file and well labels). Any corrections done prior to exporting the data (e.g., negative control correction) are indicated in text format at the top of the exported ASCII file.</p> <p>Order of Well Data—Order the data by column (e.g., A1, B1, C1) or by row (for example, A1, A2, A3) . Set this option in accordance to the way your spreadsheet or database handles well data.</p> |
| Misc. Parameters | <p>Select experiment parameters and group statistics to export with the data.</p> |
| Individual File Parameters | <p>Use the bottom section of the dialog to designate the number and type of statistic files to be created from the same data set.</p> <p>Number of Files—Click the Add Statistic button to add to the number of statistic files to be created from the same data set.</p> <p>Compute—Select a kinetic reduction type. See Measurements Tab Settings on page 171.</p> <p>Start Read—Enter the first read number to be included in the kinetic reduction.</p> <p>End Read—Enter the last read number to be included in kinetic reduction. This can equal the Start Read, if you want to extract values from a single read.</p> <p>Read Mode—Enter the read mode (or ratiometric data) from which the data processes.</p> <p>From Quadrant—You can choose a specific quadrant from which to export, or export data from all quadrants. This enables you to export data from one plate into four different spreadsheet files, if desired.</p> <p>Note: This option is only available for data created in 384-well or 1536-well formats.</p> |

Exporting Peak Pro 2.0 Group Statistics

After you run your Peak Pro 2.0 analysis, in the **File Export** dialog, use the **Peak Pro 2.0 Group Statistics** tab to configure group statistics, for example, average and standard deviation, based on the kinetic reduction defined for a select number of reads.

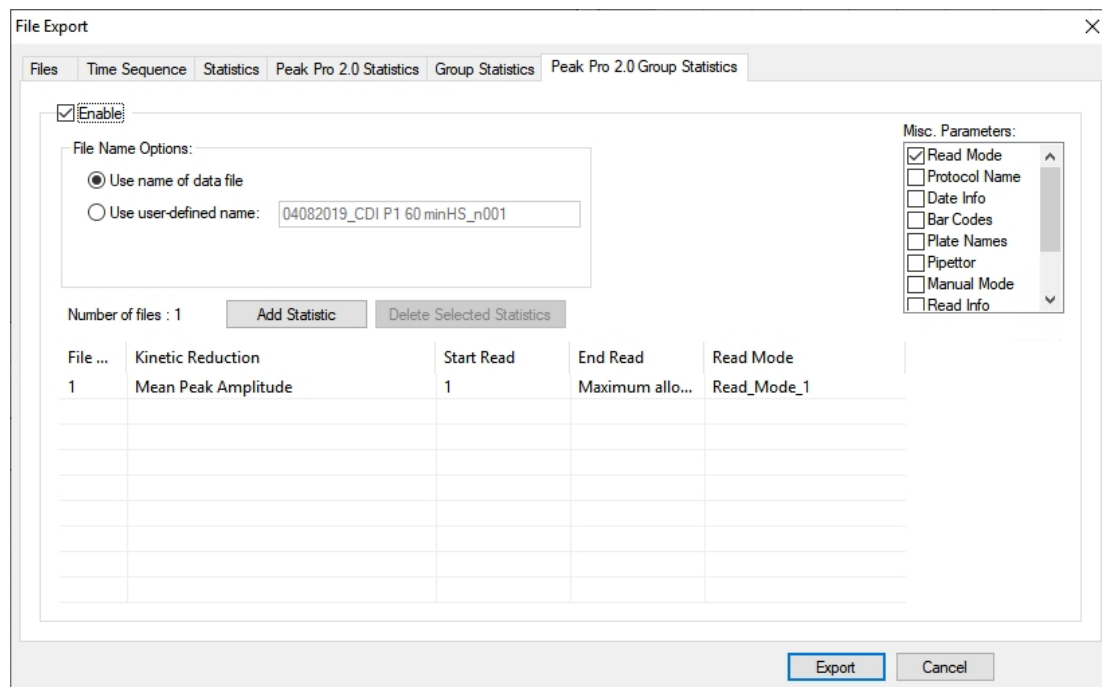


Table 7-3: Peak Pro 2.0 Group Statistics File Export Options

| Options | Description |
|-------------------------|--|
| Enable Group Statistics | Select this box to have a group statistic file created. |
| File Name Options | Choose a name for the exported group statistic file. Use name of data file —Export files with the data or protocol file name, with *.gstatn extension, where <i>n</i> is an integer ≥ 1 . Use user-defined name —Enter your own name for the export files (maximum of 25 characters). Files are given *.gstatn extension, where <i>n</i> is an integer ≥ 1 . |

Table 7-3: Peak Pro 2.0 Group Statistics File Export Options (continued)

| Options | Description |
|-------------------|--|
| Kinetic Reduction | <p>Configures the parameters used to define the kinetic reduction.</p> <p>Kinetic Reduction—Defines the reduction to be applied to the kinetic data traces exported. See Measurements Tab Settings on page 171 for details.</p> <p>Start Read—Define the first read to be used to determine the kinetic reduction.</p> <p>End Read—Define the last read used to determine the kinetic reduction.</p> <p>Read Mode—Select the read mode to apply the kinetic reduction.</p> |
| Misc. Parameters | Select experiment parameters to export with the data. |

Do only the maintenance tasks described in this guide. Maintenance procedures other than those specified in this guide must be done by Molecular Devices. See [Obtaining Support on page 226](#).

Before operating the instrument or following maintenance procedures, make sure you are familiar with the safety information in this guide. See [Safety Information on page 7](#).

The following are maintenance procedures that you can do as needed:

- [Washing Pipettor Tips and Pin Tools](#)
- [Exchanging Hardware Components on page 186](#)

Washing Pipettor Tips and Pin Tools

Pipettor tip and pin tool washing is controlled with the Wash Tips or Wash Pins process in ScreenWorks Software protocols and can be performed between fluid transfers within an experiment, or after the last fluid transfer, to prepare tips for the next experiment.

Disposable tips, as well as the 1536-tip head, and 384- and 1536-pin tools, can be washed.

The washer consists of a reservoir top of the selected pipettor format, mounted over a wash basin. Detailed instructions for exchanging the reservoir top are located in [Uninstalling Wash Reservoir Top on page 192](#). The wash basin is connected to two solvent-supply carboys and two waste carboys located on the floor beneath the computer monitor. A basin beneath the tip washer base drains to a waste carboy, to safely remove any solvent that overflows from the reservoir.

Wash solution fills the reservoir for a calibrated amount of time. Solvent is then drained from the reservoir after each wash cycle. Up to five wash cycles can be configured within a single wash process. For tips, a user-set volume of solvent is drawn up, optionally held for a time, and then expelled, up to 20 times. For pins, vertical motion of the tip block is used to agitate the wash solvent around the pins. The option is available to wash tips or pins in up to two solutions before reusing the tip washer. When additional wash solutions are required, tips or pins can be washed in a boat or reservoir, located in one of the source plate positions, using the **Mix Fluid** process.



WARNING! High volumes of volatile, flammable solvents in the reading chamber may cause explosive conditions. Use of 100% isopropanol, etc., in the tip washer is particularly discouraged without additional ventilation. Consult your facilities expert to determine the appropriate ventilation to avoid explosive conditions.

Pin tools are supplied with blotting stations that can be loaded into one of the plate positions. Blot pin steps can be configured in the protocol to remove fluid from the pins, for example, following pin washing.



Note: A waste bottle sensor override (P/N 0700-0827) is available in the FLIPR Penta System accessory kit to bypass the waste sensor and dispose waste in containers other than the dedicated waste carboys.

Exchanging Hardware Components

Use the following procedures to exchange hardware components on your FLIPR Penta System:

- Pipettor or Pin Tool Head—See [Exchanging Pipettor or Pin Tool Head](#)
- Pipettor Tips—See [Exchanging Pipettor Tips on page 194](#)
- Pin Tools—See [Exchanging Pin Tools on page 195](#)
- Alternating Position 1 Use—See [Alternating Position 1 Use on page 196](#)
- 1536 Microplate Tip Gasket—See [Exchanging the 1536 Microplate Tip Gasket on page 196](#)
- LED Modules—See [Exchanging LED Modules on page 198](#)
- Filters—See [Exchanging Filters on page 201](#)
- Plates or Tip Racks Hold-Down Devices—See [Exchanging the Plate Hold-Down Devices on page 205](#)
- Cell Reservoir—See [Exchanging the Cell Reservoir on page 204](#)

Exchanging Pipettor or Pin Tool Head

All pipettor and pin tool heads must be installed and calibrated on their intended instrument for the first time by a certified Molecular Devices Field Service Engineer. When calibrated, each head can be easily exchanged but only on the instrument for which they were calibrated.



Note: If you have any questions or concerns regarding the procedures, contact Molecular Devices Technical Support www.moleculardevices.com/service-support.

Exchanging the head type involves the following procedures:

1. [Uninstalling a Pipettor or Pin Tool Head on page 186](#)
2. [Installing the Pipettor or Pin Tool Head on page 190](#)
3. [Exchanging Pipettor Tips and Pin Tool Wash Reservoir Top on page 191](#)
4. [Resetting the System After Exchanging a Pipettor or Pin Tool Head on page 193](#)

Uninstalling a Pipettor or Pin Tool Head

The following procedure uninstalls a 384 pipettor head as an example.

To uninstall a 384 pipettor head:

1. In the ScreenWorks Software, unload all tips on the pipettor head by selecting **Instrument > Manual Operation > Unload Tips**.

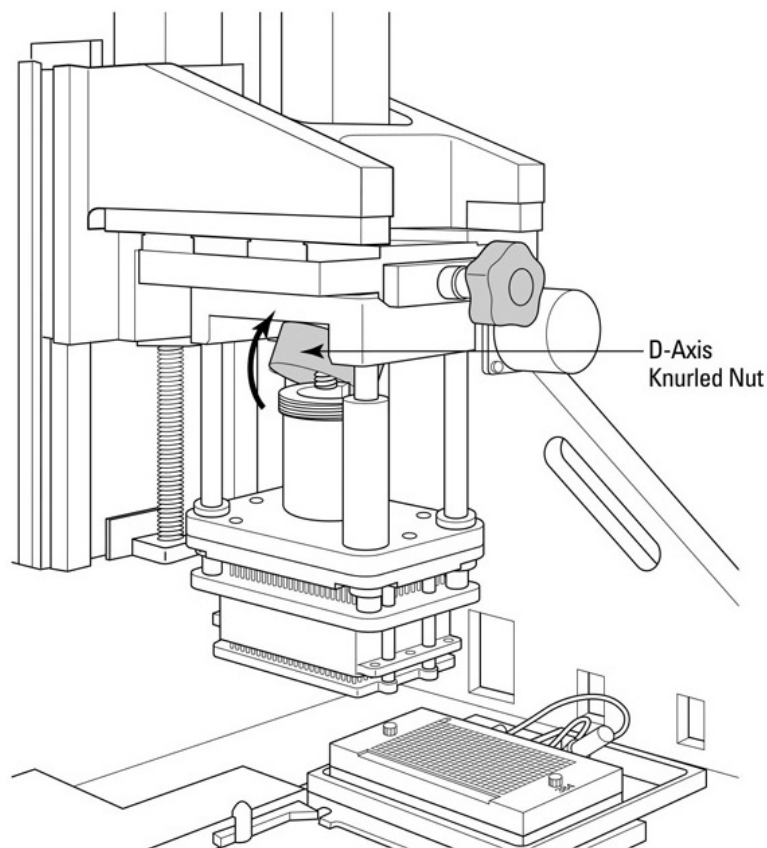
2. Select **Instrument > Manual Operation > Change Head**. This command instructs the pipettor head to move over the Read Plate position.



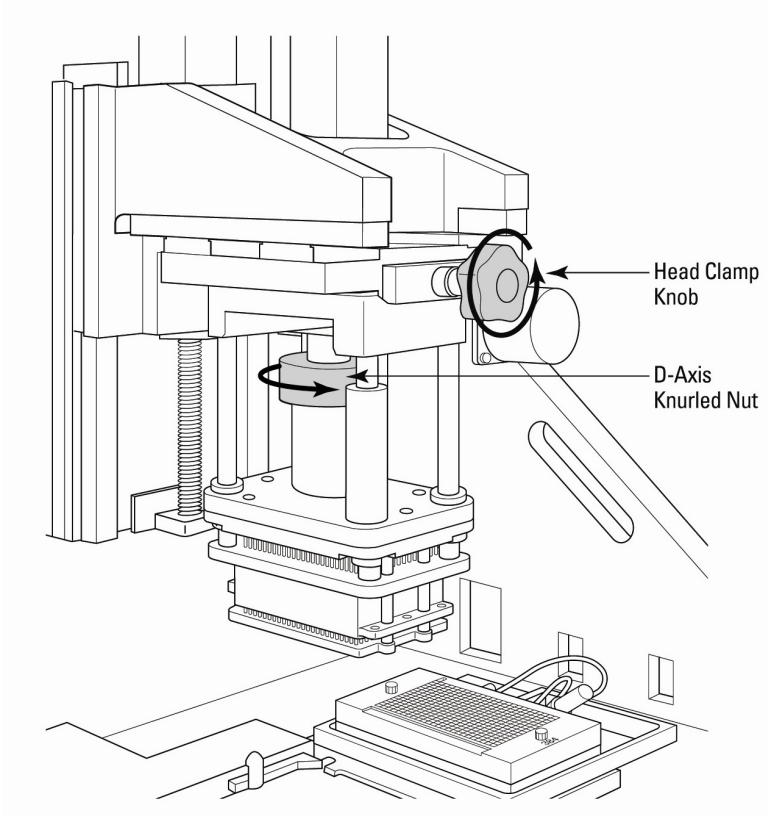
CAUTION! DO NOT select **DONE** in the dialog that opens until you have completed work on the pipettor.

3. Open the upper front door to access the pipettor head.
4. Turn the **D-Axis Knurled Nut** counter-clockwise until it loosens and can be lifted.

For 1536 pipettor heads only: After loosening the **D-Axis Knurled Nut**, install the **Pipettor Head Guard**, which seats over the 1536-pipettor head plungers, by tightening the two thumb screws. This guard protects the plungers from being damaged while the pipettor head is being handled outside the instrument.



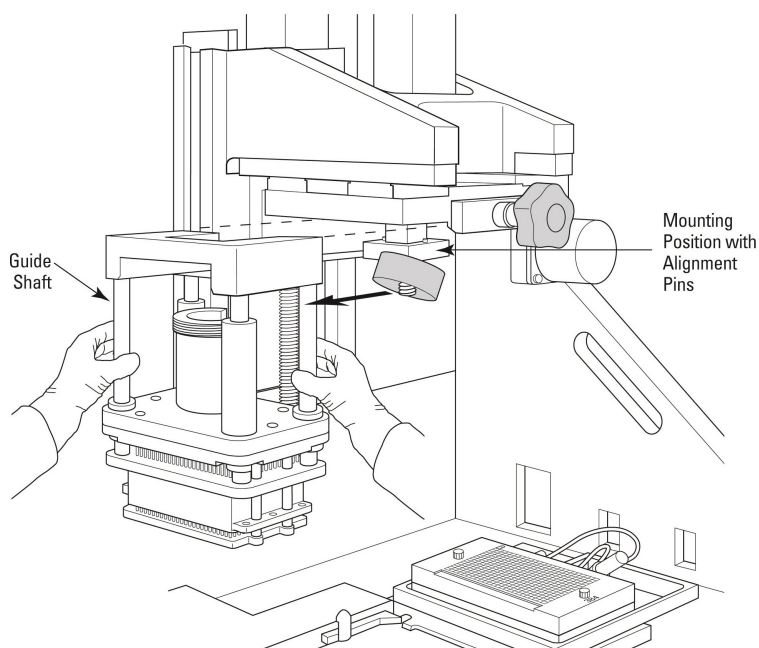
- 5. Turn the **Head Clamp Knob** counter-clockwise until the head is loose and can be lifted off its mounting position.



- To remove pipettor head, grasp the pipettor head silver **Guide Shaft**, lift upward (to unseat from **Alignment Pins** and to lower plunger pins allowing a lift gap) and then slide to the left off the pipettor mount.



Note: If the pipettor head does not easily slide off, try grabbing the D-axis knurled nut receiver and pressing it downward to give the pipettor head additional clearance from the D-axis knurled nut prior to trying again.



- Store the pipettor head in a safe location. A recommended safe location should include placing the pipettor head in its plastic storage bag to guard for particulates.



CAUTION! Do not place pipettor heads on nose cone surfaces because this can damage the nose cones. Store pipettor heads inverted, resting on the pipettor head top that seats on the pipettor mount position.

Installing the Pipettor or Pin Tool Head

The following procedure installs a 1536-pipettor head as an example.

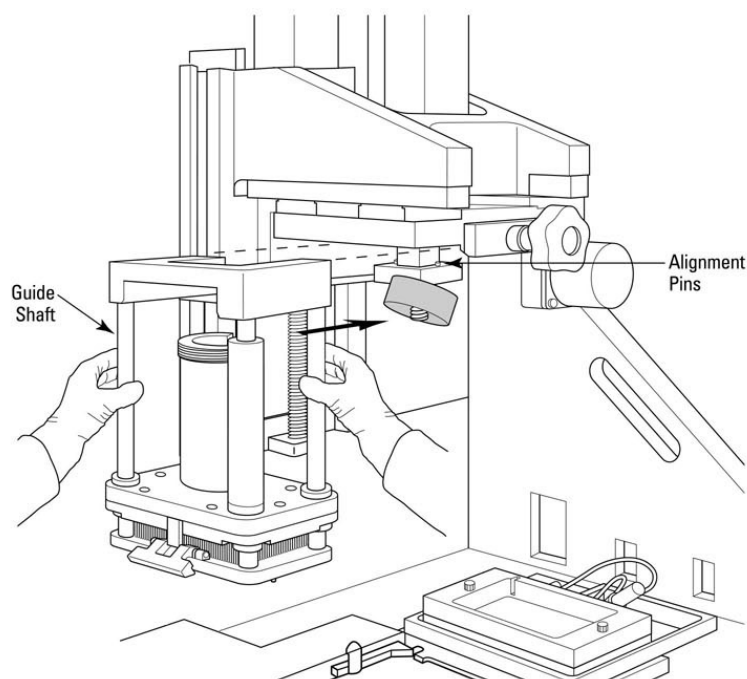


CAUTION! Only install a head that has been calibrated for your instrument by a Molecular Devices Field Service Engineer.

To install a 1536-pipettor head:

1. Hold the pipettor head by its silver **Guide Shafts** and slide it to the right over the **Pipettor Mount**.

When the pipettor head is in position, it should be seated on the **Alignment Pins** and should not slide off the pipettor mount without lifting upwards.

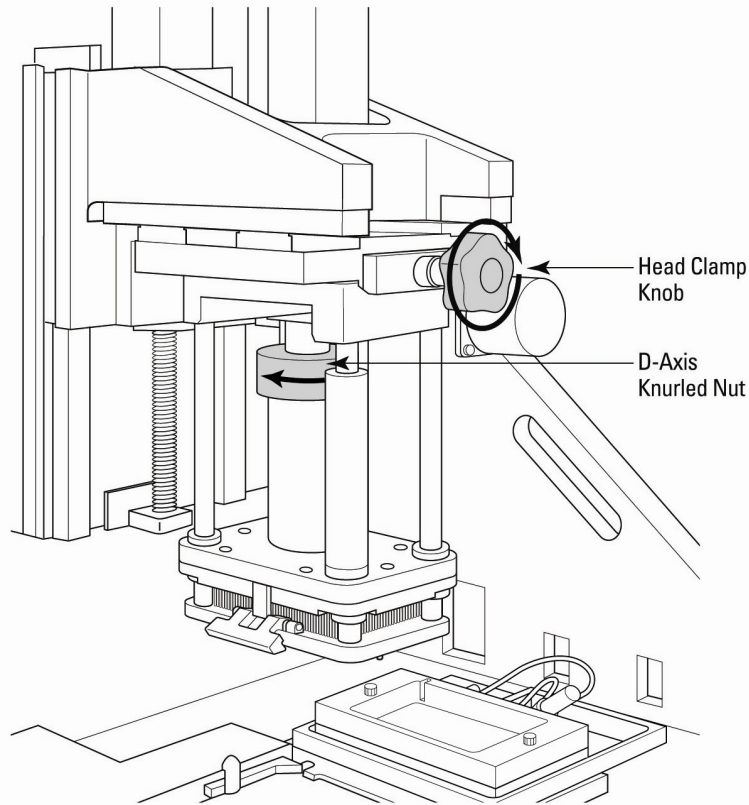


2. Tighten the **Head Clamp** knob by turning it clockwise.
3. **For 1536-pipettor heads only:** Remove the guard covering the 1536-pipettor head plungers by loosening the two thumb screws.
This allows the D-axis to move freely so it can be attached to the D-axis knurled nut in the next step.



CAUTION! If the guard is not removed before resetting the instrument, the 1536-pipettor head can be damaged.

4. Tighten the **D-Axis Knurled Nut** by turning it clockwise until tight.
If the nut does not reach its receiver threads, lift the D-axis receiver to engage the threads and tighten the nut.



5. After the head is fastened, exchange the tip wash reservoir top to match the head format. See [Exchanging Pipettor Tips and Pin Tool Wash Reservoir Top](#).

Exchanging Pipettor Tips and Pin Tool Wash Reservoir Top

After exchanging the pipettor tips or pin tool head, exchange the wash reservoir to match the new head format.



CAUTION! Failure to match head format and wash reservoir top causes damage during instrument operation.

Exchanging the wash reservoir top involves the following procedures:

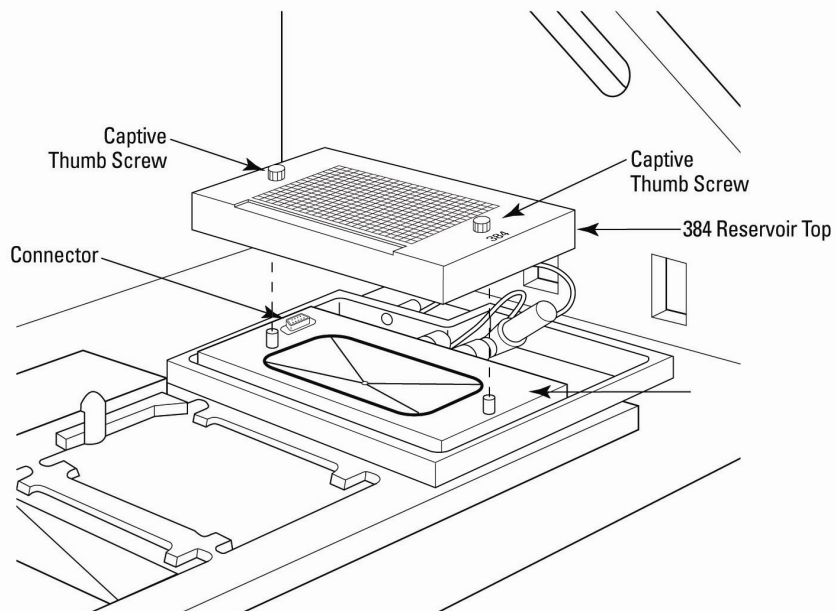
1. [Uninstalling Wash Reservoir Top on page 192](#)
2. [Installing Wash Reservoir Top on page 193](#)

Uninstalling Wash Reservoir Top

The following procedure uninstalls a 384 wash reservoir top as an example.

To uninstall a wash reservoir top:

1. Loosen both **Captive Thumb Screws** on reservoir top.



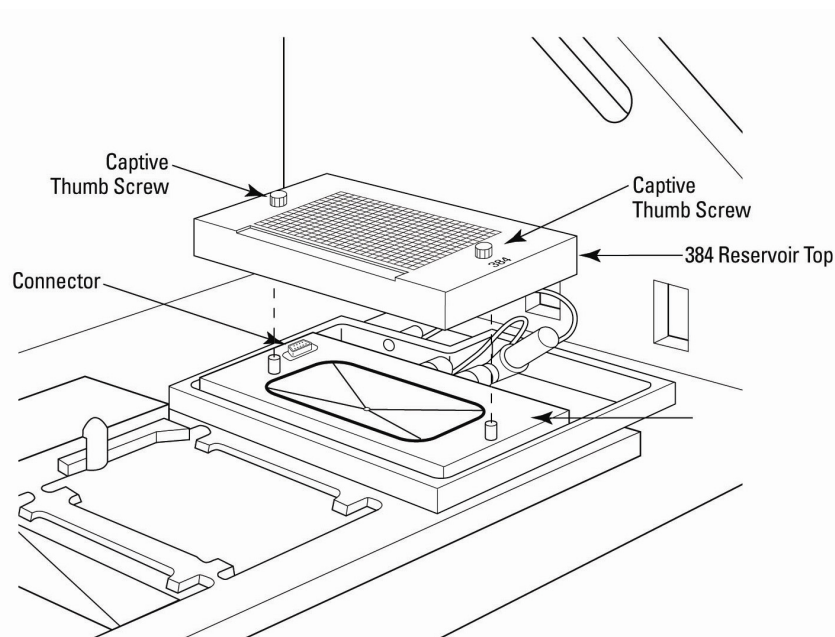
2. Lift reservoir top off of the washer base.
3. Ensure the **Blue O-Ring** on the washer base remains properly seated. If not, reinstall the o-ring into its o-ring groove.
4. Store the reservoir top in a safe location.

Installing Wash Reservoir Top

The following procedure installs a 384 wash reservoir top as an example.

To install a wash reservoir top:

1. Align the appropriate format reservoir top so the **Connector** matches the washer base connector. When aligned correctly, the reservoir top name (96, 384 or 1536) is located towards the front of the instrument.



2. When aligned, using the **Alignment Pins** to guide the placement of the top, press the top into position.
3. Tighten both **Captive Thumb Screws** to ensure the reservoir top is fastened to the base.
4. Go to [Resetting the System After Exchanging a Pipettor or Pin Tool Head](#) to reset the instrument.

Resetting the System After Exchanging a Pipettor or Pin Tool Head

To reset the FLIPR Penta System after exchanging a head and a wash reservoir top:

1. Make sure the head and tip wash reservoir top formats match before continuing.



CAUTION! If there is a head and wash reservoir top format mismatch, the instrument faults and prevents all experiment runs until corrected.

2. To finish the exchanges, click **DONE** in the software dialog that opens when you begin the head exchange. The instrument resets itself.
3. When the reset cycle is complete, the green **Assay Finished (Unlock)** light is the only indicator light illuminated on the front of the instrument.

Exchanging Pipettor Tips

The 96-pipettor head and the 384-pipettor head both use disposable plastic tips to transfer fluid.



Figure 8-1: 384-Well Tip Rack Example

The 1536-pipettor head uses a stainless steel tip block with a disposable 1536-tip gasket to transfer fluid. See [Exchanging the 1536 Microplate Tip Gasket on page 196](#).

Each pipettor head is compatible only with the tip racks of matching format. The pipettor head can only pick up a matching tip rack. For example, the 384-well pipettor head should not try to pick up pipettor tips from a 96-well tip rack.



CAUTION! Verify that there is a matching tip rack in the Tip Load position on the instrument deck before running your experiment. A mismatched pipettor head and tip rack causes a fatal error and can damage your equipment.



CAUTION! The FLIPR Penta System is only validated for use with the Molecular Devices tips. Instrument performance and troubleshooting is not guarantee with any other tips.

To automate pipettor tips loading:

- In your protocol configuration, select **Settings > Assign Plate to Position > Load Tips Position**. See [Settings Process on page 72](#) and [Creating New Protocols on page 124](#).

To load pipettor tips:

- In the ScreenWorks Software, select **Instrument > Manual Operation > Load Tip and Unload Tip**.

The pipettor tips attach to the pipettor head and the pipettor tips rack stays on the stage in Position 1.

To unload pipettor tips:

- In the ScreenWorks Software, select **Instrument > Manual Operation > Unload Tip**.

The pipettor tips are removed from the pipettor head in to the pipettor tip rack on the stage in Position 1.

For 1536-tip exchange procedures, see [Exchanging the 1536 Microplate Tip Gasket on page 196](#).

Exchanging Pin Tools

Pin tools are blocks of pins that are loaded onto a pin tool head to transfer fluid.

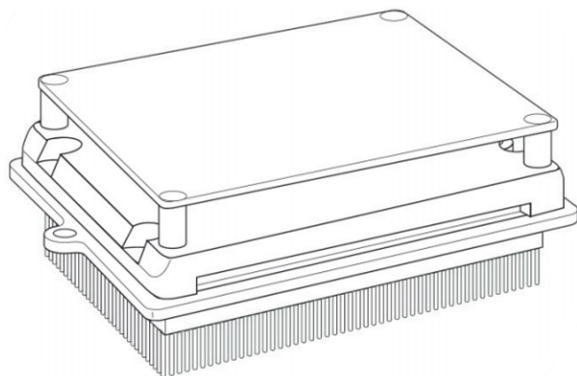


Figure 8-2: 1536-Well Pin Tool Block illustration

Two pin tool heads, 1536-well and 384-well formats, are available for the FLIPR Penta System. Each pin tool head is compatible only with the pin tool blocks of matching format. The pin tool head only picks up a matching pin tool block according to the attached magnetic ID. For example, the 384-well pin tool head does not pick up a 1536-well pin tool block. This is designed for the component safety.

To automate pipettor tips or a pin tool block loading:

- In your protocol configuration, select **Settings > Assign Plate to Position > Load Tips Position**. See [Settings Process on page 72](#) and [Creating New Protocols on page 124](#).

To load a pin tool block:

- In the ScreenWorks Software, select **Instrument > Manual Operation > Load Tip and Unload Tip**.

The pin tool block is picked up and clamped to the head using pin tool block grippers.

To unload a pin tool block:

- In the ScreenWorks Software, select **Instrument > Manual Operation > Unload Tip**.

The pin tool block is placed in Position 1 of the stage and unclamped from the pin tool head grippers.

Alternating Position 1 Use

Stage Position 1 is primarily for tip loading and unloading, but this position can also be used as a source plate position after tips are removed.



Note: Sensors indicate only the presence of an object in plate position and do not indicate whether it is a source plate or tip rack.

Use the following procedure in both manual and robotic modes.

To use a source well plate in the stage Position 1:

1. Load tips or pins manually.
2. Remove the tip rack before starting an experiment, and place a source plate in **Source 1**.

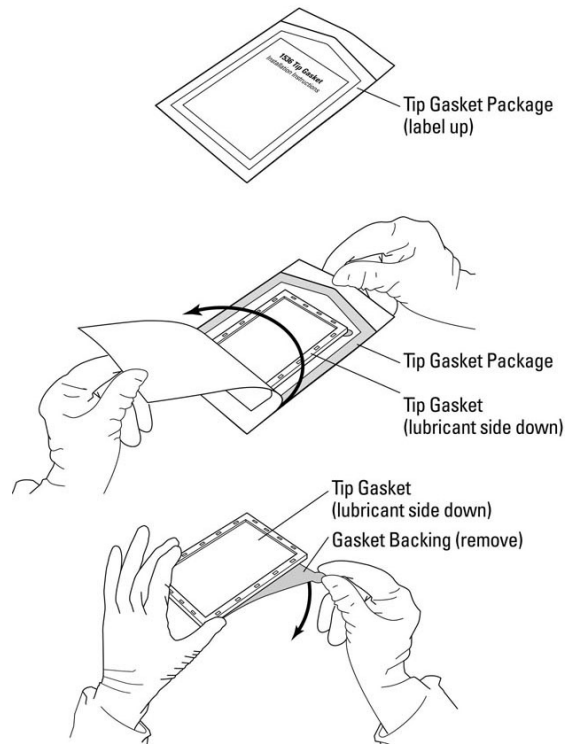
Exchanging the 1536 Microplate Tip Gasket

Gaskets should be replaced approximately every 200 cycles (1 cycle = 1 stroke down + 1 stroke up). Gasket lifetime should be monitored. It varies on the number of aspirate and dispense steps you run in your experiment.

Installing the Gasket

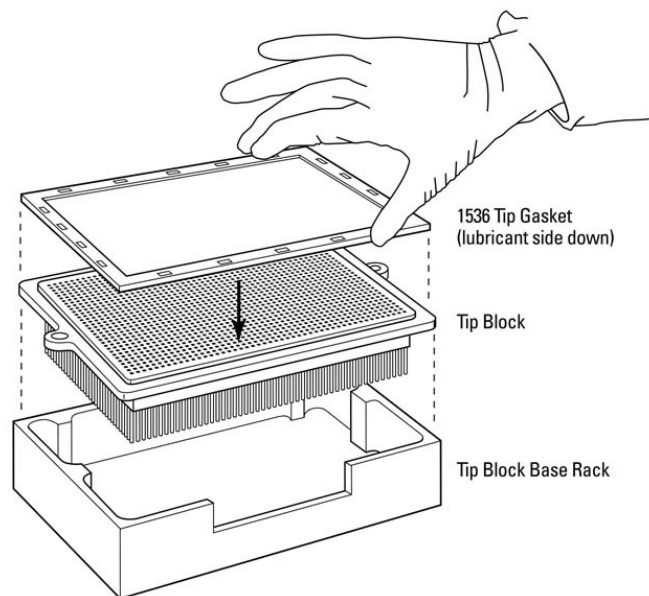
To install the 1536 tip gasket:

1. Remove the gasket from its package and discard the gasket backing.



- Place the lubricated side of the gasket down on the tip block, using the gasket recess to guide the frame position.

When aligned, the tip block and gasket are sandwiched together until the tip block grippers clamp into place.



Removing the Tip Block and Gasket

During the unloading, the tip block grippers unclamp and release the gasket and tip block into the rack assembly. When unloaded, dispose the gasket and use a new one the next time the block is loaded. Ordering information for disposable gaskets can be found in [Consumables and Accessories on page 231](#).


To load from **Position 1** for a source well plate:

- Load tip block manually and then remove the tip block before starting an experiment. When loaded, a source plate can be placed in **Source 1** in both manual and robotic modes.



Note: Sensors indicate only the presence of an object in plate position and do not indicate whether it is a source plate or tip block.

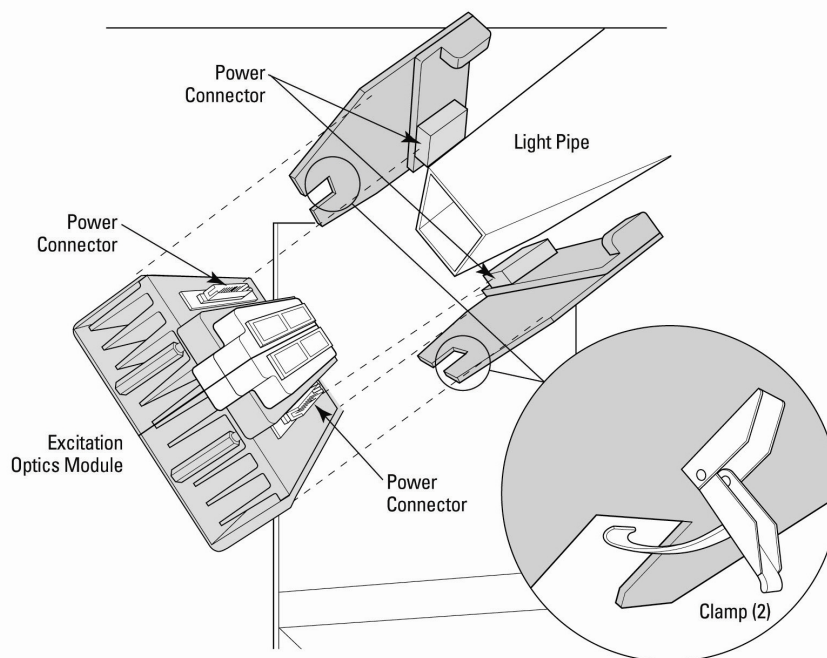
Exchanging LED Modules

 **WARNING!** During the LED exchange procedure, you have access to precision optical components. Do not touch the surfaces of the LED filters with your bare fingers or other objects because you can damage them.

Uninstalling LED Modules

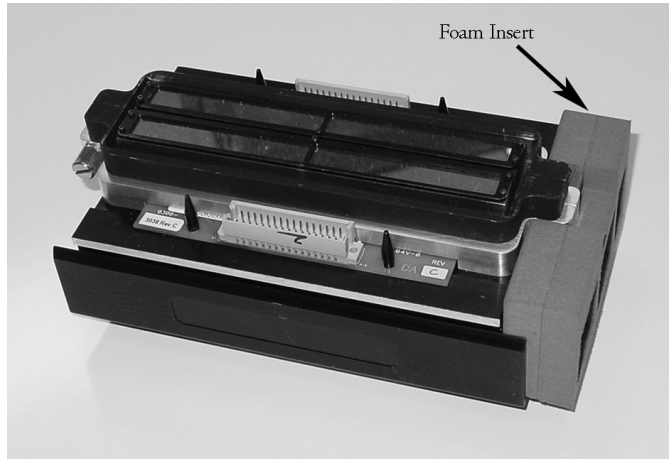
To uninstall LED modules:

1. Open the lower front door of the instrument.
When the lower door is open, power to the LEDs is disabled.
2. Loosen the two silver **Latching Clamps** securing the excitation optics module to the left light pipe.

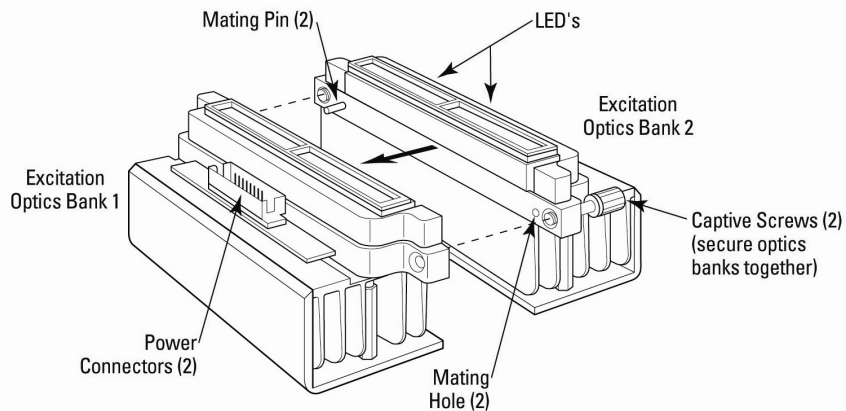


3. Pull the excitation optics module down and away from light pipe to disengage the **Power Connectors** on the LED modules from the light pipe. Be careful not to scratch the bandpass filters when pulling LED modules out of the instrument.

Remove the **Foam Insert** found on the rear side of the excitation optics module.



4. Loosen the two **Captive Thumb Screws** holding the LED modules together, and then pull the modules apart so they can be exchanged for new wavelengths.

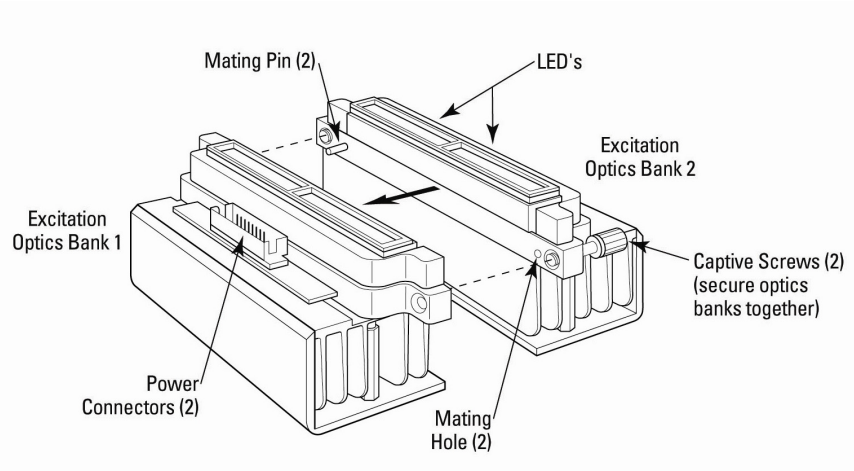


5. Repeat for the LED module mounted to the right light pipe.
6. When separated, LED modules can be exchanged using the [Installing LED Modules](#) procedure.

Installing LED Modules

To install LED modules:

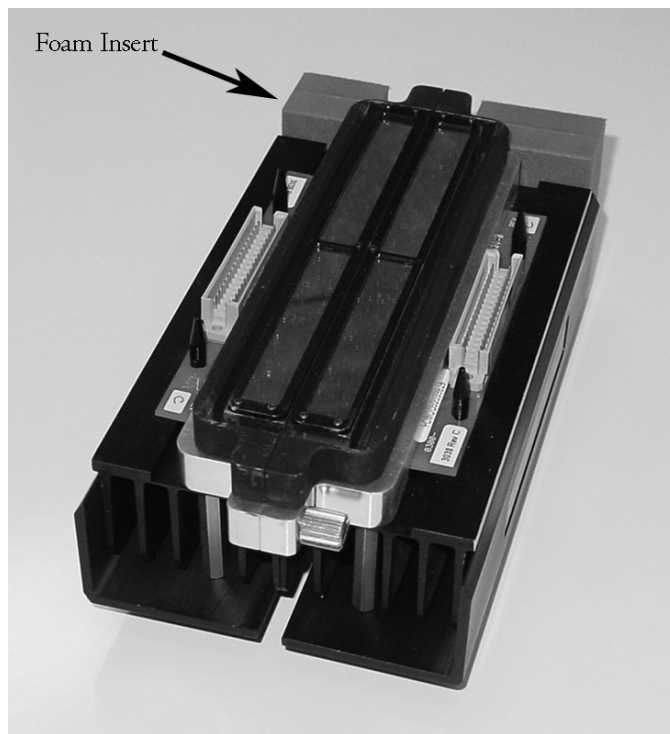
1. To create an Excitation Optics Module, you combine two LED banks of different wavelengths. Align the Mating Pins to their respective Mating Hole in the LED module that you are securing to each other.



2. When aligned, push the two LED banks of different wavelengths together and tighten the two **Captive Thumb Screws** with firm finger pressure.
3. Place the **Foam Insert** on the rear end of the excitation optics module.

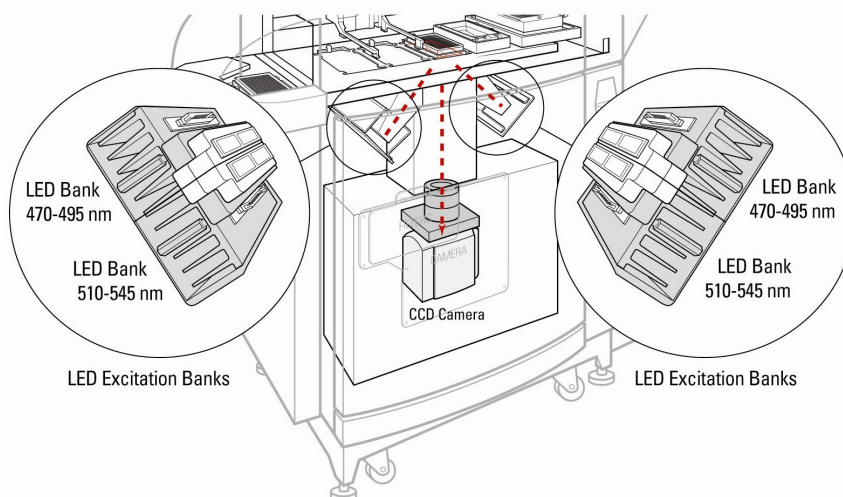


Note: Make sure that the foam piece fits inside the LED Frame.



4. Open the lower front instrument door and slide the left **Excitation Optics Module** into place. Each module is coarsely self-aligning, but observe engagement of the two **Power Connectors** and confirm they are aligned before applying mating pressure.
5. After the connectors are mated, engage and tighten the two **Latching Clamps** that secure each module to the **Light Pipe**.

- Repeat the above procedure for the right **Excitation Optics Module**, being sure the same wavelength modules occupy the corresponding position (upper/lower) on each side.



- If also changing filters, continue to [Exchanging Filters](#). If you are finished, close the lower front door.
- When the lower front door closes, and the message displays **The lower door has been opened, do you want to update optics?** Click **Yes**.
 - Filters and LEDs are identified.
 - LEDs are tested.
 - ScreenWorks Software **Instrument Status Configuration** is updated.
 - The process finishes when the progress dialog closes.

Exchanging Filters

Filters are mechanically sensed and the filter configuration can be viewed in the instrument configuration window of the FLIPR Penta System ScreenWorks Software.

If the needed filter is not available, three custom filter holders are available for purchase to use filters from an outside vendor. See [Consumables and Accessories on page 231](#). When installed, these filters appear as 1_1, 2_2, and 3_3 in the software protocol.

Uninstalling a Filter

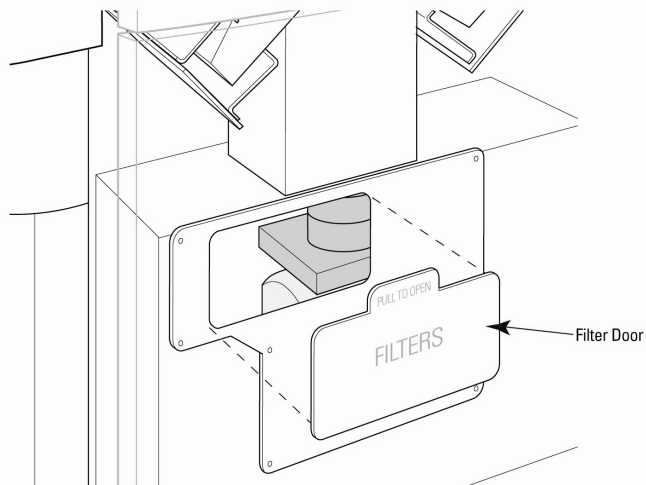
To uninstall a filter:

- Open the lower front door of the instrument.



Note: When the lower front door is open, system power is disengaged.

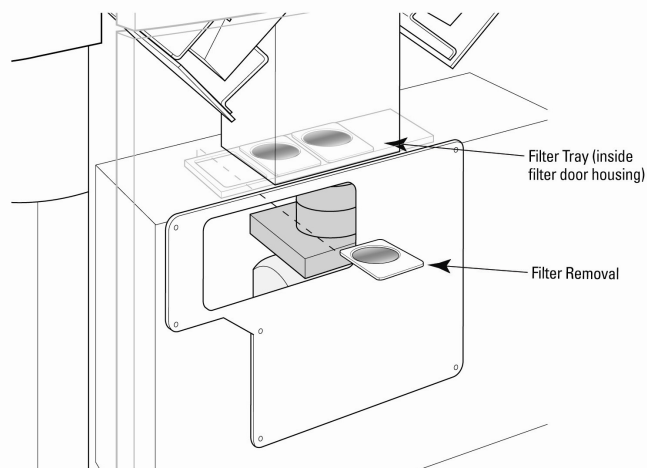
2. Remove the **Filter Cover**, located below the left LED module.



3. Using the filter tab that protrudes slightly from the filter tray, push the filter up out of its position and remove it from the instrument. Only to grasp the filter by the tab.



CAUTION! DO NOT touch the filter surface.



4. Repeat Step 3 for the remaining emission filters you want to remove.
5. After needed emission filters are removed, replace the **Filter Cover**.
6. To install new emission filters into the system, continue to [Installing an Emission Filter](#).

Installing an Emission Filter

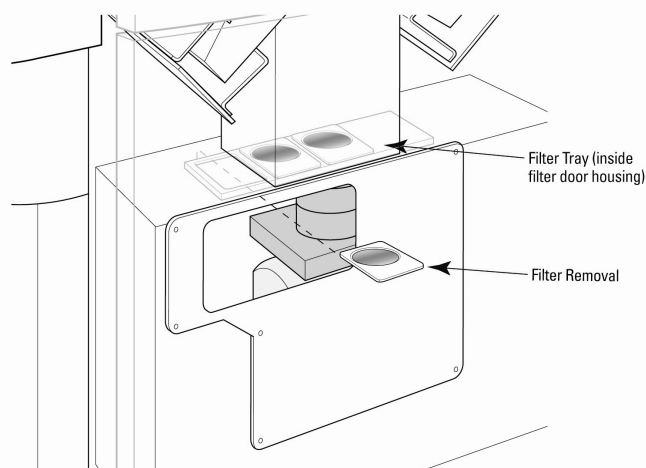
The emission filter wavelengths are automatically identified by the system and can be selected from the ScreenWorks Software menu.

To install an emission filter:

1. If needed, open the lower front door of the instrument, and remove the **Filter Cover**, located below the left LED module.
2. To install emission filters, grasp the filter you wish to install by its Tab, located where the filter table is, and place it into the appropriate filter position. Typically, this is the rightmost Position 1 in the slider. Insert the rear of the filter first, making sure it fits into the specific filter position. Push the tab down, so that the Filter Indexer clicks into the position. A notch on front of the tab should align with one of the lines on the **Filter Tray**.



CAUTION! DO NOT touch the filter surface.



3. Subsequent filters can be installed by repeating step 2 in Positions 2 and 3.



Note: If a needed filter is not available, three custom filter cassettes are available for purchase (see [Optics Consumables on page 232](#) to use filters created by an outside vendor). When installed, these filters appear as 1_1, 2_2, 3_3 in the software protocol.

4. The emission filter wavelengths are identified by the system and passed to the ScreenWorks Software where they can be selected from menus.



CAUTION! When using all 3 filter positions, do not duplicate wavelengths, because this causes a fault in the software.

5. After desired emission filters are installed, replace the **Filter Cover**.
6. When you are finished, close the lower front door.

7. When the lower front door closes, and the message displays **The lower door has been opened, do you want to update optics?** Click **Yes**.
 - Filters and LEDs are identified.
 - LEDs are tested.
 - ScreenWorks Software **Instrument Status Configuration** is updated.
 - The process finishes when the progress dialog closes.

Exchanging the Cell Reservoir

The first installation of the cell reservoir is done by a Molecular Devices certified field service engineer.

Uninstalling the Cell Reservoir

When the Cell Reservoir is configured and detected by the instrument, it is the only available choice as the Source Plate 3 selection. ScreenWorks Software displays **Cell Reservoir: OK** in the Instrument Status tab.

To use the Source 3 position for a different source plate, the Cell Reservoir must be uninstalled.

To uninstall the Cell Reservoir:

1. Remove the Reservoir Connector by loosening both Captive Thumb Screws.



2. Place the Reservoir Connector by the Tip Washer.
3. Remove the Cell Reservoir.

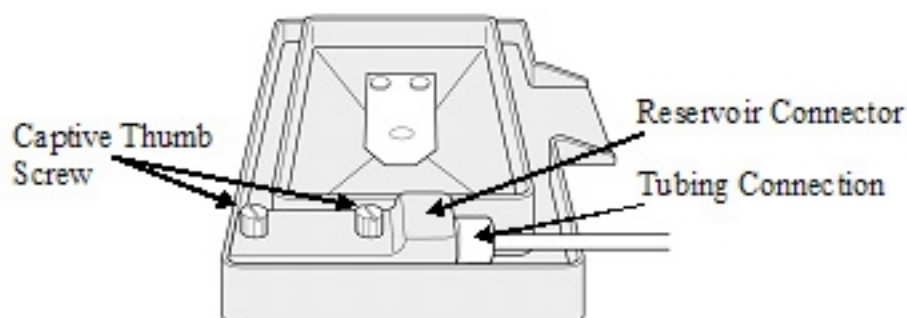
*** Tip:** You can now autoclave the Cell Reservoir.

4. Click **Instrument > Refresh Configuration**.
The Instrument Status tab displays **Cell Reservoir: NOT_OK**.
5. Select another plate for the Source Plate 3 position.

Installing the Cell Reservoir

To install the cell reservoir:

1. Ensure the tubing connection to the connector is tight.
2. Align the **Reservoir Connector** to the Cell Reservoir and tighten the two **Captive Thumb Screws**.



3. Insert the Cell Reservoir into the **Source Plate 3** position.
4. Go to the **Instrument** menu and select **Refresh Configuration**. The **Instrument Status** tab displays **Cell Reservoir: OK**.
This status indicates that only the Cell Reservoir can be selected for the **Source Plate 3** position.
5. Run a **Manual Wash** and inspect the cell reservoir area to ensure that there is no fluid leakage. See [Wash Cell Reservoir Process on page 124](#).

Exchanging the Plate Hold-Down Devices

The plate hold-down devices, part numbers 2300-1631 and 2300-1633, keep plates or tip racks from lifting after aspirating, dispensing and tip loading. Each device has an overhang that extends over the skirt of a plate to prevent upward movement when the pipettor is removed from the plate. The provided devices fit three different plate hold-down gap sizes, 2 mm, 2.5 mm, and 6 mm, which correspond to three identified plate skirt thicknesses. The 2.0 mm and 2.5 mm are incorporated in one device that rotates 90 degrees.



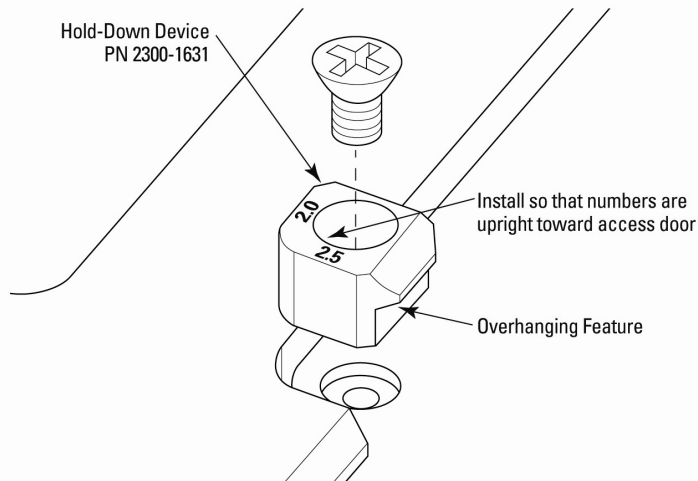
Figure 8-3: Example 2.5 mm Plate Hold-Down Device on Instrument Stage

A second plate hold-down device provided in the accessory kit, has a 6.0 mm gap. Numbers corresponding to the plate skirt thicknesses are marked on the top of the devices. The device is properly oriented for a particular flange thickness when the digits can be read as upright from the service door of the instrument.

Selecting a Plate Hold-Down Device

Most plates and tip racks use the 2.5 mm device setting. Exceptions among approved plates are the Corning/Costar plates that have a very thick 6.0 mm plate skirt and the thin 2.0 mm plate skirt of the Greiner 1536-well plates. To determine which hold-down device is most appropriate for your plate, use the lowest hold-down device that allows the plate skirt to slide under the overhanging feature freely. It is important the overhanging feature is clearly above the skirts on all plates and tip racks that are used in a particular plate position or a misalignment with the pipette tips can occur.

Test the configuration after installation.



Removing a Plate Hold-Down Device

To remove the existing plate hold-down:


1. Clear the system of plate or tips that can interfere with the position from which you are removing the plate hold-down device.
2. Use a Phillips #2 or #3 screwdriver to loosen and remove the screws from the center of the devices.
3. Remove the plate hold-down device.

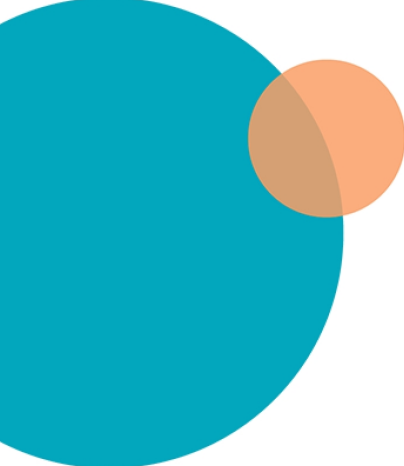
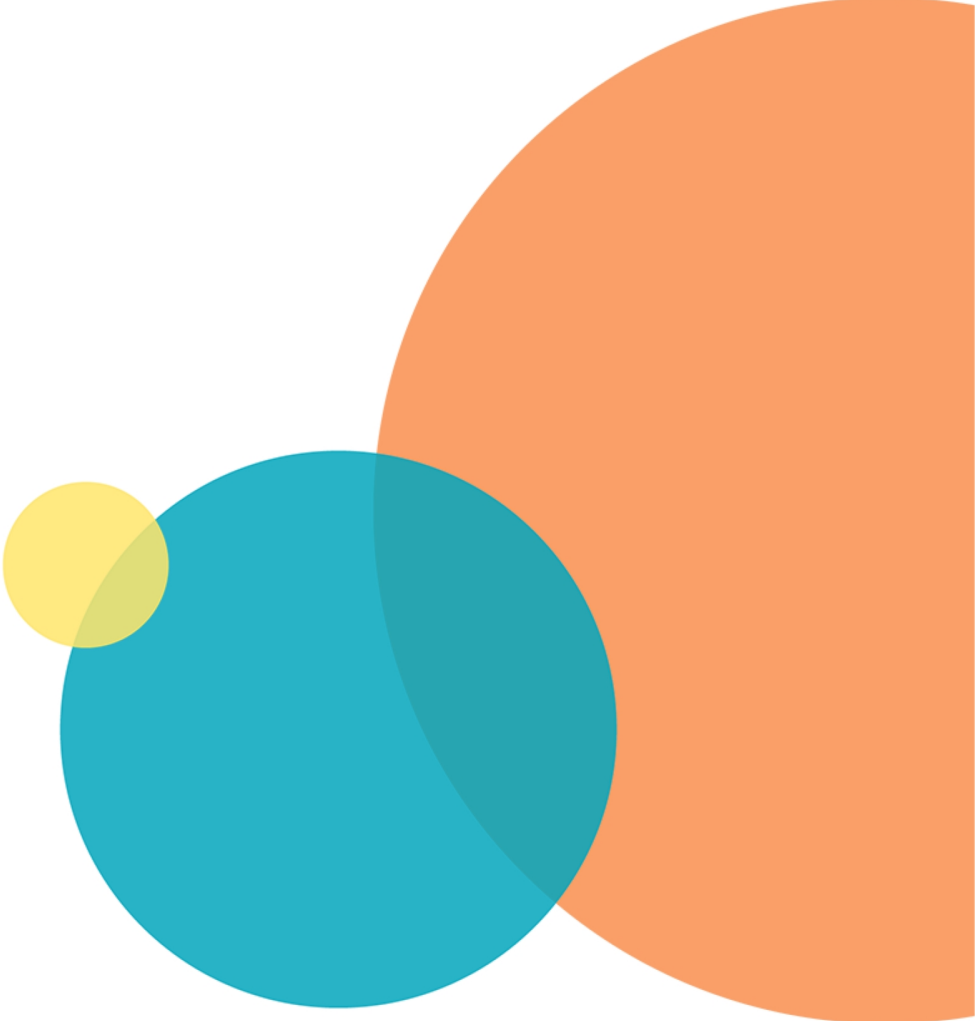
Installing a Plate Hold-Down Device

To install or reposition the plate hold-down:

1. Place the device in its proper orientation so that the numbers appear upright when viewed looking through the main user access door.

2. Insert one of the flat head screws that was originally securing the device or one included in the accessory kit through the hole in the device and into the screw threads on the left side of the plate-retaining pockets. Tighten the screw firmly. The device is self locating, therefore no additional alignment is necessary.
3. If plates or tip blocks with different flange heights are to be used in the same position, use the device specified for the largest flange thickness.
4. Test the plates or tip blocks to be used by inserting them into the plate pockets at the location in which they are used.
5. Observe that when the plate is released it is pushed to the left under the retaining device. It should come to rest against the edge locating surfaces on the left and nearest the door.
6. Pull up on the plate gently.

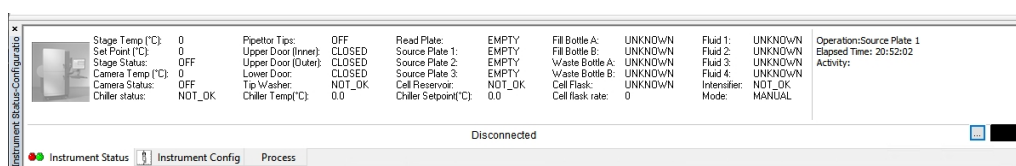
 **Tip:** A definite retaining force should be felt. A sideways movement during lifting can cause the plate to slide out from under the retaining device and be lifted upward without retention. This does not happen during instrument operation but might require attention to be avoid when lifting a plate by hand.



This chapter contains tables of symptoms and solutions to help you diagnose and repair problems with the FLIPR Penta System. In some cases, you need to contact Molecular Devices Technical Support. See [Obtaining Support on page 226](#).

Instrument Status Colors

ScreenWorks Software is designed to handle most errors that occur when running the FLIPR Penta System. The status of the system is reported in the **Instrument Status** panel. For details, see [Software Status Tabs on page 21](#).



Depending on the state of the instrument, a color icon displays at the bottom right of the **Instrument Status** window.

| If the Instrument Status Color is... | Then there is... |
|--------------------------------------|---|
| Green | Proper and open communication. |
| Yellow | Minor communication error. Clear error by selecting Clear Error from the Instrument menu. |
| Red | Instrument failure. Instrument must be reset by selecting Reset from the Instrument menu. |
| Black | No instrument communication available. This is typically shown when working in Offline mode. |

Troubleshooting Start-Up

Table 9-1: Start-Up Error Messages

| Symptom or Error Message | Possible Causes | Solutions |
|--|--|---|
| "Connecting to the instrument, please wait. Will enter offline if the instrument is not connected or if this dialog is closed. Dialog may not close immediately on pressing the close button while connection is being attempted." | ScreenWorks Software is waiting for the FLIPR Penta System to complete initialization prior to connecting. | Wait until Assay Finished (Unlocked) light is the only light illuminated on the front instrument panel before the software and instrument connect. |
| "Camera chiller subsystem not functional. Please contact Technical Support. Can't communicate with chiller." | Instrument is not communicating with the chiller. | Make sure the chiller is turned ON. Contact Technical Support. See Obtaining Support on page 226 . |

Table 9-1: Start-Up Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|--|---|--|
| 409: "Initialization halted because tips are loaded. Please verify that the appropriate tip rack (96 or 384 tip rack, 1536 tip block, or 384, 1536 pin tool holder) is loaded and select RESET". | Tips are on the pipettor head and need to be unloaded during instrument initialization/reset. | Place tip rack in source 1/tip loading position to unload tips and select Reset from the Instrument menu. |
| 410: "Initialization halted. Either the upper or lower door is open. Please shut both doors and select RESET". | The door is open during instrument initialization/reset. | Close the upper and lower doors and select Reset from the Instrument menu. |
| 411: "Front Panel Interrupt button was pushed while initializing. Initializing was not completed. Select Reset when ready to continue." | The front panel interrupt button was pushed, stopping instrument initialization. | Select Reset from the Instrument menu. |
| "Disconnected" is displayed in the bottom of the Instrument Status Window . | ScreenWorks Software is in Offline Mode . | Select Go Online from the Instrument drop-down menu. |
| | Multiple software copies are open. | Close all versions of ScreenWorks Software before starting software. |
| | Instrument timed-out. | Select Go Online from the Instrument drop-down menu. |
| Instrument Configuration does not show desired components installed on instrument. | Instrument does not have proper components installed. | Install proper components and restart system. |
| | ScreenWorks Software is in Offline Mode . | Select Go Online from the Instrument drop-down menu. |
| "Camera Temp" not in its operating temperature range EMCCD camera: -70°C (-94°F) ± 2°C HS EMCCD camera: -70°C (-94°F) ± 2°C ICCD camera: -20°C (-4°F) ± 5°C | Camera is not cooled down to temperature. | Allow 10 minutes for camera temperature to cool. Make sure the chiller is working properly. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| "Chamber Temp" not at desired temperature. | Heated stage is not warming to desired temperature. | Allow 15 minutes for the chamber to warm to the desired temperature. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| | Temperature control is not ON. | Turn on heated stage by selecting Temperature Control ON from the Instrument > Manual Operation drop-down menu. |

Table 9-1: Start-Up Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|---|---|--|
| Pipettor head is in the plate after the instrument initializes. | Instrument did not initialize properly. | Select Reset from the Instrument drop-down menu. Contact Technical Support. See Obtaining Support on page 226 . |

General Troubleshooting

Table 9-2: General System Error Messages

| Symptom or Error Message | Possible Causes | Solutions |
|--|--|--|
| 201: "Upper door open". | Upper door (outer) is not closed. | Close outer (upper) door. |
| 202: "Both upper and lower doors open". | Both upper (outer) and lower doors are not closed. | Close outer (upper) and lower doors. |
| 203: "Lower door open". | Lower door (outer) is not closed. | Close lower door. |
| 204: "Read plate position empty". | No plate is present in the Read position. | Place appropriate plate in the Read position. |
| 205: "Source 1 plate position empty". | No plate is present in the Source 1 position. | Place appropriate plate in the Source 1 position. |
| 206: "Source 2 plate position empty". | No plate is present in the Source 2 position. | Place appropriate plate in the Source 2 position. |
| 207: "Source 3 plate position empty". | No plate is present in the Source position. | Place appropriate plate in the Source 3. |
| 229: "Attempting to read a source only plate. Please assign a read plate to the read position, or if it is a read plate, go to TOOLS, PLATE LIBRARY and define the mask positions using the appropriate yellow plate." | The plate selected for the read position cannot be read because no mask has been identified. | Convert from a source only to a read plate using the plate library. |
| 100: "Command not found." | Input command is not valid. See appended string on the error message for details. | Contact Technical Support. See Obtaining Support on page 226 . |
| 101: "Invalid argument." | Input argument is not valid. See appended string on the error message for details. | Contact Technical Support. See Obtaining Support on page 226 . |

Table 9-2: General System Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|---|--|---|
| 102: "Invalid number of arguments." | Number of input arguments is not valid. See appended string on the error message for details. | Contact Technical Support. See Obtaining Support on page 226 . |
| 103: "Invalid start of command." | Start of command is not valid. See appended string on the error message for details. | Contact Technical Support. See Obtaining Support on page 226 . |
| 104: "Command invalid, system busy." | The system cannot act upon the command, as the system is busy. See appended string on the error message for details. | Contact Technical Support. See Obtaining Support on page 226 . |
| "Data buffer full". | The data buffer is full. See appended string on the error message for details. | Contact Technical Support. See Obtaining Support on page 226 . |
| 106: "Option not installed, command unavailable". | Pipettor head, LED, Emission Filter or FLIPR Cycler is not installed. | Install option you want to use. |
| | | Change protocol to use available instrument options. |
| 107: "Command failed". | General failure. | Contact Technical Support. See Obtaining Support on page 226 . |
| 108: "Command failed, system in an error state". | Minor communication error occurred. | Select Clear Error command from Instrument drop-down menu. |
| 109: "Command failed, system not initialized " | Communication failed between software and instrument. | Select Reset from the Instrument drop-down menu. |
| | | Turn the instrument power switch OFF then ON to cycle power. |
| "Password does not match". | User password is incorrect. | Contact your instrument administrator for the appropriate user name and password. |
| "Input not accepted". | Input value is not valid (e.g., select greater than 16 time sequence files in the Export menu). See appended string on the error message for details. | Change the input value so it is within the specified range. |

Table 9-2: General System Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|--|--|---|
| 302: "Temperature system not functional. Please contact Technical Support. You will not be able to run protocols with temperature control until it is repaired." | Can't read the temperature sensors for the heated stage. | Select Reset command from the Instrument drop-down menu to reset instrument to see if communication can be reconnected. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| Data files cannot be found | | Use Windows' Find utility to search for file names and determine where files were saved. Default location for ScreenWorks files is C:\Program Files\Molecular Devices\ScreenWorks\Data . |
| Hard drive fills up with data files. | Files unnecessarily saved with *.tif file. | Discard unwanted *.tif files. |
| | Data not archived. | Create free space in the hard drive. |
| "Image" option in Experiment window is grayed out. | Save images not selected in Read with TF process. | Save subsequent data with Save images selected in protocol. |
| | *.tif files not found (deleted or moved) in data folder. | *.tif files must be saved in same folder as data to retain Image button option. |
| 309: "Unable to modify or delete selected plate. It does not exist." | Plate does not exist. | Contact Technical Support. See Obtaining Support on page 226 . |
| 401: "Input/Output PCB not functional. Please contact Technical Support. You will not be able to run protocols until it is repaired." | Input/Output PCB not functional. | Contact Technical Support. See Obtaining Support on page 226 . |
| 402: "Motion reports a fatal error. Please select Reset. If this error repeats, please contact Technical Support." | Motion reports a fatal error. | Contact Technical Support. See Obtaining Support on page 226 . |
| 403: "Motion timeout waiting for done. Please select Reset. If this error repeats, please contact Technical Support." | Instrument motion timed out prior to receiving "done" command from instrument. | Select Reset from the Instrument menu. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| 404: "Motion timeout waiting for echo. Please select Reset. If this error repeats, please contact Technical Support." | Instrument motion timed out while waiting for an echo. | Select Reset from the Instrument menu. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |

Table 9-2: General System Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|---|---|--|
| 405: "Motion communication fault. Please select Reset. If this error repeats, please contact Technical Support." | Motion communication fault. | Select Reset from the Instrument menu. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| 406: "Invalid DONE response from motion. Please select Reset. If this error repeats, please contact Technical Support." | Invalid DONE response from motion. | Turn the instrument off. Wait 15 seconds. Turn instrument on. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| 407: "LCB read/write test failed. Please select Reset. If this error repeats, please contact Technical Support." | LCB read/write test failed. | Select Reset from the Instrument menu. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| 408: "Plate format file fault. Please contact Technical Support. You will be unable to use this plate until repaired." | Plate format file fault. | Select Reset from the Instrument menu. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| ScreenWorks indicates the user-defined name of the data file is invalid. | Illegal characters were used in the file name. | Eliminate illegal characters from userdefined name (for example, " , ' : ? ;). |
| | User-defined name includes too many characters. | Use 25 characters or less in the user-defined name. |
| Saturation detected. | One or more wells in the microplate saturated the camera. | A saturation warning in either the signal test dialog or on a plate prevent you from reading the plate. It is at the discretion of the user to decide whether to address the saturation warning or proceed forward. If multiple wells within a plate are saturating the camera, decrease the Excitation Intensity , Exposure Time , or Camera Gain to prevent saturation. |

Troubleshooting the Pipettor

Table 9-3: Pipettor Error Messages

| Symptom or Error Message | Possible Causes | Solutions |
|---------------------------------|--|--|
| Cannot uninstall pipettor head. | The D-Axis Knurled Nut is caught on the D-Axis receiver. | Grasp the D-Axis Knurled Nut receiver and press downward to give the pipettor head additional clearance from the D-Axis Knurled Nut. |
| | Pipettor head is caught on the alignment pins. | Lift the pipettor head up prior to moving left to lift the pipettor head off of the pipettor mount alignment pins. |

Table 9-3: Pipettor Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|--|---|--|
| Cannot install pipettor head. | The Head Clamp Knob is not tightening all the way down. | Slide the pipettor head all the way to the right over the pipettor mount. When installed properly on the alignment pins, the pipettor head should not slide left without lifting. |
| | D-Axis Knurled Nut does not reach receiver threads. | Lift D-axis receiver to engage threads and then tighten the D-Axis Knurled Nut. |
| 231: "Pipettor head not detected. Please verify proper installation and select Reset." | No pipettor head is present on pipettor mount. | Install appropriate pipettor onto instrument. |
| | Pipettor head is not installed properly. | Head clamp knob may not be secured tightly for instrument to identify pipettor head. Reinstall pipettor head. |
| | The open protocol may not use the pipettor on the instrument. | Install appropriate pipettor onto instrument. Change plate format used in the software to match the pipettor head. |
| 208: "Unable to load tips. Tips are already on pipettor." | Tips are already present on pipettor. | If you would like to change tips, select Unload Tips from the Instrument > Manual Operation drop-down menu. Once tips are exchanged, select Load Tips from the same drop-down menu. |
| | | If want to keep the present tips on, clear error and proceed with your experiment. |
| 209: "Unable to unload tips. There are no tips loaded on the pipettor." | No tips are present on the pipettor head. | Clear error. |
| 233: Unable to dispense. No tips loaded on pipettor." | No tips are present on the pipettor head. | Load tips manually on the pipettor head using the Load Tips command which is found in the Instrument > Manual Operation drop-down menu. |
| | | Change protocol to include loading of tips at the beginning of the experiment. |
| 234: "Unable to aspirate. No tips loaded on pipettor." | No tips are present on the pipettor head. | Load tips manually on the pipettor head using the Load Tips command which is found in the Instrument > Manual Operation drop-down menu. |
| | | Change protocol to include loading of tips at the beginning of the experiment |

Table 9-3: Pipettor Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|--|--|--|
| 235: “Unable to place tips in well. No tips loaded on pipettor.” | No tips are present on the pipettor head. | Load tips manually on the pipettor head using the Load Tips command which is found in the Instrument > Manual Operation drop-down menu. |
| | | Change protocol to include loading of tips at the beginning of the experiment |
| 236: “Unable to mix. No tips loaded on pipettor.” | No tips are present on the pipettor head. | Load tips manually on the pipettor head using the Load Tips command which is found in the Instrument > Manual Operation drop-down menu. |
| | | Change protocol to include loading of tips at the beginning of the experiment |
| 237: “Unable to wash. No tips loaded on pipettor.” | No tips are present on the pipettor head. | Load tips manually on the pipettor head using the Load Tips command which is found in the Instrument > Manual Operation drop-down menu. |
| | | Change protocol to include loading of tips at the beginning of the experiment |
| 246: “Unable to change head. Please unload tips and select Change Head again.” | “Change Head” action cannot be performed unless tips are removed. | Unload tips manually from the pipettor head using the Unload Tips command which is found in the Instrument > Manual Operation drop-down menu. |
| Some tips are left in the tip rack after loading the tips. | Pipettor calibration failure. | Contact Technical Support. See Obtaining Support on page 226 . |
| Some tips are left on the pipettor after unloading. | Pipettor calibration failure. | Contact Technical Support. See Obtaining Support on page 226 . |
| The tip rack is attached to the tips. | Tip box failure. | Remove tip box from tips. Choose Instrument > Manual Operation > Unload Tips to unload tips to tip loading position. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| 242: “The selected read plate has fewer wells than the pipettor head. Please select a plate with a matching format.” | The pipettor head and read plate format is incompatible (e.g., 1536 pipettor head with 96-well plate). | Load the appropriate pipettor head type. |
| | | Change protocol to include read plate format that is compatible with pipettor head. |
| 243: “The selected source 1 plate has fewer wells than the pipettor head. Please select a plate with a matching format.” | The pipettor head and Source 1 plate format is incompatible (e.g., 1536 pipettor head with 96-well plate). | Load the appropriate pipettor head type. |
| | | Change protocol to include Source 1 plate format that is compatible with pipettor head. |

Table 9-3: Pipettor Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|--|--|---|
| 244: “The selected source 2 plate has fewer wells than the pipettor head. Please select a plate with a matching format.” | The pipettor head and Source 2 plate format is incompatible (e.g., 1536 pipettor head with 96-well plate). | Load the appropriate pipettor head type. |
| | | Change protocol to include Source 2 plate format that is compatible with pipettor head. |
| 245: “The selected source 3 plate has fewer wells than the pipettor head. Please select a plate with a matching format.” | The pipettor head and Source 3 plate format is incompatible (e.g., 1536 pipettor head with 96-well plate). | Load the appropriate pipettor head type. |
| | | Change protocol to include Source 3 plate format that is compatible with pipettor head. |
| “Volume is too large.” | Volume is too large for the pipettor head or plate format you are using. See appended string on the error message for details. | Reduce the volume that you are aspirating/dispensing so it is within the specified range. |
| “Volume is too small.” | Volume is too small for the pipettor head or plate format you are using. See appended string on the error message for details. | Increase the volume that you are aspirating/dispensing so it is within the specified range. |
| “Speed is too large.” | Speed is too large for the pipettor head or plate format you are using. See appended string on the error message for details. | Reduce the speed that you are aspirating/dispensing so it is within the specified range. |
| “Speed is too small.” | Speed is too small for the pipettor head or plate format you are using. See appended string on the error message for details. | Increase the speed that you are aspirating/dispensing so it is within the specified range. |
| “Height is too large.” | Height is too large for the pipettor head or plate format you are using. See appended string on the error message for details. | Reduce the height that you are aspirating/dispensing so it is within the specified range. |
| “Height is too small.” | Height is too small for the pipettor head or plate format you are using. See appended string on the error message for details. | Increase the height that you are aspirating/dispensing so it is within the specified range. |
| “Done with plate after aspiration” is not present. | FLIPR Cycler is not present on instrument. | Must install FLIPR Cycler on instrument for command to be available. |

Table 9-3: Pipettor Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|--|--|--|
| “Done with plate after dispense” is not present. | FLIPR Cycler is not present on instrument. | Must install FLIPR Cycler on instrument for command to be available. |
| | Dialog does not appear because you are dispensing to the Read plate position. | Read plate is automatically changed at the end of an experiment when in Remote Mode . |
| “Mix Fluid After Dispense” is not present. | Command is not available when “Single aspirate – Multiple dispense” fluid transfer type is selected. | Select either Single aspirate – Single dispense or Multiple aspirate – Single dispense fluid transfer types. |

Troubleshooting the Optics

Table 9-4: Optics System Error Messages

| Symptom or Error Message | Possible Causes | Solutions |
|---|---|--|
| 240: “Attempting to use a LED module that is not installed. Please install the desired LED banks.” | Specified LED module in protocol is not installed. | Install the appropriate LED banks. |
| | | Select the appropriate LED module that is already installed on the instrument. |
| 241: “Attempting to use an emission filter that is not installed. Please install the desired emission filter.” | Specified emission filter in the protocol is not installed. | Install the appropriate emission filter. |
| | | Select the appropriate emission filter that is already installed on the instrument. |
| 210: “Unable to read plate. Calibration not valid for emission/excitation pair.” | A calibration file is not available for the LED/emission filter combination selected. | Calibrate instrument for desired LED/emission filter pair. |
| 214: “The 2 upper LED banks are not identical wavelengths. Please install identical wavelengths to both upper positions.” | The upper pair of LED banks has mismatched LED module wavelengths. | Make sure LEDs of the same wavelength are located in the upper LED bank position on either side of the read plate. |
| 215: “The 2 lower LED banks are not identical wavelengths. Please install identical wavelengths to both lower positions.” | The lower pair of LED banks has mismatched LED module wavelengths. | Make sure LEDs of the same wavelength are located in the lower LED bank position on either side of the read plate. |

Table 9-4: Optics System Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|---|--|--|
| 217: "One or more LED banks not detected. Please verify all four banks firmly installed." | Cannot detect one of the LED modules. | Make sure LED module is properly connected. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| 216: "LED module configuration/position has changes. Please recalibrate." | LED module position changed from where it was flat-field calibrated. | Calibrate instrument for current LED configuration. |
| | | Place LEDs in previous calibrated position. |
| 303: "LED modules not calibrated." | None of the installed LED banks have been flat-field calibrated. | Calibrate instrument with installed LED banks. |
| | A LED inside the LED bank burned out. | Calibrate instrument with installed LED banks. |
| | | Replace LED bank if you are unable to recalibrate LED bank. |
| 306: "Flat field calibration failed. Please verify that the correct plate is loaded. Please repeat flat field calibration." | Flat Field calibration failed; see appended string on the error message for details. | Calibrate using the correct Flat Field calibration plate in the Read position. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| 218: "Unable to define plate mask. Please install 515–575 nm emission filter." | The 515–575 nm emission filter is not installed. Thus, mask alignment cannot be performed. | Install the 515–575 nm emission filter to match the calibration file. |
| | | Calibrate instrument with new emission filter. |
| 219: "Unable to define plate mask. Please install 470–495 nm LED banks." | The 470–495 nm LED banks are not installed. Thus, mask alignment cannot be performed. | Install the 470–495 nm LED banks to match the calibration file. |
| | | Calibrate instrument with new LED bank. |

Table 9-4: Optics System Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|--|---|--|
| 307: "Unable to complete plate mask definition. Incorrect number of wells found. Please verify that the correct plate is loaded and repeat." | Mask alignment failed because it didn't find the correct number of wells (e.g., must find 384 wells for a 384 well plate). | Correct read plate format must be in Read position. |
| 308: "Unable to complete plate mask definition. Found overlapping wells. Please contact Technical Support. You will not be able to run protocols using this plate until the instrument is repaired." | Plate is tilted/popped out of position. | Make sure read plate is properly aligned in Read position. |
| 310: "Unable to perform gain curve. Please insert 515–575 nm emission filter." | 515–575 nm emission filter is not installed on system. | Install the 515–575 nm emission filter prior to performing gain curve. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| 311: "Unable to perform gain curve. Please insert 470–495 nm LED banks." | 470–495 nm LED banks are not installed on system. | Install the 470–495 nm LED banks prior to performing gain curve. |
| | | Contact Technical Support. See Obtaining Support on page 226 . |
| 304: "Gain calibration failed. Please contact Technical Support." | | Contact Technical Support. See Obtaining Support on page 226 . |
| Excitation/Emission wavelength is not available in the "edit read mode" dialog. | Instrument was not calibrated for the excitation/emission pair for which you are looking. | Calibrate instrument for the Excitation/Emission pair for which you are looking. |
| "Exposure time is too large." | The exposure time for the camera is too large. See appended string on the error message for details. | Decrease the exposure time so it is within the specified range. |
| "Exposure time is too small." | The exposure time for the camera is too small. See appended string on the error message for details. | Increase the exposure time so it is within the specified range. |
| 200: "Read interval too short." | Read interval is shorter than the sum of exposure lengths for each read interval in addition to the camera processing time. | Set a longer read interval. |

Table 9-4: Optics System Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|--|--|---|
| "First interval is too small." | The time for the first interval is too short. Interval time must always be longer than the sum of all selected read modes plus the computer processing time. See appended string on the error message for details. | Increase the read interval so it is within the specified range. |
| "First interval is too large." | The time for the first interval is too large. See appended string on the error message for details. | Decrease the read interval so it is within the specified range. |
| "First interval number of reads is too large." | The number of reads in the first interval is too large. See appended string on the error message for details. | Decrease the number of reads so it is within the specified range. |
| "Number of reads before dispense are too large." | Number of reads prior to dispense are larger than the number of reads in the first interval. See appended string on the error message for details. | Decrease the number of reads before dispense so it is within the specified range. |
| 238: "Protocol requests greater than 800 reads. This is not permitted." | Sum of all reads for all read actions and configurations must not exceed 800. | Reduce the number of reads in the experiment to less than 800 for all combined read modes. |
| 239: "Protocol requests greater than 50 images. This is not permitted." | Sum of all raw images for all read actions and configurations must not exceed 50. | Reduce the number of raw images in the experiment to less than 50 for all combined dispenses. |
| 400: "Camera is not functional. Please contact Technical Support. You will not be able to run protocols until it is repaired." | Camera is not functional. | Contact Technical Support. See Obtaining Support on page 226 . |

Troubleshooting the Yellow Plate

Table 9-5: Yellow Plate Testing Error Messages

| Symptom or Error Message | Possible Causes | Solutions |
|---|--|--|
| "Saturation detected, data may be invalid." | Incorrect settings. | Set excitation intensity to 80%, exposure length to 0.1 s and for EMCCD gain to 80, for ICCD gate to 3%. If condition repeats, lower excitation intensity. |
| | Incorrect LED and/or emission filter choice. | Check to make sure the appropriate LED and emission filters are selected. |
| | Emission optics are dirty. | Make sure no dust is on emission filters. |
| | Bottom of Read plate is dirty. | Make sure no dust or fluorescent compounds are on the bottom of the Read plate. |
| Wells are cut off in Image Viewer. | Plate not properly placed in Read position | Make sure plate is properly placed and indexed in the Read position. |
| | Dirty/misaligned optics. | Contact Technical Support. See Obtaining Support on page 226 . |
| Fluorescence counts are 0. | LED power is off. | Increase excitation intensity of LEDs. |
| | | Make sure Excitation/Emission Wavelength in the read mode does not have NONE for the excitation wavelength. |
| | Camera failure. | Contact Technical Support. See Obtaining Support on page 226 . |
| | Signal test was not initiated. | Select Test Signal in the Signal Test dialog. |

Table 9-5: Yellow Plate Testing Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|---|---|--|
| Standard deviation for the signal test is greater than 5% and fluorescence counts may or may not be in the correct range. | Instrument was calibrated with yellow boat. | If instrument was calibrated using the yellow boat, the standard deviation can be 5% or less. |
| | Camera not fully cooled down. | Wait approximately 10 minutes for the camera temperature to cool down. |
| | Optical correction out of specifications | Perform Flat-Field calibration. |
| | Plate is not aligned properly in Read position. | Make sure plate is properly aligned and indexed in the Read position. |
| | Optics is dirty or failing. | Contact Technical Support. See Obtaining Support on page 226 . |
| | Calibration plate failure— plate is dirty or scratched. | Clean calibration plate with lens paper. DO NOT use lab wipes. If cleaning does not work, replace the calibration plate. |

Troubleshooting the Tip Washer

Table 9-6: Tip Washer System Error Messages

| Symptom or Error Message | Possible Causes | Solutions |
|--|---|--|
| 221: "Tip wash reservoir format (96/384/1536) does not match pipettor head. Please install matching reservoir and select Reset." | Pipettor head and tip washer reservoir top do not match (e.g., a 384 Pipettor head with a 96-well reservoir top). | Install appropriate pipettor head to match tip wash reservoir top. |
| | | Install appropriate tip wash reservoir top to match pipettor head. |
| 222: "Tip wash reservoir top not detected. Please install matching reservoir and select Reset." | No reservoir top is installed. | Install appropriate tip wash reservoir top. |
| | Tip wash reservoir top is not properly installed. | Tip wash reservoir top connectors may not be properly connected. |
| 223: "Check Waste Bottle A - sensor indicates full." | Waste bottle A is full with tip wash solvent. | Empty tip wash solvent from waste bottle A. |
| 224: "Check Waste Bottle B - sensor indicates full." | Waste bottle B is full with tip wash solvent. | Empty tip wash solvent from waste bottle B. |
| 225: "Check TipWasher - sensor indicates overflow." | Waste bottles are full. | Empty tip wash solvent from waste bottles A and B and select Reset . |
| | Fluid level sensor malfunction. | Contact Technical Support. See Obtaining Support on page 226 . |

Table 9-6: Tip Washer System Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|---|---|---|
| 226: "TipWasher: Check Wash Fluid Bottle A - fill sensor dry." | Fill Bottle A is empty. | Place tip wash solvent into Fill Bottle A. |
| 227: "TipWasher: Check Wash Fluid Bottle B - fill sensor dry." | Fill Bottle B is empty. | Place tip wash solvent into Fill Bottle B. |
| 250: "TipWasher: Check Wash fluid. Fill sensor dry during fill after prime." | The selected wash bottle might be running low on fluid. | Check the wash bottles. |
| 251: "TipWasher: Check Wash fluid. Fill sensor dry after fill." | The selected wash bottle might be running low on fluid. | Check the wash bottles. |
| 252: "TipWasher: Check Wash Fluid Bottles A and B - sensor indicates empty." | Both of the wash bottles are detected empty. | Add fluid to both wash bottles. |
| 254: "TipWasher: Check tubing at Fill Pump - fill sensor wet after draining line." | Fluid sensor is detecting fluid when it should not. | Check for any unconnected tubing. Contact Technical Support. See Obtaining Support on page 226 . |
| 301: "Tip wash module not functional. Please contact Technical Support. You will not be able to run protocols with tip washing until it is repaired." | Can't communicate with the Tip Washer. | Select Reset command from the Instrument drop-down menu to reset instrument to see if communication can be reconnected. |

Troubleshooting the Cell Reservoir

Table 9-7: Cell Reservoir System Error Messages

| Symptom or Error Message | Possible Causes | Solutions |
|--|--|---|
| 255 "CellReservoir: Reservoir not detected. Please install it at Source 3 position and select Refresh Configuration from Instrument menu bar." | Instrument cannot detect the cell reservoir. | Check to make sure the Cell Reservoir is installed. |
| | | Check the connection of the reservoir sensor inside the instrument. |
| 256 "CellReservoir: prime sensor dry during prime/fill. Please check Fluid Bottle 1". | Fluid sensor is not detecting fluid when pumping from Fluid 1. | Check to make sure that Fluid 1 bottle is not empty. |
| 257 "CellReservoir: prime sensor dry during prime/fill. Please check Fluid Bottle 2". | Fluid sensor is not detecting fluid when pumping from Fluid 2. | Check to make sure that Fluid 2 bottle is not empty. |
| 258 "CellReservoir: prime sensor dry during prime/fill. Please check Fluid Bottle 3". | Fluid sensor is not detecting fluid when pumping from Fluid 3. | Check to make sure that Fluid 3 bottle is not empty. |

Table 9-7: Cell Reservoir System Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|---|--|--|
| 259 "CellReservoir: prime sensor dry during prime/fill. Please check Fluid Bottle 4". | Fluid sensor is not detecting fluid when pumping from Fluid 4. | Check to make sure that Fluid 4 bottle is not empty. |

Troubleshooting Data

Table 9-8: Data Processing Error Messages

| Symptom or Error Message | Possible Causes | Solutions |
|--|--|---|
| Negative control wells show a response. | Wash buffer components are different from the compound buffer. | Make sure wash buffer and compound buffer are the same. Check for a DMSO stimulation, if present. |
| Statistical results do not conform to the expected results. | Data analysis has not been optimally set up. | Refer to Data Processing Algorithms on page 249 . |
| Data parameters need to be reset (subtract bias, spatial uniformity correction, positive control, etc.) for each experiment. | *.vamp file parameters not set up correctly. | Select the Grouping or Correction buttons to adjust the parameters as necessary. |
| False negative wells. | Pipettor adjustment problem. | See Optimizing an Assay . |
| | Tip problem. | Contact Technical Support. See Obtaining Support on page 226 . |

Troubleshooting Robotic Integration

Table 9-9: Robotic Integration System Error Messages

| Symptom or Error Message | Possible Causes | Solutions |
|--|--|---|
| 220: "Unable to load plate. No plate on landing pad." | Plate present on plate landing pad, but plate presence was not detected. | Contact Technical Support. See Obtaining Support on page 226 . |
| | | If plate not positioned properly by robot, contact Technical Support for robot manufacturer. |
| | No plate was delivered to landing pad by 3rd party robot. | Contact Technical Support for robot manufacturer. |
| 221: "Unable to load plate. Plate already detected at requested location." | Plate already at requested position. | Unload plate currently in the desired position prior to delivering new plate. |
| | | If robot is attempting to load multiple plates to one position without retrieving plates, contact Technical Support for robot manufacturer. |

Table 9-9: Robotic Integration System Error Messages (continued)

| Symptom or Error Message | Possible Causes | Solutions |
|---|---|--|
| 222: "Unable to unload plate. Plate already detected at landing pad" | Plate already present on landing pad. | Remove plate on landing pad prior to unloading a new plate. |
| 223: "Unable to unload plate. No plate detected at requested position." | No plate present in position. | Clear error and proceed with experiment. |
| | Plate may have popped out of position. | Clear plate that is out of position prior to resuming experiment. |
| Bar code reader misreads the bar code and file name reads "Bad_bar_code". | Misplacement of bar code or incorrect type of bar code. | Bar code must be on the side of the plate where the bar code is reading. Place label as low as possible. Use labels with numbers on top. |

Obtaining Support

Molecular Devices is a leading worldwide manufacturer and distributor of analytical instrumentation, software, and reagents. We are committed to the quality of our products and to fully supporting our customers with the highest level of technical service.

Our Support website, www.moleculardevices.com/service-support, has a link to the Knowledge Base, which contains technical notes, software upgrades, safety data sheets, and other resources. If you still need assistance after consulting the Knowledge Base, you can submit a request to Molecular Devices Technical Support.

You can contact your local representative or Molecular Devices Technical Support at 800-635-5577 (North America only) or +1 408-747-1700. In Europe, call +44 (0) 118 944 8000.

Please have your instrument serial number or Work Order number, and your software version number available when you call.



WARNING! BIOHAZARD. It is your responsibility to decontaminate components of the instrument before you return parts to Molecular Devices for repair. Molecular Devices does not accept items that have not been decontaminated where it is applicable to do so. If parts are returned, they must be enclosed in a sealed plastic bag stating that the contents are safe to handle and are not contaminated.

Appendix A: Technical Specifications



The following table lists the technical specifications for the FLIPR Penta High-Throughput Cellular Screening System.

Table A-1: FLIPR Penta High-Throughput Cellular Screening System Technical Specifications

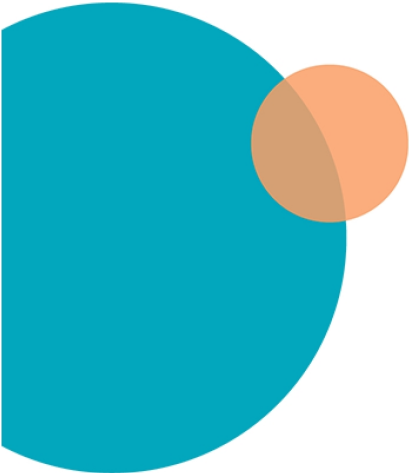
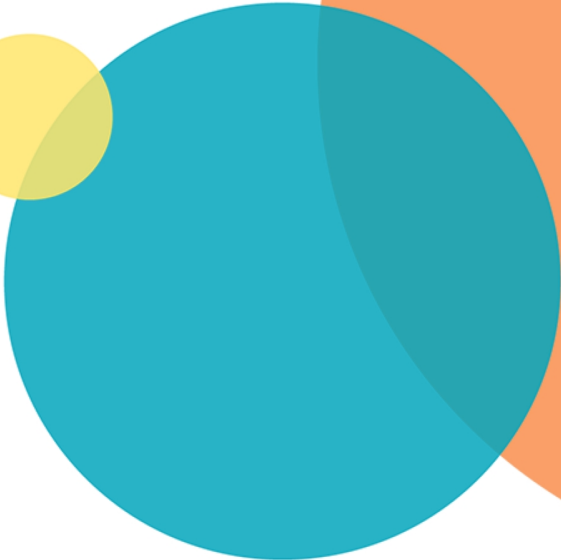
| Item | Description |
|---|--|
| Instrument Dimensions | |
| Size | Inches—39 (W) x 70.25 (H) x 27 (D) Centimeters—99 (W) x 178.44 (H) x 69 (D) |
| Weight (approx.) | 850 lbs. (386 kg) |
| Throughput | |
| Temperature control | Ambient, +25°C to 40°C |
| Throughput | 20 plates/hour (2-minute read time) |
| 96-well | 1,920 data points/hour |
| 384-well | 7,680 data points/hour |
| 1536-well | 30,720 data points/hour |
| Bar code manual mode 1 | Keyboard entry |
| Bar code manual mode 2 | Short or long side |
| EMCCD Camera Option | |
| Camera type | EMCCD |
| Detection | Fluorescence |
| Mode | Cooled CCD camera |
| Fluorescence sensitivity with $Z' \geq 0.5$ (whole plate) | 1.25 nM Fluorescein |
| Data acquisition rate | 7 Hz |
| Ratiometrics | Up to 4 EX/EM pairs per experiment |
| Data acquisition rate for ratiometric emission | 2 Hz (2EX/EM pairs) |
| HS EMCCD Camera Option | |
| Camera type | HS EMCCD |
| Detection | Luminescence and Fluorescence |
| Mode | High-Speed, High-Sensitivity CCD Camera |
| Fluorescence sensitivity with $Z' \geq 0.5$ (whole plate) | 0.625 nM Fluorescein |

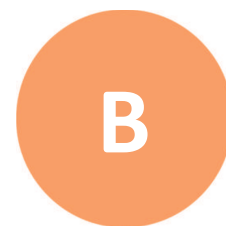
Table A-1: FLIPR Penta High-Throughput Cellular Screening System Technical Specifications (continued)

| Item | Description | |
|---|---|--------------------------|
| Luminescence sensitivity | 4.1 nM ATP (Promega Cell Titer Glow Assay) | |
| Data acquisition rate | 100 Hz | |
| Ratiometrics | Up to 4 EX/EM pairs per experiment | |
| Data acquisition rate for ratiometric emission | 2 Hz (2EX/EM pairs) | |
| ICCD Camera Option | | |
| Camera type | ICCD | |
| Detection | Luminescence and Fluorescence | |
| Mode | Intensified CCD camera | |
| Fluorescence sensitivity with $Z' \geq 0.5$ (whole plate) | 5 nM | |
| Luminescence sensitivity | 5,000 M5 AequeoZen FroZen cells in 384-well format with $Z' \geq 0.5$ | |
| Dynamic range | 4.8 log (1 measurement) 6.9 log (2 measurements) | |
| Data acquisition rate | 8 Hz | |
| Ratiometrics | Up to 4 EX/EM pairs per experiment | |
| Data acquisition rate for ratiometric emission | 2 Hz (2EX/EM pairs) | |
| Excitation LEDs and Emission Filters | | |
| Wavelength range | 335-626 nm | 400-750 nm |
| Calcium (default) | 470-495 nm | 515-575 nm |
| Membrane potential | 510-545 nm | 565-625 nm |
| Voltage sensor probes | 390-420 nm | 440-480 nm 565-625 nm |
| Fura-2 | 335-345 nm 380-390 nm | 475-535 nm |
| Instrument holds 2 pairs of LEDs, one for each side to cross-illuminate the bottom of the plate | | |
| Instrument has a 3 position filter slider that identifies the filter to the instrument | | |
| Custom filter holders are available | | |
| Fluidics | | |
| Simultaneous transfer | 96, 384 and 1536 | |
| Range-96 pipettor head | 5-200 μ L | |
| Precision @75 μ L | 3% CV | |

Table A-1: FLIPR Penta High-Throughput Cellular Screening System Technical Specifications (continued)

| Item | Description |
|---|-------------------------|
| Pipetting increments-96 pipettor head | 1 μ L |
| Range - 384-Well pipettor | 1-25 μ L |
| Precision @ 25 μ L | 4% CV |
| Pipetting increments - 384-Well pipettor | 0.5 μ L |
| Range - 1536-Well pipettor | 0.5-3 μ L |
| Precision @ 3 μ L | 6% CV |
| Precision @ 1 μ L | 10% CV |
| Pipetting increments - 1536-Well pipettor | 0.1 μ L |
| 96 pipettor head - Quadrant transfer | 384-well plate |
| 384 pipettor head - Quadrant transfer | 1536-well plate |
| Cell Suspension System | |
| Cell reservoir | Autoclavable |
| Compatible formats | 96, 384, and 1536 |
| Stir flasks available | 250 mL, 0.5 L, 1 L, 3 L |





Appendix B: Consumables and Accessories

This following catalogs consumables used with the FLIPR Penta System, including:

- [System Accessories](#)
- [Plates on page 234](#)
- [Assays on page 237](#)
- [Recommended Assay Equipment and Supplies on page 240](#)

System Accessories

Field Installations

Supplier: Molecular Devices, call +1-800-635-5577.

Table B-1: Field Installations Kits

| Item | Part Number |
|---|-------------|
| FLIPR [®] Camera Conversion Kit, ICCD or EMCCD to HS EMCCD | 5069580 |
| FLIPR [®] Cell Suspension, Field Installation Kit | 0310-5339 |
| Windows 10 Host PC with ScreenWorks [®] Software Version 5.0 | 5075062 |
| Windows 10 Host PC with ScreenWorks [®] Software Version 4.2 | 5074979 |
| FLIPR [®] FLIPR [®] Cycler | 5074848 |

Pipetting Tips and 1536 Pipetting Accessories

Supplier: Molecular Devices, call +1-800-635-5577.

Table B-2: Pipetting Consumables

| Item | Part Number |
|--|-------------|
| FLIPR [®] pipette tips, black, non-sterile, 96-well, 50 racks/case | 9000-0762 |
| FLIPR [®] pipette tips, clear, non-sterile, 96-well, 50 racks/case | 9000-0761 |
| FLIPR [®] non-sterile, black, 384-well, 50 racks/case | 9000-0764 |
| FLIPR [®] pipette tips, clear, non-sterile, 384-well, 50 racks/case | 9000-0763 |
| FLIPR [®] 1536 tip gasket, nonsterile, 40 racks/case | 9000-0746 |
| FLIPR [®] 1536 tip block | 0200-6112 |

Table B-2: Pipetting Consumables (continued)

| Item | Part Number |
|---|-------------|
| ¹ FLIPR® Pin Tool, 384 | 5075203 |
| ¹ FLIPR® Pin Tool, 1536 | 5075194 |
| 1. Contact your local Molecular Devices Sales Representative for details regarding this configure-to- order item. | |

Pipettor Heads

Supplier: Molecular Devices, call +1-800-635-5577.

Table B-3: Pipettor Head Kits

| Item | Part Number |
|---|--------------------|
| FLIPR® Pipettor Head Kit, 96 | 0200-6071 |
| FLIPR® Pipettor Head Kit, 384 | 0200-6072 |
| FLIPR® Pipettor Head Kit, 1536 | 0200-6073 |
| ¹ FLIPR® Pin Tool Kit, 384 and/or 1536 | FLIPR PIN TOOLS OC |
| 1. Contact your local Molecular Devices Sales Representative for details regarding this configure-to- order item. | |

Optics Consumables

Supplier: Molecular Devices, call +1-800-635-5577.

Table B-4: Optics Kits

| Item | Part Number |
|--|-------------|
| FLIPR® Calcium Optics Kit | 0200-6206 |
| FLIPR® Membrane Potential Optics Kit | 0200-6207 |
| FLIPR® Voltage Sensor Probes (VSP) Optics Kit | 0200-6208 |
| FLIPR® Fura-2 Optics Kit | 0200-6271 |
| FLIPR® LED Module, 335–345 nm (UV) | 0200-6272 |
| FLIPR® LED Module, 360–380 nm | 0200-6178 |
| FLIPR® LED Module, 380–390 nm (UV) | 0200-6273 |
| FLIPR® LED Module, 390–420 nm (VSP) | 0200-6135 |
| FLIPR® LED Module, 420–455 nm | 0200-6148 |
| FLIPR® LED Module, 470–495 nm (Calcium) | 0200-6128 |
| FLIPR® LED Module, 495–505 nm | 0200-6175 |
| FLIPR® LED Module, 510–545 nm (Membrane Potential) | 0200-6127 |

Table B-4: Optics Kits (continued)

| Item | Part Number |
|---|-------------|
| FLIPR® LED Module, 610–626 nm | 0200-6150 |
| FLIPR® Emission Filter, 400–460 nm | 0200-6213 |
| FLIPR® Emission Filter, 440–480 nm (VSP) | 0200-6205 |
| FLIPR® Emission Filter, 475–535 nm | 0200-6211 |
| FLIPR® Emission Filter, 515–575 nm (Calcium) | 0200-6203 |
| FLIPR® Emission Filter, 526–586 nm | 0200-6212 |
| FLIPR® Emission Filter, 565–625 nm (Membrane Potential and VSP) | 0200-6204 |
| FLIPR® Emission Filter, 646–706 nm | 0200-6214 |
| FLIPR® Custom Filter Set (3) | 0200-6221 |
| FLIPR® Single Custom Filter Holder (1) | 0200-6276 |

Cell Reservoir Consumables

Supplier: Molecular Devices, call +1-800-635-5577.

Table B-5: Cell Reservoir Consumables

| Item | Part Number |
|----------------------------------|-------------|
| FLIPR® Cell Suspension Reservoir | 0200-6222 |
| FLIPR® 250 mL Flask Assembly | 0200-6223 |
| FLIPR® 500 mL Flask Assembly | 0200-6224 |
| FLIPR® 1 L Flask Assembly | 0200-6225 |
| FLIPR® 3 L Flask Assembly | 0200-6226 |
| FLIPR® 1 L Buffer Bottle | 0200-6227 |

Plates

The following conforming plates, can be purchased from Molecular Devices.

96-Well Read Plates

Table B-6: 96-Well Read Plates

| 96-Well Read Plates | Part Number | Suggested Supplier | Phone Number |
|---|-------------|------------------------------|-----------------|
| Black, clear, tissue culture treated, sterile | 353948 | Becton Dickinson | +1-800-343-2035 |
| | 3603 | Corning/Costar | +1-800-492-1110 |
| | 655090 | Greiner (distributed by E&K) | +1-408-378-2013 |
| | 165305 | Nalge/Nunc | +1-800-766-7000 |
| Black, clear, tissue culture treated, sterile, poly-D-lysine coated | 356640 | Becton Dickinson | +1-800-343-2035 |
| | 3667 | Corning/Costar | +1-800-492-1110 |
| Black, clear, tissue culture treated, sterile, collagen coated | 356649 | Becton Dickinson | +1-800-492-1110 |
| White, clear, tissue culture treated, sterile | 353947 | Becton Dickinson | +1-800-343-2035 |
| | 3903 | Corning/Costar | +1-800-492-1110 |
| | 655098 | Greiner (distributed by E&K) | +1-408-378-2013 |
| | 165306 | Nalge/Nunc | +1-800-766-7000 |

96-Well Read Plate Masks Made by Molecular Devices

Table B-7: 96-Well Read Plate Masks Made by Molecular Devices

| 96-Well Read Plate Masks | Part Number | Plate Manufacturer |
|--------------------------|-------------|--------------------|
| 96-well slit mask | 0200-6143 | Becton Dickinson |
| | 2300-1362 | Corning/Costar |
| | 0200-6142 | Nalge/Nunc |
| | 0200-6144 | Greiner |

96-Well Source Plates

Table B-8: 96-Well Source Plates

| 96-Well Source Plates | Part Number | Suggested Supplier | Phone Number |
|---------------------------|-------------------------|---------------------------------|-----------------|
| V-bottom plate, 96-well | 353263 polypropylene | Becton Dickinson | +1-800-343-2035 |
| | 651201 polypropylene | Greiner (distributed by E&K) | +1-408-378-2013 |
| | 249944 | Nalge/Nunc | +1-800-766-7000 |
| U-bottomed plate, 96-well | 351190 polypropylene | Becton Dickinson | +1-800-343-2035 |
| | 3365 polypropylene | Corning/Costar | +1-800-492-1110 |
| | 267245 | Nalge/Nunc | +1-800-766-7000 |
| | 650201 polypropylene | Greiner (distributed by E&K) | +1-408-378-2013 |
| Deep-well plate, 96-well | 353966 polypropylene | Becton Dickinson | +1-800-343-2035 |
| | 780270 polypropylene | Greiner (distributed by E&K) | +1-408-378-2013 |
| | 278752 polypropylene | Nalge/Nunc | +1-800-766-7000 |

384-Well Read Plates

Table B-9: 384-Well Read Plates

| 384-Well Read Plates | Part Number | Suggested Supplier | Phone Number |
|---|--------------------------------------|---|-----------------|
| Black, clear, tissue culture treated, sterile | 353262 | Becton Dickinson | +1-800-343-2035 |
| | 3712 | Corning/Costar | +1-800-492-1110 |
| | 781091 | Greiner with lids (distributed by E&K) | +1-408-378-2013 |
| | MGB101-1-2 | Matrical | +1-509-343-6225 |
| | MGB101-1-3 0.17 mm-thick glass | Matrical | +1-509-343-6225 |
| | 142761 | Nalge/Nunc | +1-800-766-7000 |
| Black, clear bottom, tissue culture treated, sterile, poly-D- lysine coated | 356663 | Becton Dickinson | +1-800-343-2035 |
| | 3664 | Corning/Costar | +1-800-492-1110 |

Table B-9: 384-Well Read Plates (continued)

| 384-Well Read Plates | Part Number | Suggested Supplier | Phone Number |
|--|-------------|------------------------------|-----------------|
| Black, clear, tissue culture treated, sterile, collagen coated | 356667 | Becton Dickinson | +1-800-343-2035 |
| White, clear, tissue culture treated, sterile | 353963 | Becton Dickinson | +1-800-343-2035 |
| | 3707 | Corning/Costar | +1-800-492-1110 |
| | 781098 | Greiner (distributed by E&K) | +1-408-378-2013 |
| | 142762 | Nalge/Nunc | +1-800-766-7000 |
| Low volume black, clear, tissue culture treated, sterile | 3542 | Corning/Costar | +1-800-492-1110 |
| | 788092 | Greiner (distributed by E&K) | +1-408-378-2013 |

384-Well Source Plates

Table B-10: 384-Well Source Plates

| 384-Well Source Plates | Part Number | Suggested Supplier | Phone Number |
|------------------------|----------------------|------------------------------|-----------------|
| Flat bottom plate | 353265 polypropylene | Becton Dickinson | +1-800-343-2035 |
| | 3702 polystyrene | Corning/Costar | +1-800-492-1110 |
| | 781201 polypropylene | Greiner (distributed by E&K) | +1-408-378-2013 |
| | 265496 polystyrene | Nalge/Nunc | +1-800-766-7000 |
| U/V-bottom plate | 3657 polypropylene | Corning/Costar | +1-800-492-1110 |
| | 781280 polypropylene | Greiner (distributed by E&K) | +1-408-378-2013 |
| | 264573 polypropylene | Nalge/Nunc | +1-800-766-7000 |
| Deep-well plate | 353996 polypropylene | Becton Dickinson | +1-800-343-2035 |
| | 3965 polypropylene | Corning/Costar | +1-800-492-1110 |
| | 781270 polypropylene | Greiner (distributed by E&K) | +1-408-378-2013 |
| | MP102 polypropylene | Matrical | +1-509-343-6225 |

1536-Well Read Plates

Table B-11: 1536-Well Read Plates

| 1536-Well Read Plates | Part Number | Suggested Supplier | Phone Number |
|--|-------------|--|-----------------|
| Black/Clear, tissue culture treated, sterile | 783092 | Greiner with lids (distributed by E&K) | +1-408-378-2013 |
| White/Clear, tissue culture treated, sterile | 781098 | Greiner (distributed by E&K) | +1-408-378-2013 |

1536-Well Source Plates

Table B-12: 1536-Well Source Plates

| 1536-Well Source Plates | Part Number | Suggested Supplier | Phone Number |
|-------------------------|-------------------------|------------------------------|-----------------|
| Deep-well plate | 782270 polypropylene | Greiner (distributed by E&K) | +1-408-378-2013 |

Source Reservoirs

Table B-13: Source Reservoirs

| Source Reservoirs | Part Number | Suggested Supplier | Phone Number |
|-------------------|------------------------|--------------------|-----------------|
| Omni tray | 2428110 polystyrene | Nalge/Nunc | +1-800-786-7000 |

Assays

Calcium Flux Consumables and Accessories

Table B-14: Calcium Flux Consumables and Accessories

| Item | Part Number | Suggested Supplier | Phone Number |
|--|-------------|--------------------|------------------------------------|
| FLIPR® Calcium 6 Assay Kit: Bulk Kit | R8190 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| Explorer Kit | R8191 | | |
| Express Kit | R8195 | | |
| FLIPR® Calcium 6 QF Assay Kit: Bulk Kit | R8192 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| Explorer Kit | R8193 | | |
| Express Kit | R8196 | | |

Table B-14: Calcium Flux Consumables and Accessories (continued)

| Item | Part Number | Suggested Supplier | Phone Number |
|---|---------------------------------------|--------------------|------------------------------------|
| FLIPR® Calcium 5 Assay Kit: Bulk Kit Explorer Kit Express Kit | R8186 R8185 R8187 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| FLIPR® Calcium 4 Assay Kit: Bulk Kit Explorer Kit Express Kit | R8141 R8142 R8143 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| FLIPR® Calcium 3 Assay Kit: Bulk Kit Explorer Kit Express Kit | R8090 R8091 R8108 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| FLIPR® Cardiocytotoxicity Assay Kit: Bulk Kit Explorer Kit | R8211 (10 plates) R8210 (2 plates) | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| FLIPR® Fura-2 QBT Calcium Assay Kit: Bulk Kit Explorer Kit | R8197 R8198 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| FLIPR® Penta System Calcium Optics Kit: LED Module, 470-495 nm Emission Filter, 515-575 nm | 0200- 6206 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| FLIPR® Penta System Fura-2 Optics Kit: LED Module 333-345 nm LED Module 380-390 nm Emission Filter 475-535nm | 0200-6271 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| Hank's Balanced Salt Solution (10X stock) | 14065- 056 | Thermo Fisher | 1-800-828-6686 |
| HEPES buffer solution 1X | 9319 | Irvine Scientific | +1-800-437-5706 |
| Probenecid, crystalline | P8761 | Sigma | +1-800-325-3010 |
| Carbachol (receptormediated positive control) | C4382 | | |
| UTP, Na salt (receptormediated positive control) | U6625 | | |
| Ionomycin (positive control) | 407950 | CalBiochem | +1-800-854-3417 |

Membrane Potential Assay Kit Consumables and Accessories

Table B-15: Membrane Potential Assay Kit Consumables and Accessories

| Item | Part Number | Suggested Supplier | Phone Number |
|---|-----------------------------|--------------------|------------------------------------|
| FLIPR® Membrane Potential Assay Kit | R8034 (BLUE) R8123 (RED) | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| FLIPR® Penta System Membrane Potential Optics Kit: FLIPR® Penta System LED Module, 510–545 nm FLIPR® Penta System Emission Filter, 565–625 nm | 0200-6207 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| Hank's Balanced Salt Solution (10X stock) | 14065-056 | Thermo Fisher | 1-800-828-6686 |
| HEPES buffer solution 1X | 9319 | Irvine Scientific | +1-800-437-5706 |
| Carbachol (receptor-mediated positive control) | C4382 | Sigma | +1-800-325-3010 |

Potassium Assay Kit Consumables and Accessories

Table B-16: Potassium Assay Kit Consumables and Accessories

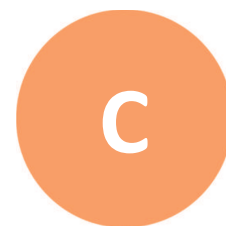
| Item | Part Number | Suggested Supplier | Phone Number |
|--|-------------------------|--------------------|------------------------------------|
| FLIPR® Potassium Assay Kit Evaluation Kit Explorer Kit Bulk Kit | R8330 R8222 R8223 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| FLIPR® Penta System Calcium Optics Kit: FLIPR® Penta System LED Module, 470-495 nm FLIPR® Penta System Emission Filter, 515-575 nm | 0200-6206 | Molecular Devices | +1-800-635-5577 +1-408-747-1700 |
| Hank's Balanced Salt Solution (10X stock) | 14065-056 | Thermo Fisher | 1-800-828-6686 |
| HEPES buffer solution 1X | 9319 | Irvine Scientific | +1-800-437-5706 |
| DMSO | D8418 | Sigma | +1-800-325-3010 |
| Probenecid | P8761 | Sigma | +1-800-325-3010 |

Recommended Assay Equipment and Supplies

The following recommended assay equipment and supplies are not available through Molecular Devices:

- 5 mL, 10 mL, 25 mL sterile serological pipettes
- Rechargeable pipettor for 2–25 mL pipettes
- Sterile tissue culture water
- Gloves
- Culture medium to grow cells
- EDTA and Trypsin/EDTA to lift cells
- Hemacytometer and counter
- Sterile test tubes 15 mL and 50 mL or smaller tubes for compounds dilutions
- 1 N NaOH solution to dissolve probenecid

Appendix C: FLIPR Penta Plate Dimensions



Note: Molecular Devices is not responsible for instrument malfunctions if the provided dimensions are incorrect.

The following dimensions derive from specifications distributed by the individual vendors.

Becton Dickinson

Table C-1: Becton Dickinson Plate Dimensions

| Plate Type | Part # | Rows | Columns | Type | Well Volume (μL) | A1 Column Offset "Y" (mm) | A1 Row Offset "X" (mm) | Bottom (MM) | Top (MM) | Well Offset (mm) |
|---------------------------------------|--------|------|---------|--------|------------------|---------------------------|------------------------|-------------|----------|------------------|
| 96-well Black/Clear | 353948 | 8 | 12 | read | 349 | 14.28 | 11.23 | 3.50 | 14.53 | 9.01 |
| 96-well Black/Clear, Collagen-Coated | 356649 | 8 | 12 | read | 349 | 14.28 | 11.23 | 3.50 | 14.53 | 9.01 |
| 96-well Black/Clear, PDL-Coated | 356640 | 8 | 12 | read | 349 | 14.28 | 11.23 | 3.50 | 14.53 | 9.01 |
| 96-well White/Clear | 353947 | 8 | 12 | read | 349 | 14.42 | 11.19 | 3.50 | 14.66 | 9.00 |
| 96-well U bottom, polypropylene | 351190 | 8 | 12 | source | 340 | 14.23 | 11.33 | 2.49 | 14.62 | 8.99 |
| 96-well V bottom, polypropylene | 353263 | 8 | 12 | source | 340 | 14.24 | 11.35 | 3.70 | 14.61 | 8.99 |
| 96-well 2mL Deep Well, polypropylene | 353966 | 8 | 12 | source | 2000 | 13.88 | 10.62 | — | 44.24 | 9.04 |
| 384-well Black/Clear | 353962 | 16 | 24 | read | 120 | 12.21 | 9.05 | 2.90 | 14.32 | 4.49 |
| 384-well Black/Clear, Collagen-Coated | 356667 | 16 | 24 | read | 120 | 12.21 | 9.05 | 2.90 | 14.32 | 4.49 |
| 384-well Black/Clear, PDL-coated | 356663 | 16 | 24 | read | 120 | 12.21 | 9.05 | 2.90 | 14.32 | 4.49 |
| 384-well White/Clear | 353963 | 16 | 24 | read | 120 | 12.26 | 9.06 | 2.90 | 14.34 | 4.49 |

Table C-1: Becton Dickinson Plate Dimensions (continued)

| Plate Type | Part # | Rows | Columns | Type | Well Volume (µL) | A1 Column Offset "Y" (mm) | A1 Row Offset "X" (mm) | Bottom (MM) | Top (MM) | Well Offset (mm) |
|------------------------------------|--------|------|---------|--------|------------------|---------------------------|------------------------|-------------|----------|------------------|
| 384-well flat bottom polypropylene | 353265 | 16 | 24 | source | 140 | 12.15 | 9.27 | 2.90 | 14.40 | 4.50 |
| 384-well Deep Well, polypropylene | 353996 | 16 | 24 | source | 400 | 11.83 | 8.83 | — | 43.69 | 4.51 |
| 1536-well Black/Clear | 353255 | 32 | 48 | read | 12 | 11.04 | 7.85 | 3.80 | 10.41 | 2.25 |
| 1536-well White/Clear | 353258 | 32 | 48 | read | 12 | 11.03 | 7.84 | 3.80 | 10.42 | 2.25 |

Corning/Costar

Table C-2: Corning/Costar Plate Dimensions

| Plate Type | Part # | Rows | Columns | Type | Well Volume (µL) | A1 Column Offset "Y" (mm) | A1 Row Offset "X" (mm) | Bottom (MM) | Top (MM) | Well Offset (mm) |
|----------------------------------|--------|------|---------|--------|------------------|---------------------------|------------------------|-------------|----------|------------------|
| 96-well Black/Clear | 3603 | 8 | 12 | read | 360 | 14.30 | 11.20 | 3.57 | 14.22 | 9.00 |
| 96-Well Black/Clear, PDL-Coated | 3667 | 8 | 12 | read | 360 | 14.30 | 11.20 | 3.57 | 14.22 | 9.00 |
| 96-well White/Clear | 3903 | 8 | 12 | read | 360 | 14.30 | 11.20 | 3.57 | 14.22 | 9.00 |
| 96-well U bottom, polypropylene | 3365 | 8 | 12 | source | 360 | 14.30 | 11.20 | 2.92 | 14.22 | 9.00 |
| 384-well Black/Clear | 3712 | 16 | 24 | read | 112 | 12.12 | 8.99 | 2.79 | 14.22 | 4.50 |
| 384-Well Black/Clear, PDL-coated | 3664 | 16 | 24 | read | 112 | 12.12 | 8.99 | 2.79 | 14.22 | 4.50 |
| 384-well White/Clear | 3707 | 16 | 24 | read | 112 | 12.12 | 8.99 | 2.79 | 14.22 | 4.50 |
| 384-well flat bottom polystyrene | 3702 | 16 | 24 | source | 112 | 12.12 | 8.99 | 2.79 | 14.22 | 4.50 |
| 384-well U bottom, polypropylene | 3657 | 16 | 24 | source | 95 | 12.12 | 8.99 | 2.65 | 14.20 | 4.50 |

Table C-2: Corning/Costar Plate Dimensions (continued)

| Plate Type | Part # | Rows | Columns | Type | Well Volume (μL) | A1 Column Offset "Y" (mm) | A1 Row Offset "X" (mm) | Bottom (MM) | Top (MM) | Well Offset (mm) |
|-------------------------------|--------|------|---------|--------|-------------------------------|---------------------------|------------------------|-------------|----------|------------------|
| 384-well block, polypropylene | 3965 | 16 | 24 | source | 180 | 12.12 | 8.99 | 2.69 | 27.80 | 4.50 |

Greiner**Table C-3: Greiner Plate Dimensions**

| Plate Type | Part # | Rows | Columns | Type | Well Volume (μL) | A1 Column Offset "Y" (mm) | A1 Row Offset "X" (mm) | Bottom (MM) | Top (MM) | Well Offset (mm) |
|--------------------------------------|--------|------|---------|--------|-------------------------------|---------------------------|------------------------|-------------|----------|------------------|
| 96-well Black/ μ Clear | 655090 | 8 | 12 | read | 392 | 14.38 | 11.24 | 3.50 | 14.40 | 9.00 |
| 96-well White/ μ Clear | 655098 | 8 | 12 | read | 392 | 14.38 | 11.24 | 3.50 | 14.40 | 9.00 |
| 96-well U bottom, polypropylene | 650201 | 8 | 12 | source | 310 | 14.38 | 11.24 | 3.70 | 14.60 | 9.00 |
| 96-well V bottom, polypropylene | 651201 | 8 | 12 | source | 340 | 14.38 | 11.24 | 3.70 | 14.60 | 9.00 |
| 96-well 2mL Deep Well, polypropylene | 780270 | 8 | 12 | source | 2420 | 14.38 | 11.24 | 3.00 | 44.00 | 9.00 |
| 384-well Black/ μ Clear | 781091 | 16 | 24 | read | 138 | 12.13 | 8.99 | 2.90 | 14.40 | 4.50 |
| 384-well White/ μ Clear | 781098 | 16 | 24 | read | 138 | 12.13 | 8.99 | 2.90 | 14.40 | 4.50 |
| 384-well flat bottom polypropylene | 781201 | 16 | 24 | source | 152 | 12.13 | 8.99 | 2.90 | 14.40 | 4.50 |
| 384-well V bottom, polypropylene | 781280 | 16 | 24 | source | 145 | 12.13 | 8.99 | 2.90 | 14.40 | 4.50 |
| 384-well Deep Well, polypropylene | 781270 | 16 | 24 | source | 240 | 12.13 | 8.99 | 2.70 | 22.00 | 4.50 |

Table C-3: Greiner Plate Dimensions (continued)

| Plate Type | Part # | Rows | Columns | Type | Well Volume (µL) | A1 Column Offset "Y" (mm) | A1 Row Offset "X" (mm) | Bottom (MM) | Top (MM) | Well Offset (mm) |
|------------------------------------|--------|------|---------|--------|------------------|---------------------------|------------------------|-------------|----------|------------------|
| 1536-well Black/µClear, Low Base | 783092 | 32 | 48 | read | 13 | 11.01 | 7.87 | 1.80 | 10.40 | 2.25 |
| 1536-well Deep Well, polypropylene | 782270 | 32 | 48 | source | 18 | 11.01 | 7.87 | 2.50 | 10.10 | 2.25 |

Matrical

Table C-4: Matrical Plate Dimensions

| Plate Type | Part # | Rows | Columns | Type | Well Volume (µL) | A1 Column Offset "Y" (mm) | A1 Row Offset "X" (mm) | Bottom (MM) | Top (MM) | Well Offset (mm) |
|----------------------------------|------------|------|---------|--------|------------------|---------------------------|------------------------|-------------|----------|------------------|
| 384-well Black/.17mm-thick glass | MGB101-1-2 | 16 | 24 | read | 120 | 12.13 | 8.99 | 2.85 | 14.40 | 4.50 |
| 384-well Black/Clear polystyrene | MGB101-1-3 | 16 | 24 | read | 120 | 12.13 | 8.99 | 2.85 | 14.40 | 4.50 |
| 384-well Deep well polypropylene | MP102 | 16 | 24 | source | 225 | 12.13 | 8.99 | 4.90 | 27.40 | 4.40 |

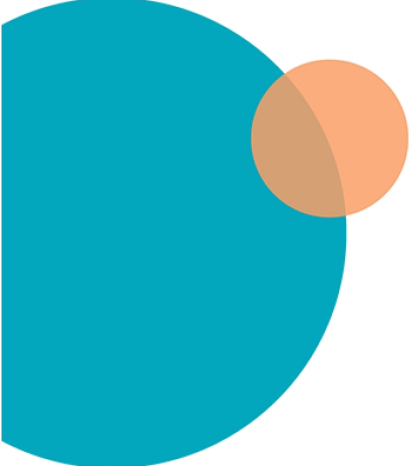
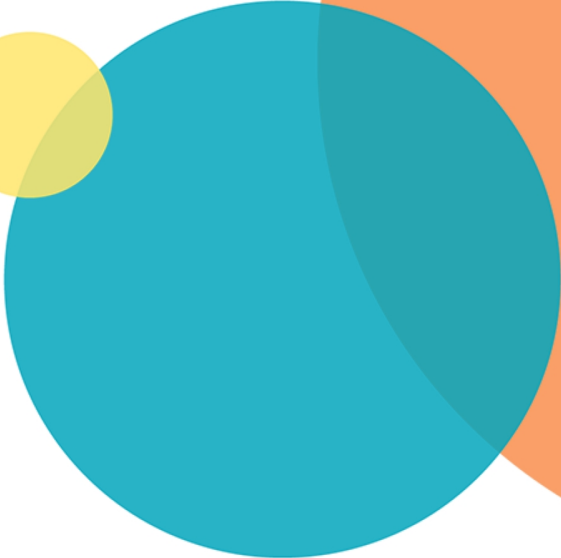
Nalge/Nunc

Table C-5: Nalge/Nunc Plate Dimensions

| Plate Type | Part # | Rows | Columns | Type | Well Volume (µL) | A1 Column Offset "Y" (mm) | A1 Row Offset "X" (mm) | Bottom (MM) | Top (MM) | Well Offset (mm) |
|---------------------|--------|------|---------|--------|------------------|---------------------------|------------------------|-------------|----------|------------------|
| 96-well Black/Clear | 165305 | 8 | 12 | read | 400 | 14.48 | 11.20 | 2.30 | 14.80 | 9.00 |
| 96-well White/Clear | 165306 | 8 | 12 | read | 400 | 14.48 | 11.20 | 2.30 | 14.80 | 9.00 |
| 96-well U bottom | 267245 | 8 | 12 | source | 500 | 14.30 | 11.30 | 2.20 | 14.30 | 9.00 |
| 96-well V bottom | 249944 | 8 | 12 | source | 450 | 14.48 | 11.20 | 3.40 | 14.40 | 9.00 |

Table C-5: Nalge/Nunc Plate Dimensions (continued)

| Plate Type | Part # | Rows | Columns | Type | Well Volume (μ L) | A1 Column Offset "Y" (mm) | A1 Row Offset "X" (mm) | Bottom (MM) | Top (MM) | Well Offset (mm) |
|--------------------------------------|--------|------|---------|--------|------------------------|---------------------------|------------------------|-------------|----------|------------------|
| 96-well 2mL Deep Well, polypropylene | 278752 | 8 | 12 | source | 2000 | 14.30 | 11.30 | 2.50 | 44.00 | 9.00 |
| 384-well Black/Clear | 142761 | 16 | 24 | read | 120 | 12.10 | 9.00 | 2.70 | 14.40 | 4.50 |
| 384-well White/Clear | 142762 | 16 | 24 | read | 120 | 12.10 | 9.00 | 2.70 | 14.40 | 4.50 |
| 384-well flat bottom polystyrene | 265196 | 16 | 24 | source | 120 | 12.10 | 9.00 | 2.70 | 14.40 | 4.50 |
| 384-well U bottom | 264573 | 16 | 24 | source | 120 | 12.10 | 9.00 | 2.70 | 14.40 | 4.50 |



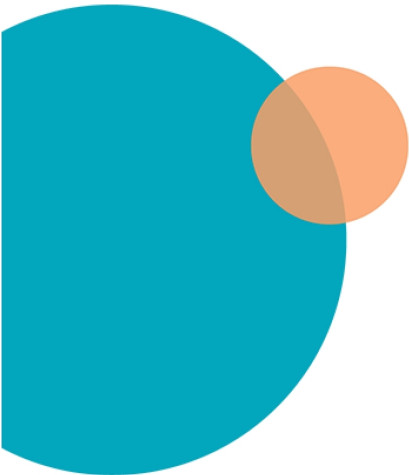
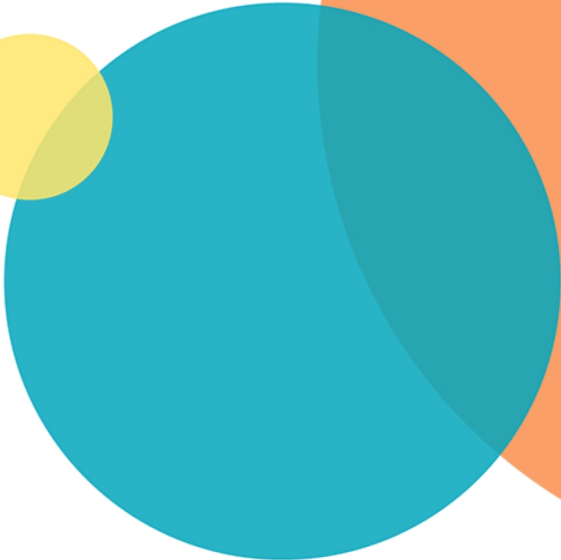


Appendix D: FLIPR Penta Instrument LED and Filter Combinations

The following are recommended LED and emission filter combinations for listed target assays.

Table D-1: FLIPR Penta Instrument LED and Filter Combinations

| Target Assay | LED (nm) | LED Part # | Emission Filter (nm) | Penta Filter Part # | Penta Kit Part # |
|--|--------------------------|------------------------|----------------------|------------------------|------------------|
| Fura-2 | 335-345 and 380-390 (UV) | 0200-6272 0200-6273 | 475-535 | 0200-6211 | 0200-6271 |
| MQAE | 360-380 (UV) | 0200-6178 | 400-460 | 0200-6213 | NA |
| Voltage Sensor Probe (VSP) | 390-420 (UV) | 0200-6135 | 440-480 | 0200-6205 | 0200-6208 |
| Tango™ GPCR Assay System (FRET) | 390-420 (UV) | 0200-6135 | 440-480 515-575 | 0200-6205 | NA |
| CFP, eCFP/YFP | 420-455 | 0200-6148 | 475-535 | 0200-6211 | NA |
| FLIPR Calcium Assay Kits | 470-495 (Blue) | 0200-6128 | 515-575 | 0200-6203 | 0200-6206 |
| Calcium, Fluo-4 | 470-495 (Blue) | 0200-6128 | 515-575 | 0200-6203 | 0200-6206 |
| Calcium, Fluo-3 | 495-505 | 0200-6175 | 526-586 | 0200-6212 | NA |
| JC-1 | 495-505 | 0200-6175 | 565-625 | 0200-6212 | NA |
| FLIPR Membrane Potential Assay Kits | 510-545 (Green) | 0200-6127 | 565-625 | 0200-6204 | 0200-6207 |
| Rhod-2 & Rhod-4 | 510-545 | 0200-6127 | 565-625 | 0200-6204 | NA |
| Alexa 633 & Bodipy | 610-626 | 0200-6150 | 646-706 | 0200-6214 | NA |
| BRET 2 luminescence assay with ICCD Camera | NA | NA | 440-480 526-586 | 0200-6205 0200-6212 | NA |
| Single Custom Filter Holder | NA | NA | NA | 0200-6276 | |
| Three Custom Filter Holders | NA | NA | NA | 0200-6221 | |





Appendix E: Data Processing Algorithms

This appendix uses results of a hypothetical experiment to describe the algorithms that can be used to adjust data. These options are all applied from the Correction dialog.

Hypothetical Experiment

Consider the results from an experiment consisting of wells A1–A9. The wells are classified into three groups:

- A1–A3: Negative Control Wells (–Ctrl). These wells contain the same dye-loaded cells and buffers as the other wells, yet are not exposed to a stimulus.
- A4–A6: Experimental Wells (Exp). These cells receive the experimental stimulus.
- A7–A9: Positive Control Wells. (+Ctrl). These cells either (1) receive a stimulus known to elicit a predetermined response or (2) demonstrate the maximal activity from an agonist.

In the following experiment, six samples, also known as pictures, readings, or frames, were taken at five second intervals. After Sample 3 was taken, buffer, with or without stimulus, was added to the wells.



Note: The longer time between readings 3 and 4. This time includes the time necessary for the 96-well pipettor to transfer liquids from the source plate to the cell plate and is automatically corrected for by the FLIPR Penta System.

The data presented is in the same format as the export data files generated using the FLIPR Penta System software. In this example, only nine wells are considered.

In [Table E-1](#), the results of the experiment are presented without Spatial Uniformity Correction, Negative Control Correction, Positive Control Scaling, or Subtract Bias Value.

Table E-1: Hypothetical Experiment

| Sample | Time | Well | | | | | | | | |
|--------|------|--------------|--------------|--------------|--------|--------|--------|--------------|--------------|--------------|
| | | A1 – Ctrl | A2 – Ctrl | A3 – Ctrl | A4 Exp | A5 Exp | A6 Exp | A7 + Ctrl | A8 + Ctrl | A9 + Ctrl |
| 1 | 0 | 8000 | 8500 | 9500 | 8200 | 9200 | 8800 | 1000 | 9500 | 8700 |
| 2 | 5 | 8200 | 8600 | 9400 | 8300 | 9000 | 9000 | 9500 | 9700 | 9000 |
| 3 | 10 | 8400 | 8500 | 9200 | 8500 | 9000 | 9100 | 9700 | 9700 | 9400 |
| 4 | 20 | 8600 | 8400 | 9200 | 45000 | 50000 | 42000 | 57000 | 52000 | 54000 |
| 5 | 25 | 8600 | 8800 | 9200 | 37000 | 42000 | 35000 | 53000 | 50000 | 50000 |
| 6 | 30 | 8800 | 8500 | 9500 | 30000 | 25000 | 29000 | 50000 | 51000 | 50000 |

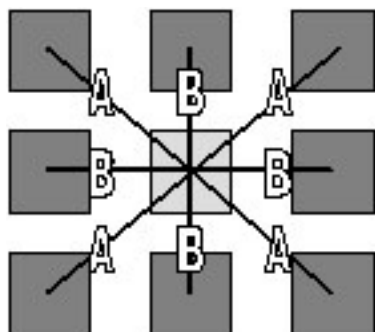
An inspection of the fluorescence counts taken at Sample 1 (time = 0) shows some inter-well variability (range 8000–10000). Variability can be the result of differences in cell density or well positions (so-called ‘edge effects’ where cells on the outside of the plate are less dense than their internal neighbors), dye loading, non-uniform illumination, buffer variations or any other effect which is constant throughout the experimental run. These variations can complicate data interpretation.

Determining Crosstalk Correction

The crosstalk correction algorithm compensates for light from one well affecting the read out from a neighboring well.

If the middle well in the drawing below was empty and the eight wells surrounding it would have signal, the middle well would show up as having a certain value based on the amount of light crossing over from each of its neighbors. The percentage of crosstalk from wells directly next to (B) the well of interest is assumed to be different from the percentage of the crosstalk from the wells diagonal (A) from the well of interest, therefore these two values are entered for the crosstalk correction. In this calculation, only the immediate neighboring wells are assumed to impact the middle well.

When the crosstalk correction is applied, the Horizontal/Vertical factor and the Diagonal Factor are applied in a matrix, which automatically subtracts the percentage of each of the neighboring well’s value from the middle well’s value. If the crosstalk correction option is selected it is the first correction applied to the data. Crosstalk correction is done in real time and no resulting data is stored. It is recommended to test this correction during assay optimization.



Determining Spatial Uniformity Correction

The spatial uniformity correction algorithm compensates, to a certain extent, for of the above-mentioned variations using a correction factor applied to all plate wells and samples.

The correction factor is derived by calculating the mean fluorescence counts of all wells at Sample 1, for example:

Table E-2: Correction Factor

| Sample | Time | Well | | | | | | | | |
|--|------|--------------|--------------|--------------|-----------|-----------|-----------|--------------|--------------|--------------|
| | | A1 – Ctrl | A2 – Ctrl | A3 – Ctrl | A4 Exp | A5 Exp | A6 Exp | A7 + Ctrl | A8 + Ctrl | A9 + Ctrl |
| 1 | 0 | 8000 | 8500 | 9500 | 8200 | 9200 | 8800 | 10000 | 9500 | 8700 |
| Mean (A1-A9) | | 8933 | | | | | | | | |
| Spatial Uniformity Correction Factor Calculation (Mean/Well) | | | | | | | | | | |
| Well- Specific Correction Factor | | | | | | | | | | |
| | | A1 | A2 | A3 | A4 | A5 | A6 | A7 | A8 | A9 |
| | | 1.12 | 1.05 | 0.94 | 1.09 | 0.95 | 1.02 | 0.89 | 0.94 | 1.03 |



Note: In actual practice, data from all 96 wells are included in the calculations, but in this example, data from only nine wells are presented, to simplify the example.

Each well-specific spatial uniformity correction factor is calculated by dividing the mean fluorescence counts of all wells by the fluorescence counts of each well (taken at Sample 1.) The table above also presents the correction factor for wells A1–A9.

All samples taken from a particular well are multiplied by their well-specific correction factor. For example, all samples from A1 are multiplied by 1.12, A2 by 1.05, and so on.

The results of applying the spatial uniformity correction factor are presented in the table below. Note the decreased variability range of wells A1–A9 in Sample 1 (8900–8976) as compared to the same data prior to applying the correction algorithm (8000–10000).

Table E-3: Spatial Uniformity Correction factor

| Sample | Time | Well | | | | | | | | |
|--------|------|--------------|--------------|-------------|--------|--------|--------|--------------|--------------|--------------|
| | | A1 – Ctrl | A2 – Ctrl | 3 – Ctrl | A4 Exp | A5 Exp | A6 Exp | A7 + Ctrl | A8 + Ctrl | A9 + Ctrl |
| 1 | 0 | 8960 | 8925 | 8930 | 8938 | 8924 | 8976 | 8900 | 8930 | 8961 |
| 2 | 5 | 9194 | 9030 | 8836 | 9047 | 8730 | 9180 | 8455 | 9118 | 9270 |
| 3 | 10 | 9408 | 8925 | 8648 | 9265 | 8730 | 9282 | 8633 | 9118 | 9682 |
| 4 | 20 | 9632 | 8820 | 8648 | 49050 | 48500 | 42840 | 50730 | 48880 | 55620 |
| 5 | 25 | 9632 | 9240 | 8648 | 40330 | 40740 | 35700 | 47170 | 47000 | 51500 |

Table E-3: Spatial Uniformity Correction factor (continued)

| Sample | Time | Well | | | | | | | | |
|--------|------|-----------|-----------|----------|--------|--------|--------|-----------|-----------|-----------|
| | | A1 – Ctrl | A2 – Ctrl | 3 – Ctrl | A4 Exp | A5 Exp | A6 Exp | A7 + Ctrl | A8 + Ctrl | A9 + Ctrl |
| 6 | 30 | 9856 | 8925 | 8930 | 32700 | 33950 | 29580 | 44500 | 47940 | 51500 |

If the spatial uniformity correction factor is applied to plates with empty wells, non dye-loaded cells, or a panel of cells containing different dyes and/or dye concentrations, the well-specific fluorescence counts are skewed by the correction factor. However, the EC₅₀ of the agonists tested are not affected.

Determining Negative Control Correction

The negative control correction algorithm corrects for changes in fluorescence that occur in all wells over the course of the experiment. Causes for these changes in fluctuations in fluorescence include dye leakage from cells, fluid addition artifacts, changes in illumination power, dye photo-bleaching, and temperature drifts.

The negative control correction factor is derived by first calculating the mean fluorescence of each of the samples of the negative control wells.

| Sample | Time | Well | | | Mean | Correction Factor |
|--------|------|-----------|-----------|----------|------|-------------------|
| | | A1 – Ctrl | A2 – Ctrl | 3 – Ctrl | | |
| 1 | 0 | 8960 | 8925 | 8930 | 8938 | 1.00 |
| 2 | 5 | 9194 | 9030 | 8836 | 9017 | 0.99 |
| 3 | 10 | 9408 | 8925 | 8648 | 8994 | 0.99 |
| 4 | 20 | 9632 | 8820 | 8648 | 9033 | 0.99 |
| 5 | 25 | 9632 | 9240 | 8648 | 9173 | 0.97 |
| 6 | 30 | 9856 | 8925 | 8930 | 9237 | 0.997 |

The mean of Sample 1 is divided by the mean of each of the samples to give the sample-specific correction factor. Each sample is multiplied by its sample-specific correction factor.

| Sample | Time | Well | | | | | | | | |
|--------|------|-----------|-----------|----------|--------|--------|--------|-----------|-----------|-----------|
| | | A1 – Ctrl | A2 – Ctrl | 3 – Ctrl | A4 Exp | A5 Exp | A6 Exp | A7 + Ctrl | A8 + Ctrl | A9 + Ctrl |
| 1 | 0 | 8960 | 8925 | 8930 | 8938 | 8924 | 8976 | 8900 | 8930 | 8961 |
| 2 | 5 | 9194 | 8940 | 8748 | 8957 | 8643 | 9088 | 8370 | 9027 | 9177 |
| 3 | 10 | 9314 | 8836 | 8562 | 9172 | 8643 | 9189 | 8547 | 9027 | 9585 |
| 4 | 20 | 9536 | 8732 | 8562 | 48560 | 48015 | 42412 | 50223 | 48391 | 55064 |
| 5 | 25 | 9343 | 8963 | 8389 | 39120 | 39518 | 34629 | 45755 | 45590 | 49955 |
| 6 | 30 | 9560 | 8657 | 8662 | 31719 | 32932 | 28693 | 43165 | 46502 | 49955 |

For example, A1 Sample 1 is multiplied by 1.00.



Note: The negative control correction factor wells contain the same cells and dyes as the experimental and positive control wells.

Determining Positive Control Scaling

The positive control algorithm compares the percent change in fluorescence counts of the positive control wells with all wells. This algorithm facilitates comparisons of results between data runs, such as different plates, by controlling factors such as cell density, cell response, laser power, or exposure time. This algorithm also makes EC₅₀ comparisons easier.

Calculate the means of all samples of the positive control wells.

| Sample | Time | Well | | | Mean | Difference from Sample 1S |
|--------|------|-----------|-----------|-----------|-------|---------------------------|
| | | A7 + Ctrl | A8 + Ctrl | A9 + Ctrl | | |
| 1 | 0 | 8900 | 8930 | 8961 | 8930 | 0 |
| 2 | 5 | 8370 | 9027 | 9177 | 8858 | -72 |
| 3 | 10 | 8547 | 9027 | 9053 | 9053 | 123 |
| 4 | 20 | 50223 | 48391 | 55064 | 51226 | 42296 |
| 5 | 25 | 45755 | 45590 | 49955 | 47100 | 38170 |
| 6 | 30 | 43165 | 46502 | 49955 | 46541 | 37610 |

The difference in fluorescence counts between Sample 1 and all of the samples is calculated. The greatest difference is determined. 100 is divided by the greatest difference in fluorescence counts to give the positive control correction factor, for example, $100/42296 = 0.0024$. All samples are multiplied by the correction factor 0.0024.

| Sample | Time | Well | | | | | | | | |
|--------|------|-----------|-----------|-----------|--------|--------|--------|-----------|-----------|-----------|
| | | A1 – Ctrl | A2 – Ctrl | A3 – Ctrl | A4 Exp | A5 Exp | A6 Exp | A7 + Ctrl | A8 + Ctrl | A9 + Ctrl |
| 1 | 0 | 21 | 21 | 21 | 21 | 21 | 21 | 21 | 21 | 21 |
| 2 | 5 | 21 | 21 | 21 | 21 | 20 | 21 | 20 | 21 | 23 |
| 3 | 10 | 22 | 21 | 20 | 22 | 20 | 22 | 20 | 21 | 23 |
| 4 | 20 | 23 | 21 | 20 | 115 | 113 | 100 | 119 | 114 | 130 |
| 5 | 25 | 22 | 21 | 20 | 92 | 93 | 82 | 108 | 108 | 118 |
| 6 | 30 | 23 | 20 | 20 | 75 | 78 | 68 | 1021 | 110 | 118 |



Note: All wells, at samples 1–3, range from 20 to 23%. The positive control wells at Sample 4 range from 114 to 130% while the negative control wells remain at around 20% throughout the experiment.

Determining Subtract Bias

The subtract bias algorithm determines the change in fluorescence over background fluorescence. The user determines the sample number chosen as background. This algorithm, by default, is always switched on and subtracts from Sample 1. The choice in the sample number to use as background depends on the assay conditions. For example, the addition of agonist can dilute the fluorescence signal in the supernatant, which would appear as a drop in fluorescence. In this case, the best choice for background is the sample taken at the bottom of this drop in fluorescence.

In our example, the percent positive fluorescence at Sample 1 is around 20% for all wells. Subtracting the background at Sample 1 would make the data easier to interpret.

Sample 1 (or any sample chosen by the operator) from each well is subtracted from the samples from the same well.

| Sample | Time | A1 – Ctrl | A2 – Ctrl | A3 – Ctrl | A4 Exp | A5 Exp | A6 Exp | A7 + Ctrl | A8 + Ctrl | A9 + Ctrl |
|--------|------|--------------|--------------|--------------|-----------|-----------|-----------|--------------|--------------|--------------|
| 1 | 0 | | | | | | | | | |
| 2 | 5 | | | | | | | | | 1 |
| 3 | 10 | 1 | | | 1 | | | | | 1 |
| 4 | 20 | 1 | | | 94 | 92 | 79 | 98 | 93 | 109 |
| 5 | 25 | 1 | | | 71 | 72 | 61 | 87 | 87 | 97 |
| 6 | 30 | 1 | | | 54 | 57 | 47 | 81 | 89 | 97 |

The results show only numbers above zero. The subtract bias algorithm places the data in a readily interpretable form. The positive control wells at Sample 4 range from 93–109% positive. In contrast, the negative control wells remain close to zero percent positive throughout the experiment. The experimental wells show two wells with about 90% activity and one with 79% activity at Sample 4.

Determining Response Over Baseline

The response over baseline correction algorithm compensates, to a certain extent, for variations within a single well.

The correction factor is derived by calculating the mean fluorescence counts in each well between the baseline start and end samples.

| Sample | Time | Well | | | | | | | | |
|--------|------|--------------|--------------|--------------|-----------|-----------|-----------|--------------|--------------|--------------|
| | | A1 – Ctrl | A2 – Ctrl | A3 – Ctrl | A4 Exp | A5 Exp | A6 Exp | A7 + Ctrl | A8 + Ctrl | A9 + Ctrl |
| 1 | 0 | 8000 | 8500 | 9500 | 8200 | 9200 | 8800 | 10000 | 9500 | 8700 |
| 2 | 5 | 8200 | 8600 | 9400 | 8300 | 9000 | 9000 | 9500 | 9700 | 9000 |
| 3 | 10 | 8400 | 8500 | 9200 | 8500 | 9000 | 9100 | 9700 | 9700 | 9400 |
| Mean | | 8200 | 8533 | 9367 | 8333 | 9067 | 8967 | 9733 | 9633 | 9033 |

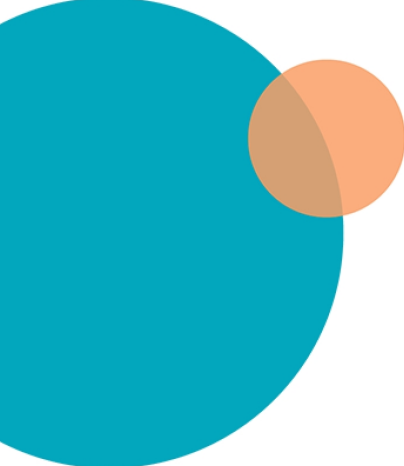
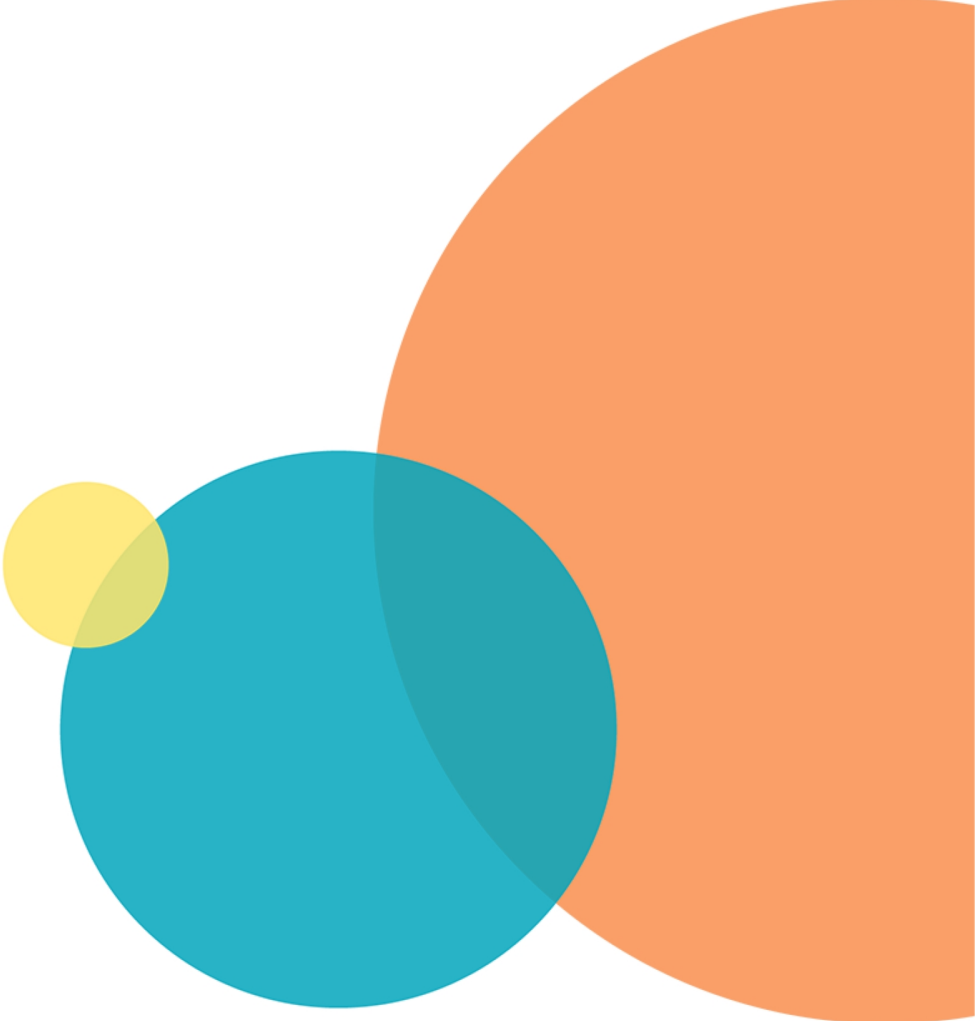


Note: In actual practice, data from all 96-wells are included in the calculations, but in this example, data from only nine wells are presented, to simplify the examples.

All samples taken from a particular well are divided by their well-specific mean baseline correction factor. For example, all samples from A1 are divided by 8200, A2 by 8533, and so on.

The results of applying the response over baseline correction factor are presented in the table below. Note the data is displayed as a fold increase of the response compared to the baseline with Sample values ranging in the single digits (0–6).

| Sample | Time | Well | | | | | | | | |
|--------|------|--------------|--------------|--------------|-----------|-----------|-----------|--------------|--------------|--------------|
| | | A1 – Ctrl | A2 – Ctrl | A3 – Ctrl | A4 Exp | A5 Exp | A6 Exp | A7 + Ctrl | A8 + Ctrl | A9 + Ctrl |
| 1 | 0 | 0.98 | 1.00 | 1.01 | 0.98 | 1.01 | 0.98 | 1.03 | 0.99 | 0.96 |
| 2 | 5 | 1.00 | 1.01 | 1.00 | 1.00 | 0.99 | 1.00 | 0.98 | 1.01 | 1.00 |
| 3 | 10 | 1.02 | 1.00 | 0.98 | 1.02 | 0.99 | 1.00 | 0.98 | 1.01 | 1.00 |
| 4 | 20 | 1.05 | 0.98 | 0.98 | 5.4 | 5.51 | 4.68 | 5.86 | 5.40 | 5.98 |
| 5 | 25 | 1.05 | 1.03 | 0.98 | 4.44 | 4.63 | 3.90 | 5.45 | 5.195 | 5.54 |
| 6 | 30 | 1.07 | 1.00 | 1.01 | 3.60 | 2.76 | 3.23 | 5.14 | 5.29 | 5.54 |



Appendix F: Using AquaMax Sterilant

F

AquaMax Sterilant can be used to clean and sterilize the fluid paths in the Cell Reservoir and Cell Suspension module.

Principle of Use

Accumulated residues and contaminating microorganisms have the potential for causing significant adverse effects on the operation of liquid handling equipment. Proteins in cell culture media, salts and minerals in buffers and other reagent components used in liquid handling equipment can build up on fluid path surfaces, clogging channels and probes resulting in incomplete or inaccurate dispensing and aspiration of liquids. In addition, microorganisms such as bacteria, fungi and molds can also accumulate in the fluid path creating similar fouling problems. Simply flushing the fluid path with water or ethanol/water solutions is often ineffective in the removal of residue.

AquaMax Sterilant was developed for the purpose of removing organic and inorganic material, and killing microorganisms, including endospores, in liquid handling instrument fluid paths. In addition, AquaMax simultaneously removes endotoxin from the fluid paths. Endotoxin is a substance produced by Gram negative bacteria that is known to cause a variety of significant cellular reactions *in vitro*, including cell activation and death. Thus minimizing the amount of endotoxin in fluid paths is necessary in order to avoid detachment and adverse effects on mammalian cells during the course of subsequent experiments. A two-hour treatment with AquaMax substantially removes endotoxin adhering to interior fluid path surfaces.

Contents

- 4X Concentrated Solution A: detergent in buffered aqueous solution (2 x 1 L bottles)
- Solution B: sodium hypochlorite in aqueous solution (2 x 14 mL bottles) [16 500 mL DW4 washes]

Materials Required but Not Provided

- Deionized water
- 250 mL graduated cylinder
- 1 L beaker, bottle or flask
- 2 mL pipettor

Storage

The unopened kit package can be stored at room temperature (between 10°C/50°F and 40°C/104°F).

After opening the kit, store 4X Concentrated Solution A and 1X Solution A (prepared by mixing 1 volume of 4X Concentrate Solution A with 3 volumes of deionized water) between 10°C/50°F and 40°C/104°F.

After opening the kit, store the tightly capped bottle of Solution B between 2°C/35.6°F and 8°C/46.4°F. Avoid freezing any of the solutions.

Reagent Preparation

500 mL Sterilant is sufficient to run the automated cleaning cycle on the AquaMax DW4 liquid handling system.

1X Solution A can be stored for 30 days in a capped bottle between 10°C/50°F and 40°C/104°F.

Preparation of 500 mL 1X Solution A

- Combine 125 mL of 4X Concentrate Solution A and 375 mL deionized water in a 1 L beaker, bottle or flask and mix thoroughly.

For other needed volumes:

- Add 1 part of 4X Concentrate Solution A and 3 parts deionized water.

Preparation of Complete Sterilant:

- Within 1 hour of use, add 0.2% (final v/v) Solution B to 1X Solution A and mix thoroughly.

For 500 mL of complete Sterilant:

- Use 1 mL of Solution B.

Reagent Use

Follow the recommended cleaning procedures in the user guide for the specific liquid handling instrument to be cleaned.

For stubborn cleaning situations, the Sterilant must sit static in the fluid path for 30 minutes half-way through the cleaning cycle before continuing to completion.

For endotoxin removal, the Sterilant must sit static in the fluid path for 2 hours half-way through the cleaning cycle before continuing to completion.

Rinse the fluid path with at least twice the volume of deionized water immediately after cleaning.

Warnings, Precautions and Limitations

- For research use only.
- When handling these reagents, gloves and eye protection must be worn.

- Do not ingest.
- Use in a well ventilated environment.
- For more safety information, see the Material Safety Data Sheet for Solutions A and B.
- As supplied, Solution A might not be sterile.

The following information is being provided in compliance with Worker and Community Right-To-Know Regulations:

Solution A, 4X Concentrate

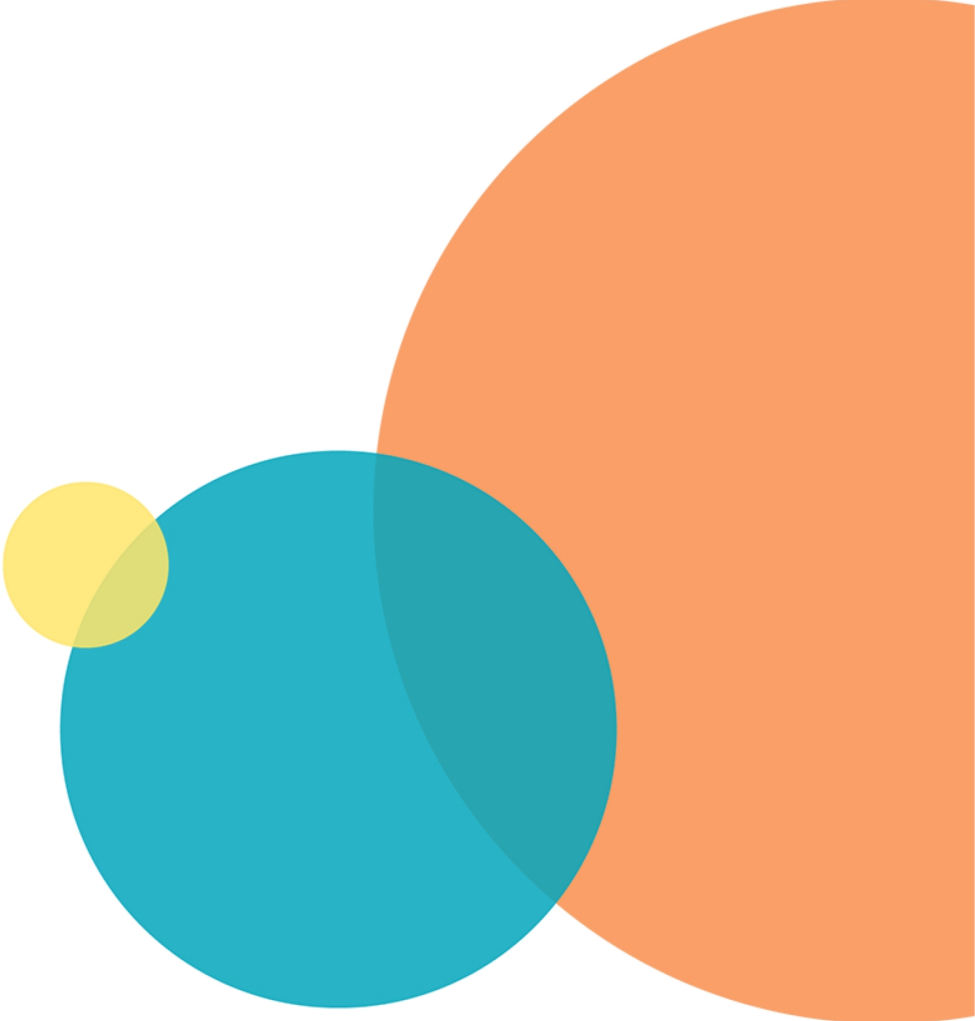
Chemical Composition

The chemical composition of this solution is a trade secret and proprietary property of Molecular Devices.

Solution B

Chemical Composition, CAS #

Sodium Hypochlorite, Solution, 7681-52-9



Appendix G: Robotic Integration



This document describes the interface between a FLIPR Penta System and a robotic instrument. The intention of the document is to provide a general overview of the interface between the instrument and robot and a detailed description of the communication protocol between the ScreenWorks Software, which run on the host computer of the FLIPR Penta System instrument, and the robotic control software.

Information contained in this document might change as new development occurs—we recommend checking with your Molecular Devices sales representative for updated information.

Conventions

- Bold text is used for commands
- Italic text is used for parameters
- Parameters in bracket [parameter1] are optional
- Parameters are separated by comma
- <CR> is the ACSII code for 'carriage return' which indicates end of the command or end of the response
- <TAB> represents a tab
- <SP> represents a space
- | represents OR

Interface Versioning

As of version 1.1 of the automation interface, the interface is versioned separately from the ScreenWorks Software and FLIPR Penta System firmware. The interface now provides a command so it can be queried for its version number. This command is described in detail later in this document.

Instrument Overview

Basic Function

FLIPR Penta System is a fluorescence microplate reader used primarily for kinetic live cell-based fluorescence and luminescence assays. A typical assay involves the following steps:

- Introduction of a read plate containing cells in solution or compounds into the instrument.
- Acquisition of a baseline read (or series of reads) to determine the background fluorescence or luminescence signal from the read plate.
- Introduction of a potentially activating (or potentially inactivating) compound into the well or cells in suspension.
- Monitoring the changes in the fluorescence or luminescence signal from the read plate for a period of time immediately following the introduction of the compound.
- Further additions of different compounds to the well are typical in an assay on a FLIPR Penta System. Between additions, the pipette tips used for fluid transfer may be replaced or washed.

Cell response in a typical assay typically begins within 2 seconds of fluid addition and is monitored for 1–3 minutes. Fluorescence or luminescence signals are monitored at wavelengths and frequencies selected by the user within the physical constraints of the instrument.

Hardware Introduction

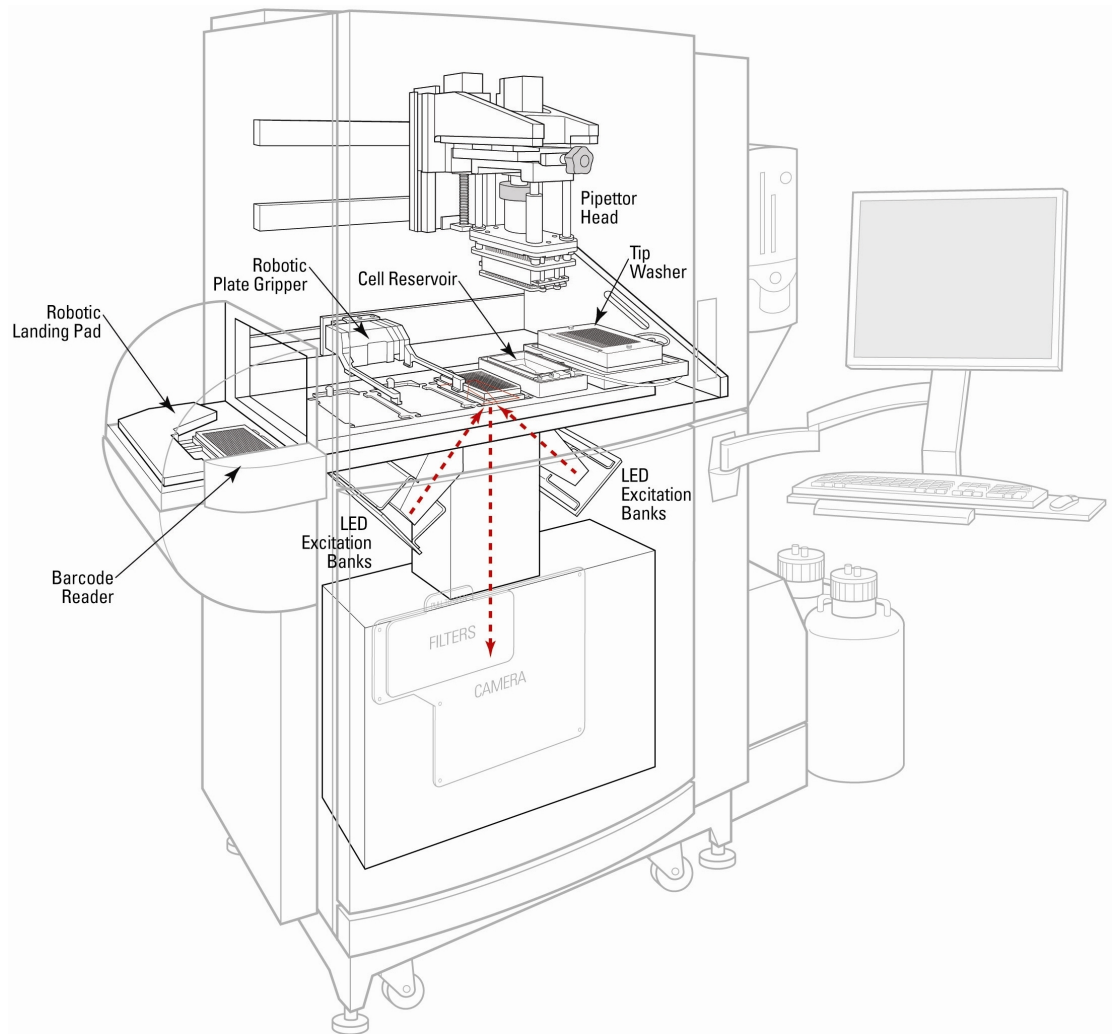
To accomplish the assays described, the instrument includes the following subsystems:

- Embedded control system — The instrument contains an embedded control system for use in controlling the instrument. This control system is accessed only by the ScreenWorks Software, which is supplied with the instrument.
- Plate reader — The optics subsystem includes an illumination source with excitation filters, emissions filters in a three-position filter changer, and a CCD camera. This system is used to read both fluorescence and luminescence assays.
- Pipettor — In order to accomplish kinetic assays, the instrument contains a pipettor which can add fluids to the read plate during a series of reads. The pipettor, equipped with a 96-well, 384-well, or 1536-well head, can be used for general purpose fluid transfer between any of the plates in the instrument. Each head type uses pipette tips, or a pin tools. For 96- and 384-well pipettor heads, this is a disposable plastic pipette tip. For 1536-well this is a metal tip block. Alternatively, 384-well and 1536-well pin tool heads have replaceable pin tools.
- Tip and pin washer — Each instrument includes a washer which can be used to wash tips or pins between fluid transfers. The washer consists of two parts—a washer control module and a wash reservoir. Installed on the right side of the instrument, the sensors, and control valves. These components connect to the wash reservoir located in position 5 of the system stage.

- FLIPR Cycler Plate Handler — An optional plate handler is available that is capable of moving plates from a location outside of the instrument to any of the first four stage positions and vice versa. The second component of the FLIPR Cycler Plate Handler is an automated door system that allows plates to be exchanged while an assay is in progress.

Layout

The basic layout of the instrument subsystems is demonstrated in the following illustration:



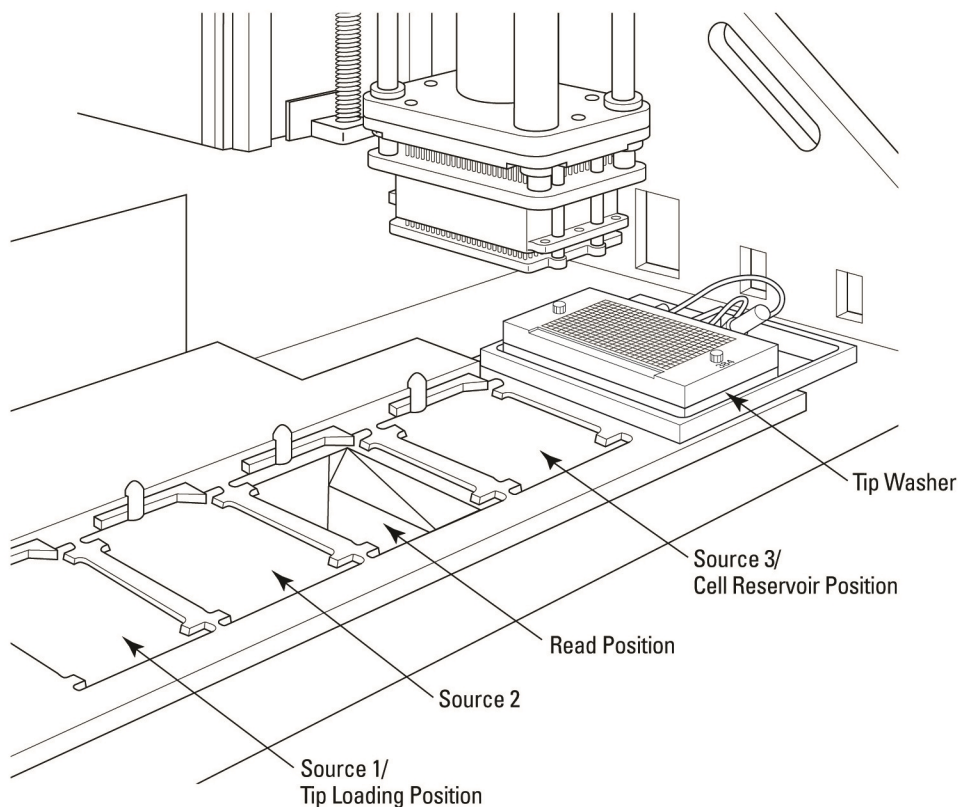
As the illustration indicates, the top instrument compartment contains the pipettor and plate platform (stage). Plates are placed in the read position of the 5-position platform in order to be read by the plate reader, which is located in the optics compartment below the platform.

Plate Layout

The stage has 4 positions where plates and/or tips may be placed. These may be:

- Position 1: Source Plate 1 or Tip-loading
- Position 2: Source Plate 2
- Position 3: Read Plate
- Position 4: Source Plate 3 or Cell Reservoir

A tip wash reservoir for tips or pins is available in position 5.



Users may place either tip racks or plates in the tip-loading position (position 1) and Cell Reservoir or plates in Cell Reservoir position (position 4). The wash reservoir is not disposable but is replaced by users when the head type is changed.

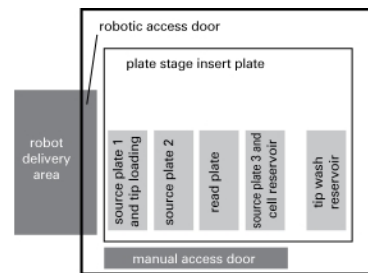
Plate-Handling System

Users may manually place plates or tip racks in the instrument or deliver them automatically using a robot or stacker.

In manual mode, users access the plate stage through a large, manually operated door. This door opens on a vertical hinge, mounted to the left side of the door. An observation panel is built into this door and may be removed by users to attach to the top instrument compartment. In this arrangement, users can observe assays in progress (data quality can be poor or simulated) in order to debug assay setup issues.

At the start of an assay, the door latches to prevent users from accidentally interrupting an assay in progress or injuring themselves. The door also latches for most other instrument actions not associated with an assay, for example, load tips, unload tips, and signal test. This latch remains engaged throughout the assay or operation. In manual mode, this latch automatically disengages at the end of the assay.

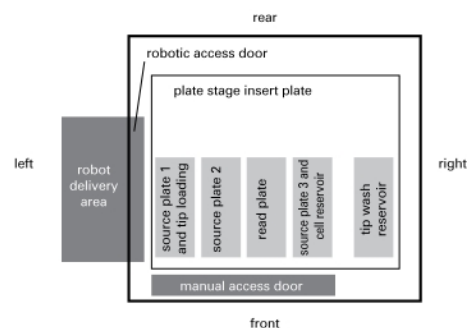
For robotic integration, plates are delivered to a single location outside of the instrument. The placement area for robotic delivery of plates to the instrument is to the left of the instrument. The locations for manual and robotic access to the instrument are shown below:



A further illustration of this arrangement shows the relation of the plate handler to the pipettor system.

Terminology

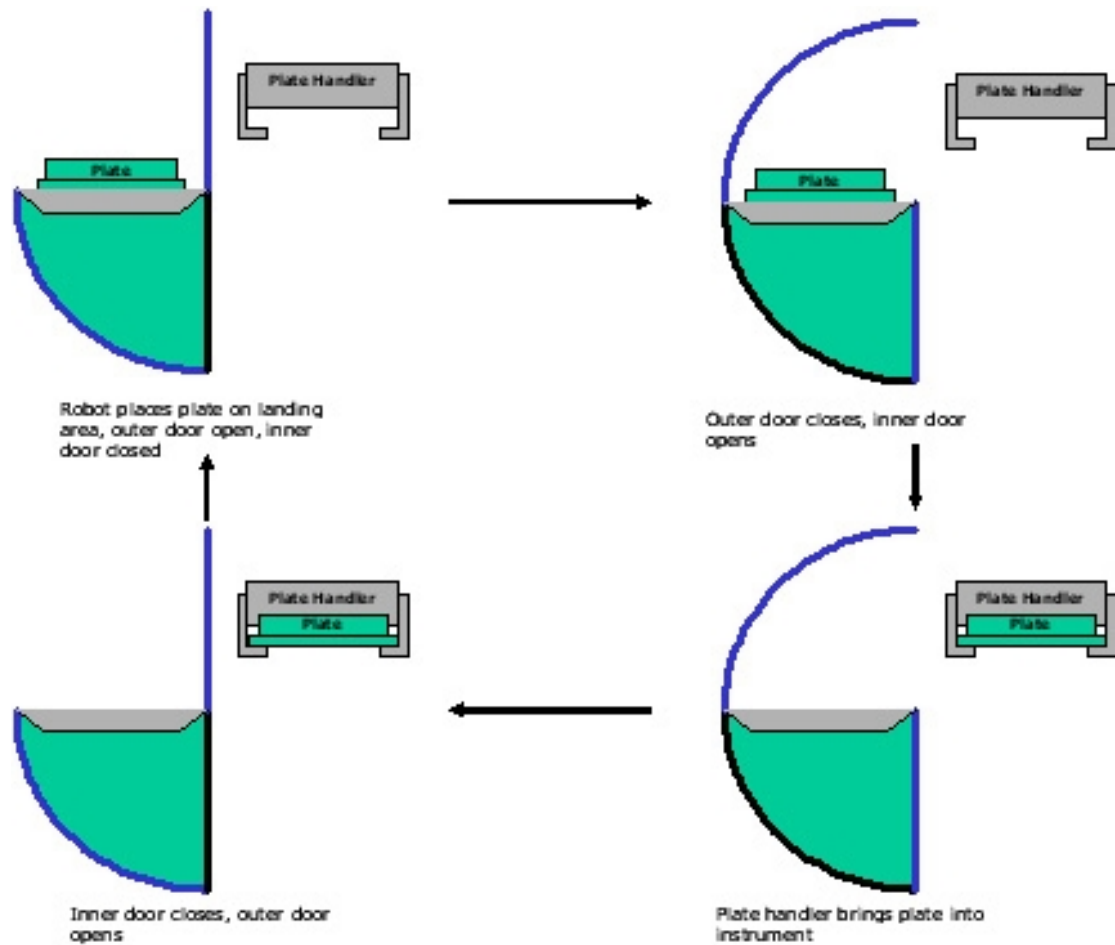
The manual access door side of the instrument is called the front of the instrument. From this, the following illustrates the named sides of the instrument:



Robotic Plate Loading

Plates are shuttled into the instrument by a plate handling subsystem called the FLIPR Cycler Plate Handler. To lessen dead time between assays and to handle more source plates in an assay than can be accommodated on the plate stage, the robotic access door(s) and FLIPR Cycler plate handler are able to exchange plates and tips while an assay is running.

The following diagram illustrates the steps involved in loading plates in robotic mode:



A reverse process is used when unloading plates in remote mode.

Optics Access Door

The plate reader optics are mounted to the plate stage directly below the read plate position. A door is provided to the optics compartment for users to change emission filters and LED modules. This door is also latched when an assay is in progress to prevent users from interrupting assays or injuring themselves.

Washer Placement

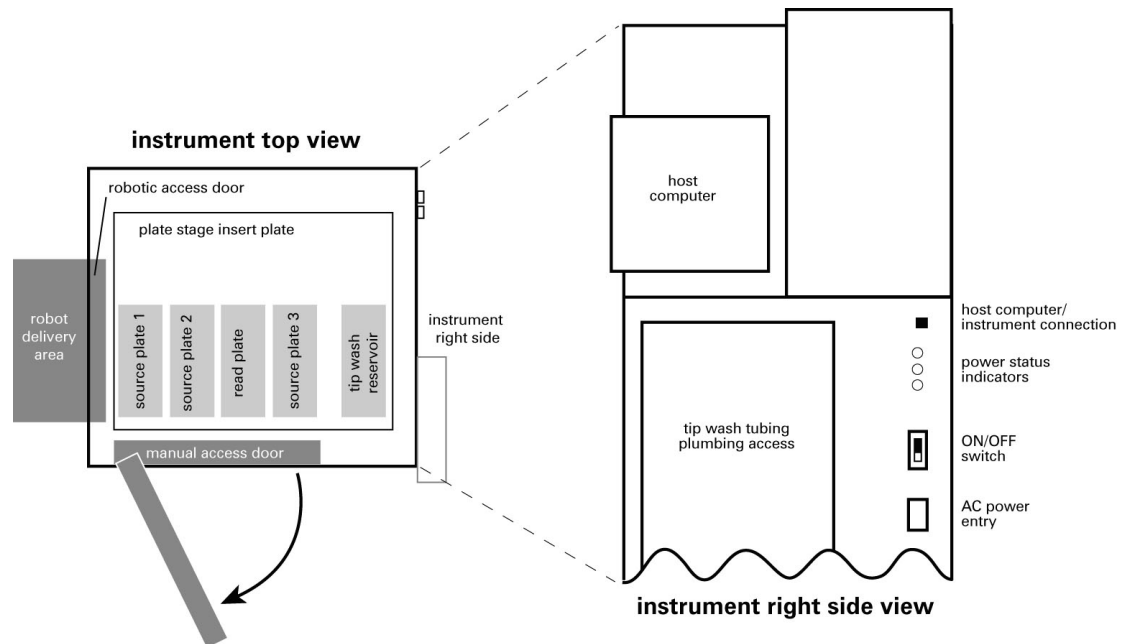
A tip/pin washer is included with the system and is placed on the lower right side of the instrument, towards the front. Extensive work has been done to isolate plumbing for the washer from other optical or power components. Access for tubing to the washer is required in any installation.

Cell Suspension Placement

The Cell Suspension module option is available for the FLIPR Penta System. When installed, the Cell Suspension module is mounted on the lower right side of the instrument and the Cell Reservoir is inside of the instrument. Access to the module is necessary during testing.

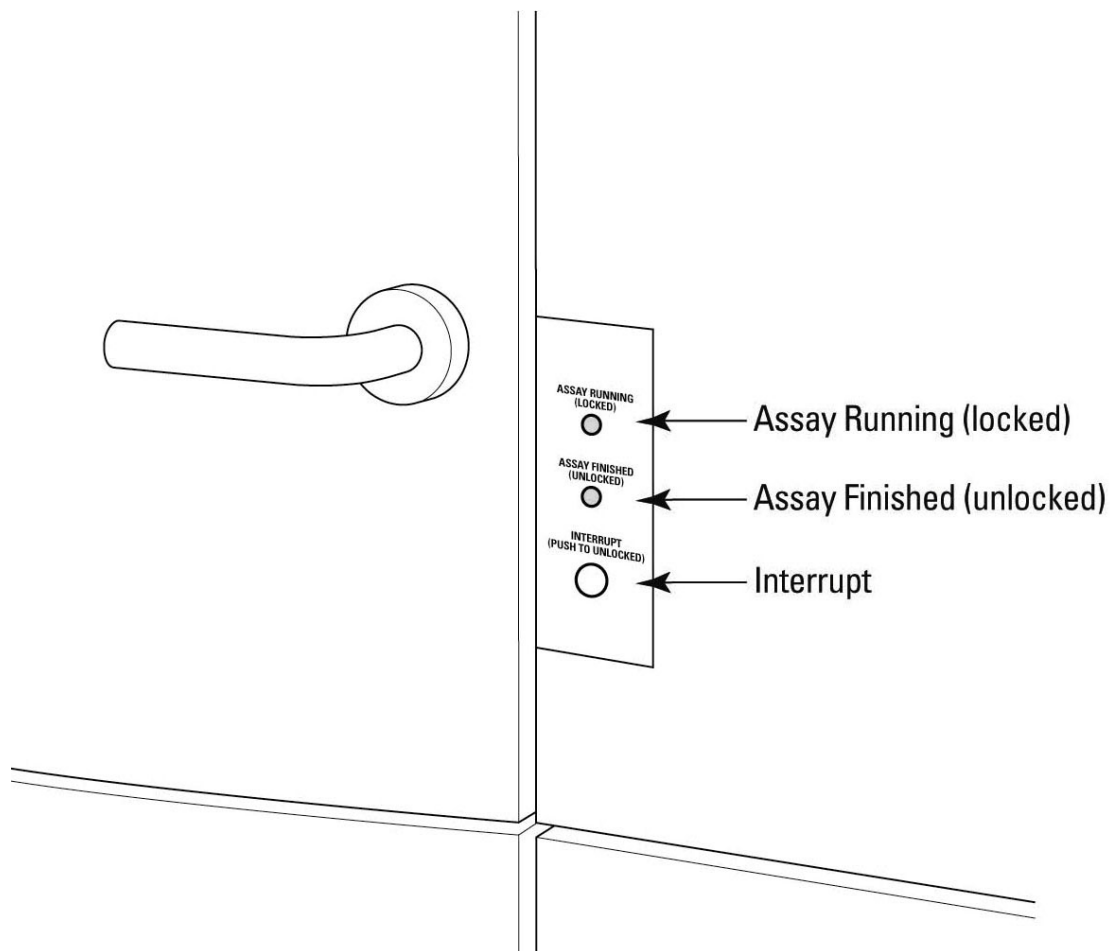
Other Instrument Access Areas

The following diagram illustrates other important access areas in the instrument.



On the right rear of the instrument, AC power enters. Directly above this entry is the main power switch. The power switch should be accessible in both robotic and manual modes. There are three indicator lamps near the power switch. Access to view these indicators can be useful for debugging instrument power problems. The main communication connection from the instrument to the host computer is located immediately above these power indicators.

The instrument status panel, located next to the upper door handle, indicates whether or not the instrument is running and safe to open. It includes an emergency Interrupt button to stop any processes.



The panel has two lights as well as the Interrupt button. From the top of the panel these are:

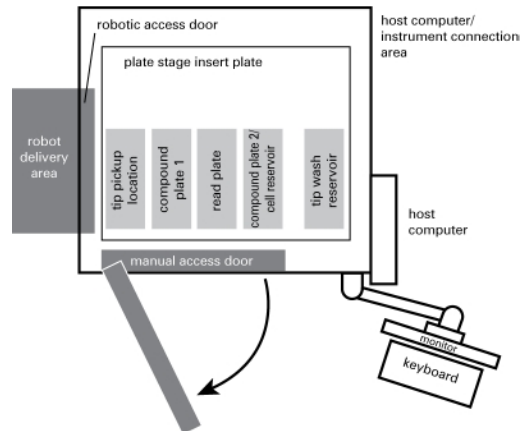
- Assay Running (Locked)—Yellow light
The FLIPR Penta System is performing a task. The upper and lower doors are locked and cannot be opened until the task finishes or is halted using the Interrupt button.
- Assay Finished (Unlocked)—Green light
No tasks are being run and it is safe to open the upper and lower instrument doors.

The Interrupt button is an override button to halt all tasks, so you can access the instrument. If pressed the yellow light flashes until the system has reached a safe state to open the doors, then the green light comes on.

CAUTION! The Interrupt button immediately ends the experiment and should only be used in emergencies. Interrupting the instrument while an assay in progress almost always results in the loss of the plate being run. If this button is pressed, the yellow ‘busy’ light flashes until the instrument is safely unlatched. The system may need to be reinitialized by selecting Reset from the Instrument menu prior to resuming normal instrument function.

Monitor and Keyboard Placement

The host computer provided with FLIPR Penta System is mounted to the instrument's right side. This computer provides the only direct user interface and path to control the instrument. Monitor and keyboard for this computer are mounted to an adjustable arm on the instrument's front right. The following image illustrates this installation.



The host computer, keyboard, and monitor may be moved to a different location at the user's discretion. Wherever the host computer is placed, it must be directly connected to the instrument. It is not possible to communicate between the host computer running the ScreenWorks Software and the instrument across a network. A short communication cable is provided between the computer and instrument. In order to move the computer, the user may need to provide an alternate cable. This cable must be a shielded crossover Category 5 cable.

Required Access Areas

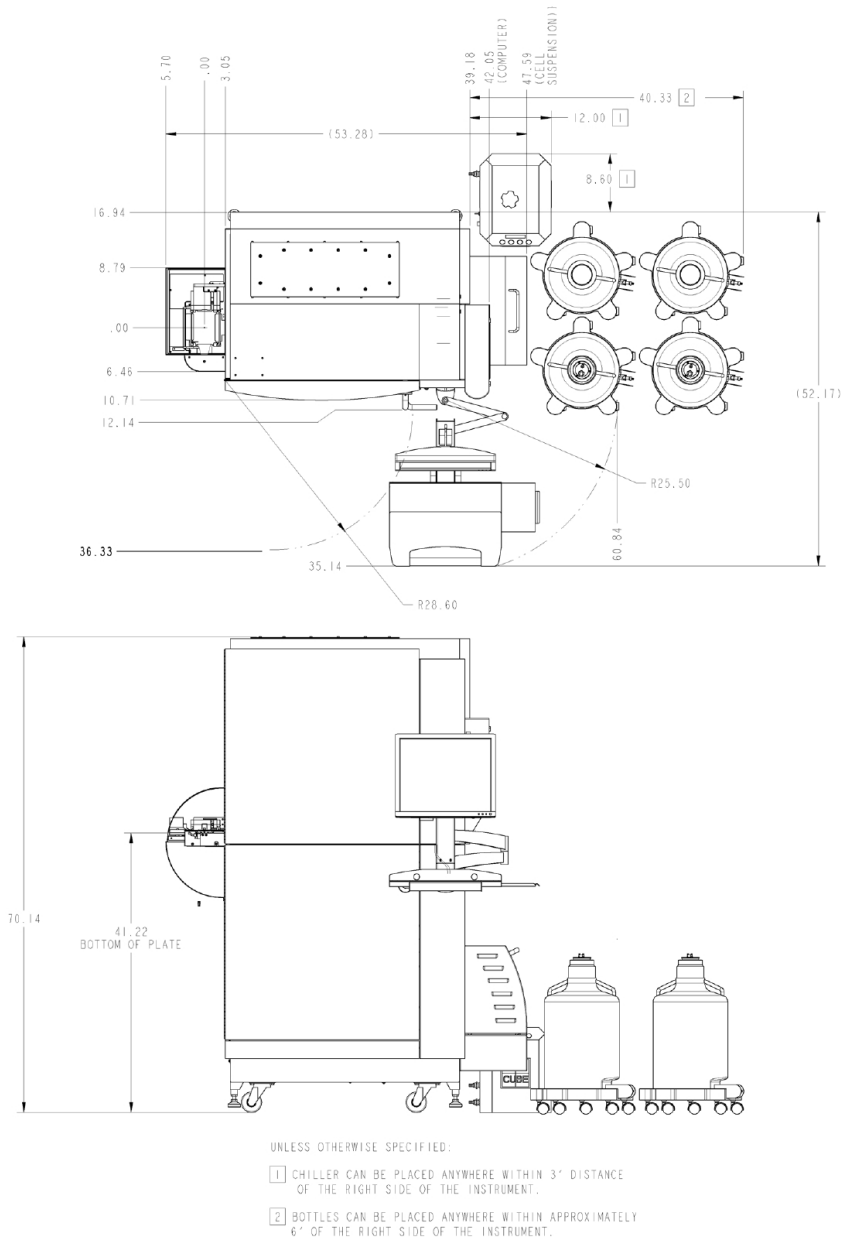
The all inclusive system dimensions with the FLIPR Cycler Plate Handler and Cell Suspension modules are approximately 53 inches (1346 mm) wide by 29 inches (737 mm) deep by 70 inches (1787 mm) tall. The FLIPR Penta System is designed with rolling castors so it can be readily moved to make necessary adjustments and perform maintenance. Leveling feet are also installed on the lower instrument chassis. These feet are typically used for stabilizing the instrument when integrated with a robot, but can also be used to establish a uniform instrument deck level in situations where the lab floor is not flat. When running an experiment, please make sure the instrument's feet are lowered and leveled.

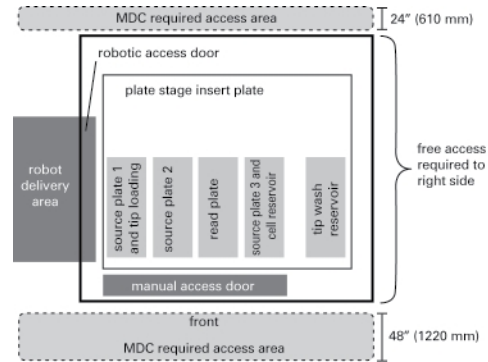
The computer and monitor are mounted to the right front side of the instrument with the included clamp, requiring a minimum lab space of 85 inches (2159 mm) wide by 61 inches (1549 mm) deep for maneuverability. A minimum 28 inch (711 mm) square footprint for tip wash bottles and chiller is required to the right side of the instrument. This space is also recommended to allow access to the main power-switch, washer tubing and communication connections.

The cabinet should have a user access space of 48 inches (1219 mm) in front for users to exchange tips, plates and reservoirs. An additional 24 inches (610 mm) behind and 10 inches (254 mm) to the left should be left available for servicing the instrument.

⚠ WARNING! The FLIPR Penta System can weigh as much as 860 lbs (390 kg). Ensure adequate personnel are present when installing or moving the system. Follow all necessary safety precautions and use proper moving techniques.

The following drawings illustrate these requirements:





FLIPR Penta System Control Architecture

General Description

The FLIPR Penta System consists of a host computer and a stand-alone instrument. ScreenWorks Software is installed on the host computer supplied with the instrument. This software is the primary user interface. ScreenWorks Software is used to setup protocols, run assays, analyze and export data. ScreenWorks Software communicates to the standalone instrument with a dedicated connection using a communication protocol proprietary to Molecular Devices. The status of the instrument and assay is monitored by ScreenWorks Software, which then displays data as it is collected.

A stand-alone instrument has all the hardware and firmware components required to perform an assay and report the results. It does not have the ability to display the results of experiments directly to a user. The user may not directly communicate with the stand-alone instrument in order to obtain data or to configure and run assays.

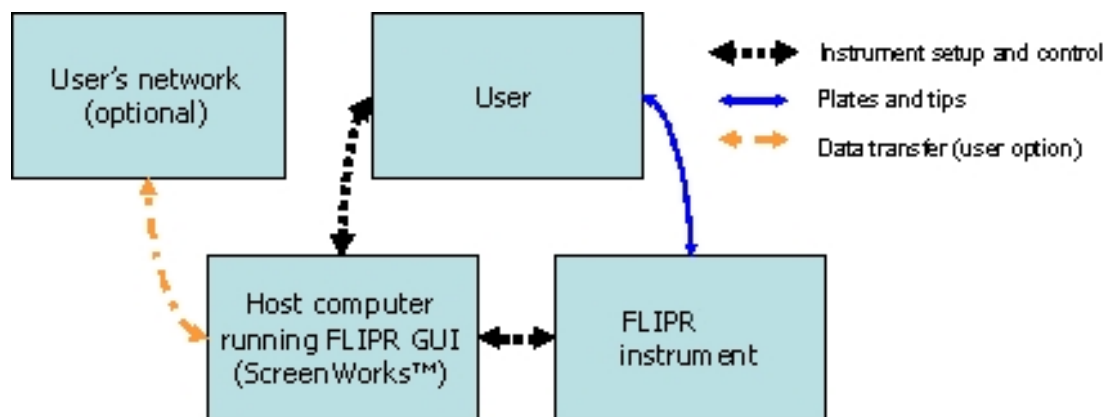
The host computer supplied with the FLIPR Penta System is configured with two Network Interface Cards (NICs). The first NIC is used to communicate with the instrument using a fixed IP address - this connection cannot be used for any other purpose. The second NIC is provided to connect the host to the Local Area Network (LAN) or to other devices (for example, router, hub, scheduling workstation).

ScreenWorks Software and the instrument have two operating modes. The first of these is Manual mode. The second mode is Remote mode. ScreenWorks Software must be placed in **Remote Mode** in order to be controlled externally by third-party software via the automation interface.

Manual Mode

By default, the instrument starts in Manual mode.

In Manual mode, configure assays in ScreenWorks Software and manually load plates and tips into the instrument. The following simple diagram illustrates this interaction:



In Manual mode, a second network connection on the host computer is not monitored. It is provided primarily to connect to your network to transfer data from the FLIPR Penta System host computer for storage or analysis elsewhere.

Remote Mode

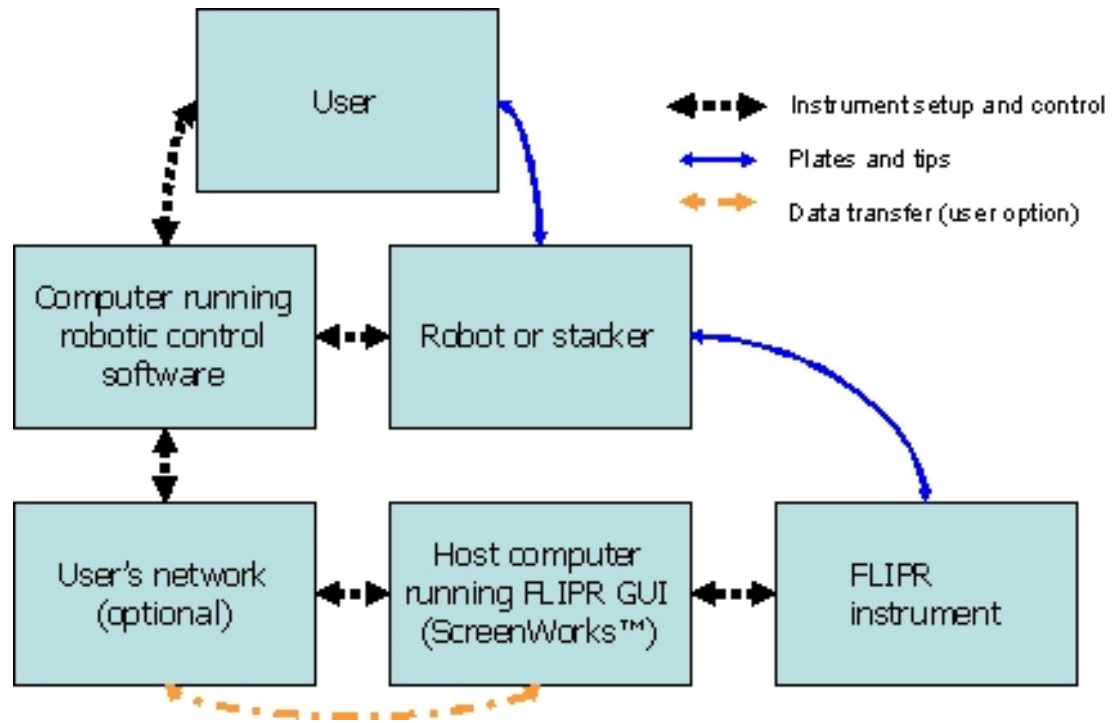
By transitioning into Remote Mode, ScreenWorks Software can be configured to communicate with external devices such as robots and stackers supplied by third parties. These devices can be used to supply plates and tips to the FLIPR Penta instrument.

When in remote mode, third-party software is given full control over the internal plate handler, FLIPR Cycler plate handler. The FLIPR Penta System software does not initiate any plate handling events. The general philosophy here is to put the third party software in charge and provide status information to allow it to make decisions on when the FLIPR Penta System should load or unload plates and start assays. Third-party software monitors the assay through the ScreenWorks Software automation interface and provides additional plates and tips to the instrument as needed. External devices might not communicate directly with the FLIPR Penta System instrument but only with the automation interface.

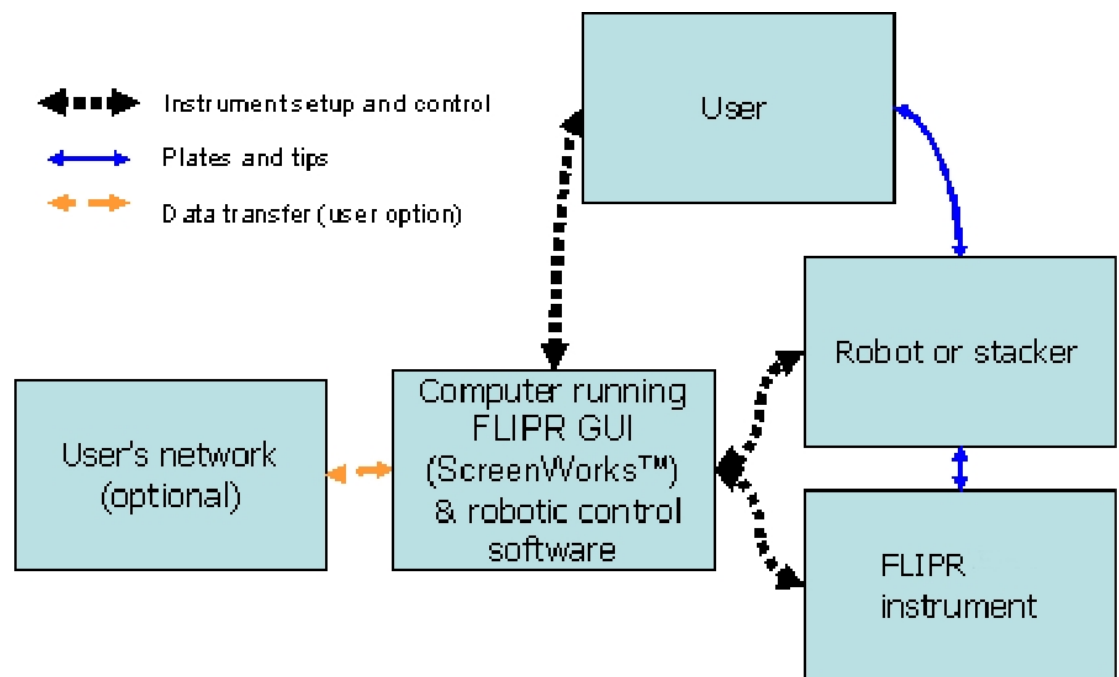
Interact with the third-party software to select a protocol or preload a protocol in ScreenWorks Software before entering remote mode. The automation interface does not provide the functionality to create assay protocols. Interact with ScreenWorks Software while assays are running remotely to QA data and monitor progress.

The control system to communicate between ScreenWorks Software and robotic controller can be configured in several ways. The following illustrations are provided to illustrate the range of options available. These examples are not exhaustive.

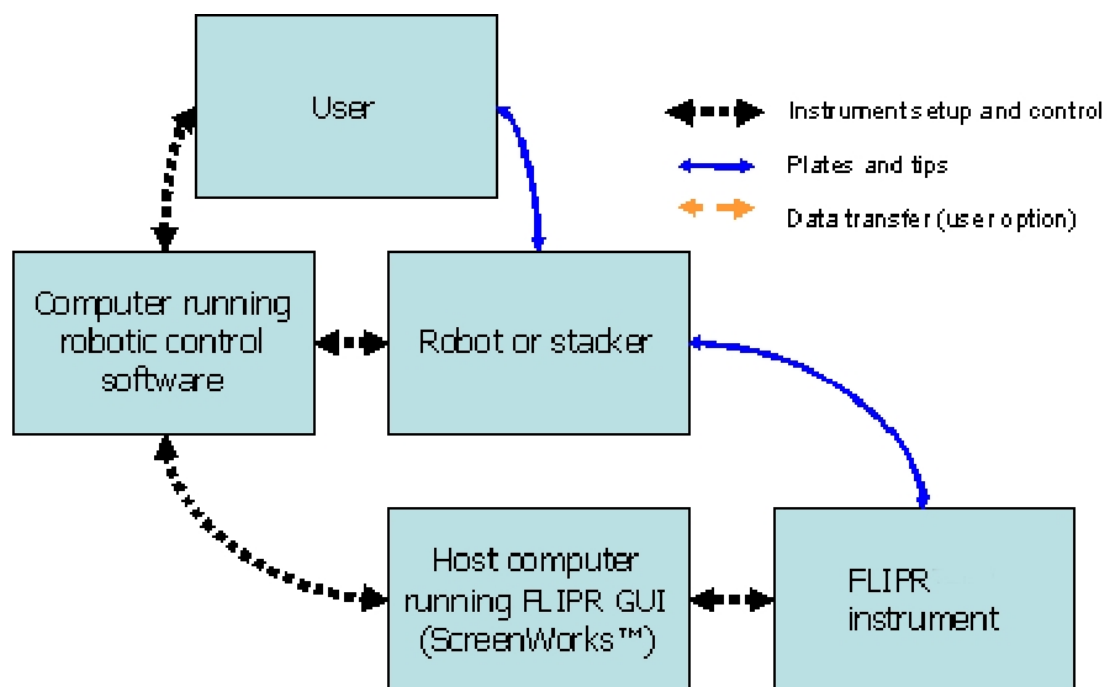
Third-Party Software via Network



Third-Party Software on the FLIPR Penta System Host Computer



Direct Connection From a Third-Party Software Host to the FLIPR Penta System host



Communication Protocol and Address

Communication between third-party software and ScreenWorks Software uses the TCP/IP protocol. The third-party software can be resident on another computer outside of the FLIPR Penta System or on the instrument host computer.

If the third-party software is remote to the host computer, the ScreenWorks Software receives communication through the IP address of the host computer on which it is running.

If the third-party software is run on the instrument host computer, the ScreenWorks Software receives communication through IP address 127.0.0.1 (localhost).

In all cases, ScreenWorks Software sends and receives commands through TCP/IP port 7.

General Remote Mode Use

This section provides a brief overview of running a FLIPR Penta System in remote mode. Later sections describe some detailed issues in more detail. This section provides some context for the introduction of the interface command set in the following section.

Instrument Startup

When the instrument is powered on, it performs a series of initialization steps which can take several minutes. These steps include:

- Configuring the camera
- Initializing the pipettor
- Initializing the plate handler and automatic door (if included)
- Determining and checking the installed LEDs
- Determining the installed emission filters
- Initializing the tip washer
- Communicating with the chiller

By default, the instrument is in manual mode upon startup. In this mode, the main manual access door is unlocked, the outer automated door is closed, and the FLIPR Cycler plate handler is moved outside the instrument, above the automation landing pad. As long as the instrument is in manual mode, the outer automated door and plate handler do not move from these positions, except during an instrument reset.

Software Startup

To use the instrument, the host computer must be started and ScreenWorks Software must be run. ScreenWorks Software establishes communication with the instrument and determine the instrument status and configuration. The instrument status and configuration can be determined by looking at the **Instrument Status** toolbar. This toolbar has two tabs. The **Instrument Status** tab shows the current instrument status including the status of assays in progress. The **Instrument Config** tab shows which LED modules, emission filters, and heads are installed.

Protocol Creation

The user must create assay protocols in manual mode. They do not need to load these protocols prior to placing the instrument into remote mode, but any protocol that is used in remote mode must be created and saved with a valid protocol file name prior to entering remote mode.

Users should test all protocols in manual mode prior to attempting to run them in remote mode.

Instrument Setup

Prior to placing the instrument into remote mode, the user may want to load tips or plates into the instrument. These actions can be done in remote mode but in certain conditions it is advisable to do this through the main manual door prior to entering remote mode.

If tips are used for an extended period of time and the tips/source 1 position is used as a source position, it is best to load the tips manually before placing the instrument into remote mode.

If boats or reservoirs are used as source plates, these should be loaded by hand in manual mode. The FLIPR Cyclor plate handler must not be used to load reservoirs because the plate handler moves rapidly and large wells spill liquid within the instrument.

Instrument Mode Setting

Prior to running an assay in remote mode, users must place the instrument into remote mode by selecting **Instrument > Set Remote Mode** from the main menu.

After the user requests transition to remote mode, the instrument locks the main manual doors, moves the plate handler inside of the instrument, closes the inner automated door, and opens the outer automated door. This is the default position for remote mode. At the end of any plate handling operation, the instrument returns to this position. The outer manual doors remain locked as long as the instrument is in remote mode.

The instrument assumes that any plates in the instrument when it transitions to remote mode are acceptable for use until a protocol has been run, which indicates that the plate is no longer needed.

Loading Protocols

In remote mode, protocols are loaded using the **openprotocol** command. This command loads the requested protocol and closes any other files open at the time that the command is sent. Protocols open regardless of whether they are valid. This command is described in more detail later in this document.

Running Protocols

After the instrument has been placed in remote mode and a protocol is opened, the protocol may be run by sending the **runprotocol** command. When a protocol begins to run in remote mode, any data files currently open close. This command is described in more detail later in this document.

Instrument Status Updates

The instrument can be queried at any time for instrument status, even when in manual mode. The status command is described in detail later in this document.

Plate Status Tracking

The most important information in the status command is the status of the plates in the instrument. From the status response, four possible information items may be determined about the plates in each of the four plate positions.

The first item that can be determined is the plate assigned to the plate position for the protocol to be run. This is the plate name assigned in ScreenWorks Software for a particular plate type and assigned to that position for the current protocol. This information could be used by the robot to confirm that the correct plate type is delivered to the required position. ScreenWorks Software and the FLIPR Penta System instrument have no method to determine whether the plate delivered is of the required type and always assumes that the plate type delivered is correct.

The status response also indicates whether a plate is present in each possible location. This indication is a direct readout of the status of the plate presence sensors in the instrument. This readout does not depend on whether a plate is required in the position or if the plate present is the correct plate.

The **REMOVEPLATES** line in the status signals that the plates in the indicated positions have completed. Plates complete in different ways. If the plates are source plates, ScreenWorks Software indicates that they are finished when an aspiration from the plate is completed and that aspiration is checked to show that the plate is finished after the aspiration. For tip racks, they are marked as complete as soon as tips have been loaded or unloaded to/from them. Read plates are marked as complete at the end of the last read in the assay in which they are used.

The **NEEDPLATES** line indicates plates that are needed for the assay in progress to continue. The instrument indicates that it needs a plate if it gets to an operation that requires a plate in a particular location and that location is either empty or contains a plate that was previously marked as finished.

Bar Code Tracking

Bar codes can be tracked in an assay in two ways.

First, the bar code can be supplied to the software by the robot controller when the robot controller instructs the ScreenWorks Software to load a plate. The software records the bar code values of the plates delivered in the appropriate data file in which they are used.

Second, the instrument attempts to read a bar code for all plates it loads using the plate handler. If this value is not overridden by one supplied by the robot controller, this bar code is recorded in the data file when the plate is used.

Regardless of the method used to deliver bar code information to the instrument, all bar code information might not be present when an assay is started, since all required plates might not be present in the instrument. Since the software can be configured to name the data file based on the bar codes of the plates used, it is possible that the data file name changes as an assay is completed.

Setting Up Protocols for Remote Control Use

Protocols must be created in manual mode before entering remote mode.

Protocols should also be tested in manual mode to ensure that they work as expected.

Setting up protocols is described in the user guide and you should be familiar with this process before reading further in this section. See [Setting Up an Assay Protocol on page 149](#). This section describes particular items you should be aware of when setting up protocols to be run in remote mode.

Settings Screen

Read Mode Settings

In the settings screen, configure the read modes. The settings for the read mode are dependent on the assay dye, plate type, cell type, and condition. These settings are determined during assay development and no changes are required when switching into remote mode.



Note: It is not possible to automatically vary these settings based on the particular cell plate loaded.

Assign Plate Settings

Assign plates to the required positions for the assay in the **Assign Plate to Position** group box. The assignment of plates can be critical to the smooth functioning of the assay.

Only one type of plate can be assigned to a position throughout an assay. For example, it is not possible to assign a normal plate to a position for the first part of an assay and then assign a deep-well plate to that position for the rest of the assay.

The read plate is always assigned to the read position. There are no special suggestions in remote mode for this assignment.

The assignment of source plates and tips should be carefully considered. The pipettor and the plate handler are both used in remote assays and must share the space above the plate locations. The instrument monitors the locations of the pipettor and the plate handler and moves them if necessary to prevent collisions. The pipettor and plate handler are never allowed to be above the same plate position at the same time. Plates should be arranged to minimize potential interactions between the plate handler and the pipettor.

The following guidelines can be helpful in assigning plate positions for automation:

Open reservoirs with very large wells cannot be handled by the plate handler. These plates must be placed in the instrument manually prior to entering manual mode. Replacing these plates requires the user to halt remote control and place the instrument in manual mode to manually replace or refill the reservoir. Deep well plates are a good alternative to using these plates.

If tips are loaded in the experiment, they must be assigned to the Source Plate 1 location. When tips are assigned to this position, the position cannot be used for source plates. Unless the tips are going to be changed for every plate, it is a good idea to load the tips prior to placing the instrument into remote mode. Alternatively, the user could create a tip configuration protocol which load and wash the tips. This assay can be run and then the empty tip rack can be removed by the robot to allow the source 1 position to be used for other plates. As of v1.1 of the automation interface, tip loading and tip washing is available.

Plates that are switched frequently should be placed closer to the left side of the instrument (near the automated landing pad) than plates which are not switched frequently.

Deep well plates (or reservoirs) which can be used for a series of assays before being replaced can be placed in the source 3 location, as long as it is not being used for Cell Reservoir. This minimizes the interaction between the pipettor and the plate handler when switching plates during the majority of assays.

It is only possible to RESET the instrument if a tip rack is present in the instrument. A RESET is required if the instrument enters a fault condition. If automated recovery from fault conditions is desired, the tips/source 1 position should be used for tips only and an empty tip rack should be left in this position at all times.

Data File Name Settings

Configure the desired data file name in the **Data File Name** group box.

If the **Include Bar Code** box is selected, the bar codes for all of the plates which have been checked as requiring them are included in the data file name. Bar codes are separated by an underscore (_). If a bar code cannot be read, or one has not been provided during the **loadplate** command, the software records **Bad_bar_code**.

The settings for the data file name are ignored in remote mode when the data file name is included with the **runexperiment** command line. Care should be taken when using this command to ensure that data file names are always unique to prevent the overwriting of data files.

If the ScreenWorks Software is allowed to create the data file name, it always ends in **_###.fmd**, where **###** is the plate number with this base name. Using this option ensures that no data is lost due to file overwriting.

Folder Settings

This area is used to assign the directories used in saving the data and export files. There is no difference between settings for these directories in remote control mode and in manual mode.

Running in remote control can produce large amounts of data. ScreenWorks Software does not allow new assays to begin if less than 50MB of hard disk space is available. Monitor hard disk capacity and maintain adequate space to save data.

Saving data across a network connection is not recommended in any case. If a protocol is created that saves data across a network connection, you might be unable to run the protocol if the network connection becomes unavailable. If a protocol that saves across the network is running when a network connection is interrupted, data can be lost and the software can encounter problems.

Temperature Control Settings

This area allows the user to set the temperature control set point for the protocol. The instrument must already be at this set point prior to running the assay.

If a protocol requiring temperature control is loaded and the **runexperiment** command is sent, the software confirms that the instrument has achieved this temperature. If the instrument has not reached the temperature, the software responds with an error. If the temperature has been reached, the experiment runs.

Auto Print Options Settings

In remote mode, problems with the printer can disrupt the continued operation of the system. No automatic printing is recommended in remote mode.

Analysis Screen

There are no special notes regarding the Analysis screen in remote control. Setup this screen as you would for any manual assay.

It is very important to take care to set up groups, display, and reduction settings prior to running a large number of plates. These settings should be set in the protocol so that they are automatically replicated in each data file as it is created. It is very tedious to go back to each data file and change these settings manually.

Transfer Fluid Processes

In general, settings for transferring fluids in remote control are the same as in manual mode, with three exceptions. These are in aspiration height over multiple plates, and the **Pause In Well** edit box.

Aspiration Height over Multiple Plates

When running in remote mode, a user might want to use a common addition source plate over a series of plates. The protocol(s) for this can be done in a couple of ways. The simplest way to do this is to create a single protocol which aspirates from the lowest allowed point in the source well. When doing this, you have to determine the maximum amount of liquid in the well so that the liquid displaced during insertion of the tip into the well is not pushed out of the well. This limits the amount of liquid that can be used in the source well. Alternatively, you can choose to create a series of protocols at successively lower aspiration start levels.

To provide an example of this situation, consider a user using a 384-well plate with a 250 μL well. This source plate is used to add a common agonist to all wells. In the assay, 20 μL of this agonist is added. Leave the plate in the instrument for a series of plates before replacing it. The FLIPR Penta System allows protocols that begin aspiration as low as 8 μL for this plate. Inserting a 384-well tip all the way to the bottom of this plate displaces 50 μL of liquid.

In the simple method described above, do not start with more than 190-200 μL of liquid in the plate because any more can spill liquid when the tip is fully inserted. The single protocol begins aspiration at 8 μL height. Run this assay 9 times before replacing the plate, leaving 10-20 μL of liquid in the wells after the source plate is removed.

In the more complex method, fill the wells to 230–240 μL . Then, set up an assay that would start aspiration at \sim 240 μL in height, a second assay to start at \sim 220 μL , and a third assay to start at \sim 200 μL . Continue to create separate assays at different heights or use the third assay for the remaining plates. In this case, 11 plates can be run before the plates need to be changed.

The simple method is simpler to setup and run. The more complex method allows source plates to be switched less often. It also decreases the proportion of wasted common reagents. Finally, it can decrease carryover of compounds sticking to the outside of the tips, especially if the protocols are created to retrieve liquid from progressively lower heights all the way through the well height.

Pause in Well Edit Box

The **Pause in Well** edit box on the **Edit Dispense** screen allows you to require that the pipettor remain in the well after dispensing. This pipettor operation leaves the tips in the well for the time specified. When the tips are in the well, it might not be possible to exchange plates from any position, and certainly not from the read or source 3 positions. For this reason, use of this option is not recommended in remote mode.

Wash Tips Processes

No special requirements are needed for the **Wash Tips** process.



Note: The pipettor and the plate handler do not share the tip reservoir position. For this reason, plate handling can proceed unhindered during wash operations.

Mix Fluid Processes

The **Mix Fluids** process does not require any special changes during protocol creation.



Note: The same warning that applies to **Pause in Well** for dispensing also applies here.

Also, mixing in any well can interfere with plate handling, so use of this function might require some experimentation.

Read Processes

No special requirements are needed for **Read** processes.

Finish with Source Processes

This notifies the plate handler to remove the plates in remote mode.

Wash Cell Reservoir Processes

No special requirements are need for the **Wash Cell Reservoir** process.

Command Set

Command Syntax

All commands and parameters must be sent in lower case and end with **<CR>**.

All responses are given in lower case and end with **<CR>**.

Response from ScreenWorks Software has one of the following forms:

```
ok<CR>
errorcode, error string<CR>
status<CR>
```

- **ok<CR>** indicates the command was started successfully or command is executed successfully.
- **errorcode, error string<CR>** indicates command failed because of either an indicated condition or an instrument error. Error codes are followed by a comma, a space, and then a string describing the error of up to 300 characters. Error strings are entirely lower case.
- **status<CR>** is the formatted status text which was requested.

Version <CR>

Command Description

This command returns the version string for the automation interface command set.

This command is only supported for versions 1.1 and newer. Older versions return the error *c69, badly formed or unrecognized command* in response to this command.

Command Parameters

There are no parameters associated with this command.

Example Command Line

```
version<CR>
```

Example Response

```
1.1<CR> or c69, badly formed or unrecognized command.<CR>
```

Suggested Steps in Command Usage

This should be the first call made to the instrument after making a TCP/IP connection to ScreenWorks Software. To ensure problem free operation of your application for FLIPR Penta System customers, the robotic integrator must consider limiting use of the application to interface versions they have tested.

Status <CR>

Command Description

The **status** command returns the status of the instrument. The status response is an ASCII text string which includes:

- Information about the instrument type and firmware version.
- Whether the instrument is functioning normally or an error occurred. If an error has occurred, this field contains an error code and an error string of less than 300 characters. The error code and error string is separated by a comma and space.
- The instrument status field reports a summary of status information including whether the instrument is busy, faulted, error, resetting, aborted, or idle. It also reports whether the instrument is remote or manual mode. In manual mode, only status information is available through this interface. All other commands do not function. Finally, this line reports whether the software is online (connected to the instrument) or not.
- Whether an experiment is running or not—do not open or start another experiment until the first is finished.
- Name of the plates assigned to each position.
- Whether plates are present or not on plate locations.
- Whether plates need to be loaded or removed.
- Whether the plate handler is busy or idle.
- Whether the landing pad has a plate or not.
- Camera temperature. Status returns “not_ok” while camera CCD temperature is cycling and/or out of range as defined by factory firmware settings, “off” if camera temperature control is disabled, or “ok” otherwise. The factory setpoint depends on the installed camera:
 - EMCCD camera: -70°C (-94°F) $\pm 2^{\circ}\text{C}$
 - HS EMCCD camera: -70°C (-94°F) $\pm 2^{\circ}\text{C}$
 - ICCD camera: -20°C (-4°F) $\pm 5^{\circ}\text{C}$
 Also returns actual CCD temperature and setpoint.
- Chamber temperature. Status reports “not_ok” while chamber is not at setpoint temperature (typical factory tolerance of $\pm 0.5^{\circ}\text{C}/\pm 41^{\circ}\text{F}$, “fault” if sensor does not return a reading, “off” if chamber temperature control has been disabled, or “ok” otherwise.
- Waste/fill bottle status.
- Tips on | off.
- Outer auto door, open/closed—if it is closing it is reported as closed.

Example Command Line

```
status<CR>
```

Response Format

```
instrument <SP>instrument name<TAB>
version<SP>mm/dd/yyyy<TAB>
function<SP>ok|errorcode|fatalcode, [errorstring]<TAB>
inst_status<SP>busy|fault|error|resetting|abort|idle, remote|manual, o
ffline|online<TAB>
exp_running<SP>yes|no<TAB>
readplate<SP>platename|none, have|empty<TAB>
sourceplate1<SP>platename|none, have|empty<TAB>
sourceplate2<SP>platename|none, have|empty<TAB>
sourceplate3<SP>platename|none, have|empty<TAB>
loadplates<SP>re[, s1, s2, s3, ti]|none<TAB>
removeplates<SP>re[, s1, s2, s3, ti]|none<TAB>
platehandler<SP>idle|busy_load|busy_remove<TAB>
plate_on<SP>yes|no<TAB>
camera_temp<SP>ok|not ok|off, temp, setpoint<TAB>
chamber_temp<SP>ok|not_ok|fault|off<TAB>
waste_bottlea<SP>ok|full<TAB>
waste_bottleb<SP>ok|full<TAB>
fill_bottlea<SP>ok|empty<TAB>
fill_bottleb<SP>ok|empty<TAB>
tips<SP>on|off<TAB>
outer_auto_door<SP>open|closed<CR>
```

Example Response

```
instrument<SP> tetra<TAB>
version<SP> sep 30 2004<TAB>
function<SP> d70, initialization halted because tips are loaded. please
verify that the appropriate tip rack (96,384,1536)is loaded and select
reset <TAB>
inst_status<SP> idle, remote, online<TAB>
exp_running<SP>no<TAB>
readplate<SP> default384, have<TAB>
sourceplate1<SP>default384, have<TAB>
sourceplate2<SP>none, have<TAB>
sourceplate3<SP>none, have<TAB>
loadplates<SP>none<TAB>
removeplates<SP>none<TAB>
platehandler<SP>idle<TAB>
plate_on<SP>no<TAB>
camera_temp<SP>ok, -35, -35<TAB>
chamber_temp<SP>off, 0, 0<TAB>
waste_bottlea<SP>ok<TAB>
```

```
waste_bottleb<SP>ok<TAB>
fill_bottlea<SP>ok<TAB>
fill_bottleb<SP>ok<TAB>
tips<SP>on<TAB>
outer_auto_door<SP>closed<CR>
```

Suggested Steps in Command Usage

- Send **status** command.
Response is the instrument status.

Loadplate<TAB>Location, Last Plate[, BAR CODE] <CR>

Command Description

Loads plate or tips to the specified location from landing pad. The plates or tips must have been placed on the instrument landing pad prior to the command being issued. Up to 13 plates (1 source plate, 12 compound plates) can be used in single experiment.

Additional parameters are available to provide the bar code for the delivered plate and identify the final plate for an assay.

The bar codes of up to 5 plates can be recorded in the data file.

Command Parameters

There are four parameters available for use with this command.

The parameters are location, bar code, and last plate. The location parameter is required. The bar code and last plate parameters are optional. If the bar code parameter is used then the plate order parameter must also be used.

Location parameter must be one of following:

- re—which indicates read plate location.
- s1—which indicates source plate 1 location.
- s2—which indicates source plate 2 location.
- s3—which indicates source plate 3 location.
- ti—which indicates tips loading location.

The tips and source 1 location are the same position. Either the “ti” or “s1” parameter can be used to indicate this position.

Bar code parameter must be:

- String up to 20 characters.
- String can have letters, numbers, underscore, and spaces.

The instrument always attempts to read a bar code while loading a plate. If a bar code is supplied through the loadplate command, any bar code actually found on the actual plate loaded is ignored.

The last plate parameter indicates whether or not the plate being loaded is the last plate in the current experiment. This parameter must be “yes” or “no”. The last plate parameter is no longer used by the FLIPR Penta System and ScreenWorks Software but continues to be required for compatibility reasons.

Example Command Line

```
loadplate<TAB>re, no, abcd1234<CR>
```

Example Response

```
ok<CR> or c10, error string<CR>
```

Suggested Steps in Command Usage

1. Check if instrument status is acceptable, such as no error state, no plate on landing pad, no plate in target location, plate handler is idle.
2. Put plate on landing pad.
3. Send **loadplate** command.
Response is **OK**, confirming command started execution or Condition code related to command, that confirming command could not start execution.
4. Monitor plate handler status for completion, and error status for any other kind of error.

Command Specific Error Responses

- Instrument is not functionalCondition code is C00
- Invalid plate locationCondition code is C05
- Plate handler busyCondition code is C10
- Landing pad does not have a plateCondition code is C20
- Plate location already has a plateCondition code is C15

Removeplate<TAB>location <CR>

Command Description

Removes plate or tips from the specified location. If source plate3 location is set for tips in the current experiment, the instrument assumes tips are present in that position.

Command Parameters

There is one parameter, location, for this command.

Location parameter must be one of following:

- re—which indicates read plate location.
- s1—which indicates source plate 1 location.
- s2—which indicates source plate 2 location.
- s3—which indicates source plate 3 location.
- ti—which indicates tips loading location.



Note: The tips and source 1 location are the same position. Either the “t1” or “s1” parameter may be used to indicate this position.

Example Command Line

```
removeplate<TAB>s1<CR>
```

Example Response

```
ok<CR> or c20, error string <CR>
```

Suggested Steps in Command Usage

1. Check if instrument status is acceptable (no error state, no plate on landing pad, plate present in target location, plate handler is idle).
2. Send **removeplate** command.
Response is **OK**, that confirming command started execution or Condition Code related to command, that confirming command could not start execution.
3. Monitor plate handler status for completion, and error status for any other kind of error.

Command Specific Error Responses

- Instrument is not functionalCondition code is C00
- Invalid plate locationCondition code is C05
- Plate handler busyCondition code is C10
- Landing pad has a plateCondition code is C21
- Plate location doesn't have plateCondition code is C16

Openprotocol<TAB>File Name<CR>

Command Description

This command opens the specified protocol. Any documents open before this command is sent close automatically. Only one protocol can be open at the time.

Command Parameters

There is one parameter, file name, for this command.

File name parameter is the complete file name of the protocol to be opened, including path and file type identifier (*.fmp).

Example Command Line

```
openprotocol<TAB>c:\screenworks\protocols\myprotocol.fmp<CR>
```

Example Response

```
ok<CR> or c30, error string <CR>
```

Suggested Steps in Command Usage

1. Make sure experiment is not running.
2. Send **openprotocol** command.

Response is **OK**, that confirming command executed or Condition code related to command, that confirming command did not execute.

Command Specific Error Responses

- Experiment is runningCondition code is C25
- Protocol not foundCondition code is C30
- Protocol could not be openCondition code is C35
- Invalid protocol nameCondition code is C36

Findprotocols<TAB>Folder<CR>

Command Description

This command returns all the protocols in the specified folder. The command returns a comma-delimited list containing all of the *.fmp type files saved in the indicated directory. Each of these files are NOT checked to confirm that the protocols that they contain are valid for the current instrument configuration.

Command Parameters

There is one parameter, folder, for this command.

The folder parameter identifies the folder for which a list of all *.fmp type files are returned. The folder must be complete and in the correct Windows format, including the drive identifier and a closing backslash(\).

Example Command Line

```
findprotocols<TAB>c:\screenworks\protocols\<CR>
```

Example Response

```
[myprotocol1.fmp,myprotocol2.fmp,myprotocol3.fmp]<CR> or c45, error string<CR>
```

Suggested Steps in Command Usage

- Send **findprotocol** command at any time.
Response is a list of protocols separated by comma or Condition code related to command, that confirming command could not execute.

Command Specific Error Responses

- Folder does not existCondition code is C45
- Invalid folder nameCondition code is C46
- Current document is not valid protocolCondition code is C50
- No protocols found in dirCondition code is C56

Runexperiment<TAB>[Data File Name]<CR>

Command Description

This command runs the currently open protocol, if it is valid for the current instrument configuration.

Running any protocol creates a data file and shifts the software focus to this data file for the duration of the experiment. When running in Remote mode, the data file is closed automatically when the experiment is done and the focus automatically returns to the protocol file.

Command Parameters

There is one optional parameter, data file name, for this command.

If a data file name parameter is used with this command, the data file created by running the current protocol is named data file name. The filename is appended to the **Data** folder path specified in the protocol Settings screen in the ScreenWorks Software. Never specify data file name containing a full path.

If the data file name parameter is not given, the data file created is saved with the auto save name as indicated in the Settings screen in the ScreenWorks Software.

Example Command Line

```
runexperiment<TAB>mydata10312003.fmd<CR>
```

Example Response

```
ok<CR> or c50, error string <CR>
```

Suggested Steps in Command Usage

1. Check if instrument is functioning fine.
2. Make sure the experiment is not running.
3. Open protocol first by sending **openprotocol** command.
4. Send **runexperiment** command.

Response is **OK**, that confirming command started execution or Condition code related to command, that confirming command could not start execution.

5. Monitor status for completion, and error status for any other kind of error.

Command Specific Error Responses

- Instrument is not functionalCondition code is C00
- Experiment is runningCondition code is C25
- Current document is not valid protocolCondition code is C50
- Invalid data file nameCondition code is C47

Stopexperiment<CR>

Command Description

This command stops the current action being performed by the instrument.

The **stopexperiment** command stops the experiment immediately. This can halt the instrument pipettor in the midst of a motion. In this case, the instrument immediately enters a fault state and an instrument reset is required. In addition, the pipettor might not be in motion but might have tips that contain fluid or are in a source or target plate. In these cases, it might be possible to recover without resetting by washing the tips. Regardless, an instrument reset is recommended following the use of any **stopexperiment** command.

Command Parameters

There are no parameters associated with this command.

Example Command Line

```
stopexperiment<CR>
```

Example Response

```
ok<CR> or c55, error string <CR>
```

Suggested Steps in Command Usage

1. Check if instrument is functioning fine.
2. Make sure experiment is running.
3. Send **stopexperiment** command.

Response is **OK**, confirming that command execution has begun, or a Condition code related to the command indicating that execution could not begin.

4. Monitor status for completion, and error status for any other kind of error.

Command Specific Error Responses

- Instrument is not functionalCondition code is 00
- No experiment is runningCondition code is 55

Clearerror<CR>

Command Description

Cleans error in case recoverable error occurred.

Command Parameters

There are no parameters associated with this command.

Example Command Line

```
clearerror<CR>
```

Example Response

```
ok<CR>
```

Suggested Steps in Command Usage

1. Check instrument status.
2. Check if error is occurred.
3. Send **clearerror** command.

Response is **OK** confirming the software has attempted to clear the error.

4. Monitor status for error is cleared.

Loadtips<CR>

Command Description

This command loads the tips onto the pipettor head from position s1.

This command executes when commanded if tips are not already loaded and any plate is in the tips position. The instrument is not capable of confirming that the plate in the tip load position is actually a tip rack. It is the responsibility of the user to ensure that the plate in that position is actually a tip rack. Attempting to load tips when a plate, rather than a tip rack, is in the tip load position results in an instrument fault and can result in damage to the plate, the instrument, or both.

Command Parameters

There are no parameters associated with this command.

Example Command Line

```
loadtips<CR>
```

Example Response

```
ok<CR>
```

Suggested Steps in Command Usage

1. Make sure an experiment is not running.
2. Check the instrument status to make sure tips are not already loaded.
3. Place tiprack in tipload position via the **loadplate** command.
4. Send **loadtips** command.

Response is **OK**, confirming that command execution has begun, or a Condition code related to the command indicating that execution could not begin.

5. Monitor status for completion, and error status for any other kind of error.

Command Specific Error Responses

- Can't load tips, tips on pipettorCondition code is C58

Unloadtips<CR>

Command Description

This command unloads the pipette tips from the pipettor head into an empty tip rack placed into the tip load position.

This command executes when commanded if tips are loaded and any plate is in the tips position. The instrument is not capable of confirming that the plate in the tip load position is actually an empty tip rack. It is the responsibility of the user to ensure that the plate in that position is actually an empty tip rack. Attempting to load tips when a plate or full tip rack, rather than an empty tip rack, is in the tip load position likely results in an instrument fault and can result in damage to the plate, the instrument, or both.

Command Parameters

There are no parameters associated with this command.

Example Command Line

```
unloadtips<CR>
```

Example Response

```
ok<CR>
```

Suggested Steps in Command Usage

1. Make sure an experiment is not running.
2. Check the instrument status to make sure tips are currently loaded.
3. Ensure that an empty tip rack is in tip load position.
4. Send **loadtips** command.

Response is **OK**, confirming that command execution has begun, or a Condition code related to the command indicating that execution could not begin.

5. Monitor status for completion, and error status for any other kind of error.

Command Specific Error Responses

- Can't unload tips, tips not on pipettorCondition code is C59

Cyclecameratemper<CR>

Command Description

This operation cycles the camera temperature by setting the camera setpoint to 20°C (68°F), waiting for the camera to reach this setpoint, resetting the camera to -45°C (-49°F), and then waiting for the camera to return to this temperature.

This command can be useful when transitioning from fluorescence to luminescence experiments. Cycling the camera temperature allows the camera CCD chip to recover from the high light levels present during fluorescence experiments and improves data quality in subsequent luminescence experiments.

This command requires up to 15 minutes to complete execution.

Command Parameters

There are no parameters associated with this command.

Example Command Line

```
cyclecameratemper<CR>
```

Example Response

```
ok<CR>
```

Suggested Steps in Command Usage

1. Make sure an experiment is not running.
2. Send **cyclecameratemper** command.
Response is **OK**, confirming that command execution has begun, or a Condition code related to the command indicating that execution could not begin.
3. Monitor status for completion, and error status for any other kind of error.

Tempcontrolonoff<TAB>Temp<CR>

Command Description

This command sets the temperature setting for the instrument chamber. After sending this command, the robotic controller should monitor the instrument temperature to confirm that the instrument reaches the setpoint. Some protocols do not execute until the required temperature is reached.

Example Command Line

```
tempcontrolonoff<TAB>12<CR>
```

Example Response

ok<CR>

Suggested Steps in Command Usage

1. Make sure an experiment is not running.
2. Send **tempcontrolonoff** command.
Response is **OK**, confirming that command execution has begun, or a Condition code related to the command indicating that execution could not begin.
3. Monitor status until the instrument reports that it has reached the requested temperature.

Command Specific Error Responses

- Temperature Control System is not functioningC60

Washtips<TAB>Fluid Type, Wash Cycles, Volume/Stroke, Aspirate Speed, Pump Speed, Strokes, Hold Time, Dispense Speed<CR>

Command Description

This command is the equivalent of selecting **Instrument> Manual Operation > Wash Tips** from the main menu of the ScreenWorks Software. This selection washes the pipette tips with the parameters provided.

Command Parameters

There are eight required parameters for this command. All eight parameters must be supplied for this command to properly execute. For some of the parameters, acceptable values can vary depending on the head type currently installed on the instrument. The following table lists the parameters, their purpose, and the appropriate range for the commands.

| Parameter | Description | Acceptable values or range | | |
|---------------|--|----------------------------|---------------|----------------|
| | | 96-well head | 384-well head | 1536-well head |
| Fluid type | Indicates which wash fluid bottle to be used. | a, b | | |
| Wash Cycles | Number of wash reservoir fill/drain cycles to complete. The requested # of strokes is repeated for each cycle. | 1,2,3,4,5 | | |
| Volume/stroke | Volume to be drawn into the pipette tip in each stroke. | 5–206 (double) | 1–28 (double) | 1–3 (double) |

| Parameter | Description | Acceptable values or range | | |
|----------------|--|--|---|---|
| | | 96-well head | 384-well head | 1536-well head |
| Aspirate speed | Pipettor aspiration speed in micro liters per second. | 5.00, 10.00, 20.00, 30.00, 40.00, 50.00, 75.00, 100.00, 125.00 | 0.50, 1.00, 2.00, 5.00, 7.00, 10.00, 15.00, 20.00, 25.00, 30.00 | 0.25, 0.50, 0.75, 1.00, 1.25, 1.50, 2.00, 3.00, 5.00, 10.00 |
| Pump speed | Speed to run the pumps which fill and drain the wash reservoir. Two speeds, fast and slow, are available. Slow is recommended for use with wash solvents which tend to bubble. | fast, slow | | |
| Strokes | Number of times which the pipettor fills and empty with the requested volume in each wash cycle. | 1–20 (integer) | | |
| Hold time | Pause time in seconds at the top of the stroke of the pipettor | 0–15 | | |
| Dispense speed | Pipettor dispense speed in micro liters per second. | 5–200 (double) | 1–50 (double) | 1–10 (double) |

Example Command Line

```
washtips<TAB>a,2,28.0,20.00,fast,5,0.0,20.0<CR>
```

Example Response

```
ok<CR>
```

Suggested Steps in Command Usage

1. Make sure an experiment is not running.
2. Ensure that tips are loaded.
3. Send **washtips** command.

Response is **OK**, confirming that command execution has begun, or a Condition code related to the command indicating that execution could not begin.

4. Monitor status until the wash is complete.

Command Specific Error Responses

- Tips not loaded C68

Configuration<CR>

Command Description:

The **configuration** command returns the configure of the instrument. The configuration response is an ASCII text string which includes:

- Information about the instrument type and firmware version.
- Pipettor head's type.
- Tip washer's type.
- Bar code reader installed or not.
- Plate handler installed or not.
- Cell Reservoir installed or not.
- Camera type installed.
- Chiller installed or not.
- LEDs installed.
- Emission filters' wavelength.

Example Command Line

```
configuration<CR>
```

Response format

```
instrument<SP>name<TAB>
version<SP>version#<TAB>
pipettor<SP>96|384|1536|384pintool| 1536pintool|no<TAB>
tip_washer<SP>96|384|1536|no<TAB>
bar_code<SP>installed|not installed<TAB>
plate_handler<SP>installed|not installed<TAB>
cell_reservoir<SP>installed|not installed<TAB>
camera_installed<SP>emccd|iccd<TAB>
chiller_installed<SP>installed|not installed<TAB>
wave1<SP>no led|470-495 nm|510-545 nm|390-420 nm|420-455
nm|610-626 nm|360-380 nm|525-550 nm|525-570 nm|570-595
nm|590-614 nm|620-648 nm|495-505nm|360-380 nm <TAB>
wave2<SP>no led|470-495 nm|510-545 nm|390-420 nm|420-455
nm|610-626 nm|360-380 nm|525-550 nm|525-570 nm|570-595
nm|590-614 nm|620-648 nm|495-505 nm|360-380 nm <TAB>
emfilt1<SP>no filter|515-575 nm|565-625 nm|440-480 nm|475-535
nm|646-706 nm|400-460 nm|570-630 nm|590-650 nm|615-675
nm|634-694 nm|668-728 nm|526-585 nm|400-460 nm<TAB>
emfilt2<SP>no filter|515-575 nm|565-625 nm|440-480 nm|475-535
nm|646-706 nm|400-460 nm|570-630 nm|590-650 nm|615-675
nm|634-694 nm|668-728 nm|526-585 nm|400-460 nm<TAB>
emfilt3<SP>no filter|515-575 nm|565-625 nm|440-480 nm|475-535
nm|646-706 nm|400-460 nm|570-630 nm|590-650 nm|615-675
nm|634-694 nm|668-728 nm|526-585 nm|400-460 nm<TAB>
<CR>
```

Example Response

```
Instrument<SP>tetra<TAB>
version<SP>2.1.0.0<TAB>
pipettor<SP>96<TAB>
tip_washer<SP>96<TAB>
bar_code<SP>installed<TAB>
plate_handler<SP>installed<TAB>
cell_reservoir<SP>installed<TAB>
camera_installed<SP>emccd<TAB>
chiller_installed<SP>installed<TAB>
wave1<SP>470-495 nm<TAB>
wave2<SP>510-545 nm<TAB>
emfilt1<SP>515-575 nm<TAB>
emfilt2<SP>515-575 nm<TAB>
emfilt3<SP>515-575 nm<CR>
```

Suggested Steps in Command Usage

- Send configuration command.
Response is instrument configuration.

Statusex<CR>

Command Description

The **statusex** command is an extension of status command. It returns the status of the instrument in more details than the status command. The **statusex** response is an ASCII text string which includes all status command response information plus:

- Tip washer status
- Cell Reservoir status
- Camera chiller status
- Camera intensifier status

Example Command line

```
statusex<CR>
```

Response format

```
instrument <SP>instrument name<TAB>
version<SP>mm/dd/yyyy<TAB>
function<SP>ok|errorcode|fatalcode, [errorstring]<TAB>
inst_
status<SP>busy|fault|error|resetting|abort|idle, remote|manual, offline|online<TAB>
AB>
exp_running<SP>yes|no<TAB>
readplate<SP>platenam | none, have|empty<TAB>
sourceplate1<SP>platenam | none, have|empty<TAB>
```

```

sourceplate2<SP>platename|none,have|empty<TAB>
sourceplate3<SP>platename|none,have|empty<TAB>
loadplates<SP>re[,s1,s2,s3,ti]|none<TAB>
removeplates<SP>re[,s1,s2,s3,ti]|none<TAB>
platehandler<SP>idle|busy_load|busy_remove<TAB>
plate_on<SP>yes|no<TAB>
camera_temp<SP>ok|not_ok|off|busy|fault,temp,setpoint<TAB>
chamber_temp<SP>ok|not_ok|fault|off|busy<TAB>
waste_bottlea<SP>ok|full<TAB>
waste_bottleb<SP>ok|full<TAB>
fill_bottlea<SP>ok|empty<TAB>
fill_bottleb<SP>ok|empty<TAB>
tips<SP>on|off<TAB>
outer_auto_door<SP>open|closed<TAB>
tip_washer<SP>ok|not_ok|off|busy|fault<TAB>
cell_reservoir<SP>ok|not_ok|off|busy|fault,stir_rate<TAB>
chiller_status<SP>ok|not_ok|off|busy|fault,temp,setpoint<TAB>
intensifier_status ok|not_ok|off|busy|fault<CR>

```

Example Response

```

instrument<SP> tetra<TAB>
version<SP>2.1.0.0<TAB>
function<SP>ok<TAB>
inst_status<SP>idle,remote,online<TAB>
exp_running<SP>no<TAB>
readplate<SP>default96,have<TAB>
sourceplate1<SP>default96,have<TAB>
sourceplate2<SP>default96,have<TAB>
sourceplate3<SP>default96,have<TAB>
loadplates<SP>none<TAB>
removeplates<SP>none<TAB>
platehandler<SP>idle<TAB>
plate_on<SP>no<TAB>
camera<SP>temp ok,0,0<TAB>
chamber_temp<SP>off,0,0<TAB>
waste_bottlea<SP>ok<TAB>
waste_bottleb<SP>ok<TAB>
fill_bottlea<SP>ok<TAB>
fill_bottleb<SP>ok<TAB>
tips<SP>on<TAB>
outer_auto_door<SP>open<TAB>
tip_washer<SP>ok<TAB>
cell_reservoir<SP>not_ok,0<TAB>
chiller_status<SP>ok,18.0, 18.0<TAB>
intensifier_status<SP>ok<CR>

```

Cellflaskcontrol<TAB>Rate<CR>

Command Description

This command sets the stir-rate for the cell-flask control. After sending this command, the robotic controller should monitor the instrument stir-rate to confirm that the instrument reaches the setpoint.

Command Parameters

There is one required parameter, rate, for this command.

If a value of zero is used for the rate parameter, the cell flask control stops stirring. If a valid value for rate is used, the cell flask control turns on and stirs at this rate. Valid values for rate are integers between 1–25.

Example Command Line

```
cellflaskcontrol<TAB>22<CR>
```

Example Response

```
ok<CR>
```

Suggested Steps in Command usage

1. Make sure an experiment is not running.
2. Send **cellflaskcontrol** command.
Response is **OK**, confirming that command execution has begun, or a Condition code related to the command indicating that execution could not begin.
3. Monitor status until the instrument reports that it has reached the requested stir-rate.

Command Specific Error Responses

- C76–Cell flask stirring rate parameter was not sent
- C77–Parameter is out of range. Range is ...

Washcellreservoir<TAB>Fluid Type, Fill Speed, Drain Destination, Drain Speed, Wash Cycles, Hold Time, Volume Level<CR>

Command Description

This command is the equivalent of selecting **Instrument > Manual Operation > Wash Cell Reservoir** from the main menu of the ScreenWorks Software. This selection washes the Cell Reservoir with the parameters provided.

Command Parameters

There are seven required parameters for this command. All seven parameters must be supplied for this command to properly execute. The following table lists the parameters, their purpose, and the appropriate range for each parameter.

| Parameter | Type | Description | Acceptable values or range |
|-------------------|---------|---|---|
| Fluid type | string | Indicates the fluid bottle to be used. | Fluid 1, fluid 2, fluid 3, fluid 4, cell flask. |
| Fill speed | integer | Speed to run the pumps which fill the Cell Reservoir. Range from 1–10 where 1 is very slow, and 10 is very fast. | 1–10 |
| Drain destination | string | Indicate the bottle to drain fluid to. | Fluid 1, fluid 2, fluid 3, fluid 4, cell flask, waste a, waste b. |
| Drain speed | integer | Speed to run the pumps which drain the Cell Reservoir. Range from 1–10 where 1 is very slow, and 10 is very fast. | 1–10 |
| Wash cycles | integer | Number of wash reservoir fill/drain cycles to complete. | 1,2,3,4,5 |
| Hold time | float | Pause time in seconds, leaving the fluid in the reservoir for desired period of time before draining it. | 0–300 |
| Volume | string | This indicate the level of fluid to be filled in the Cell Reservoir. | low, high |

Example Command Line

```
washcellreservoir<TAB>fluid 2, 6, waste a, 8, 1, 0, high<CR>
```

Example Response

```
Ok<CR>
```

Suggested Steps in Command usage

1. Make sure an experiment is not running.
2. Ensure that tips are loaded.
3. Send **washcellreservoir** command.
Response is **OK**, confirming that command execution has begun, or a Condition code related to the command indicating that execution could not begin.
4. Monitor status until the wash is complete.

Command Specific Error Responses

- C77–Parameter is out of range. The range is ...
- C78–The number of parameters for wash Cell Reservoir commands is invalid.
- C79–Invalid volume parameter. It must be 'low' or 'high'.
- C72–Hold time is too small.
- C73–Hold time is too large.

Error Handling and Reporting

Interface Errors

Interface errors are errors that are generated by issuing a command. These errors are detected upon receipt of the command by ScreenWorks Software and are reported in the response to the command. These errors consist of a single alpha character and a two or three-digit code. Specific errors are listed with the command that can generate them in [Table G-1](#).

Table G-1: Interface Errors

| Error Code | Description |
|------------|---|
| C01 | Instrument is not in Remote Mode. Only 'status' command can be sent while instrument is in Manual Mode. |
| C05 | Invalid plate location. |
| C10 | Plate handler is busy. |
| C15 | Plate location already has a plate. |
| C16 | Plate location has no plate. |
| C20 | Landing pad has no plate. |
| C21 | Landing pad has plate. |
| C25 | Experiment is running. |
| C35 | Protocol could not be opened. |
| C36 | Invalid protocol name. |
| C45 | Folder does not exist. |
| C46 | Invalid folder name. |
| C47 | Invalid data file name. |
| C50 | Current document is not valid protocol.S |
| C55 | No experiment is running. |
| C56 | No protocols found in dir. |
| C57 | The temperature parameter (int) was not sent. |

Table G-1: Interface Errors (continued)

| Error Code | Description |
|------------|---|
| C58 | Can't load tips, tips on pipettor. |
| C59 | Can't unload tips, no tips on pipettor. |
| C60 | The temperature control system is not functioning. |
| C61 | The aspirate speed is out of range. |
| C62 | The pump speed can only be Slow or Fast. |
| C63 | Invalid number of wash strokes. |
| C64 | Invalid number of parameters for the wash tips command. |
| C65 | Invalid number of wash cycles. |
| C66 | Invalid protocol parameters discovered at runtime. |
| C67 | Invalid wash fluid parameter. |
| C68 | Can't wash tips, tips are not on. |
| C69 | Unrecognized or badly formatted command. |
| C70 | Volume stroke too large. |
| C71 | Volume stroke too small. |
| C72 | Hold time is too small. |
| C73 | Hold time is too large. |
| C74 | Dispense speed is too large. |
| C75 | Dispense speed is too small. |
| C76 | Cell flask stirring rate parameter was not sent. |
| C77 | Parameter is out of range. |
| C78 | The number of parameters for wash Cell Reservoir commands is invalid. |
| C79 | Invalid volume parameter. It must be 'low' or 'high'. |

Table G-2 documents the errors one might encounter when sending the **runexperiment** command. The errors explain how the currently opened protocol does not match instrument configure or instrument is not ready to run current protocol.

Table G-2: runexperiment Command Errors

| Error Code | Description |
|------------|--|
| E100 | Invalid LED filter combination! Calibrate it or choose valid combination. |
| E110 | Plate does not exist in plate library. Choose existing plate or add it into plate library. |
| E111 | Plate is not present or is not in the proper location. |

Table G-2: runexperiment Command Errors (continued)

| Error Code | Description |
|------------|---|
| E112 | Door(s) are opened. Close the doors before start your experiment. |
| E113 | Load Tips position not checked! Check Load Tips position for auto load tips or manually load tips before running experiment. |
| E114 | The chamber temperature is not in range. |
| E115 | Experiment wash tips and there is no tip washer. |
| E116 | Bottle A is empty! Refill it or use another bottle. |
| E117 | Bottle B is empty! Refill it or use another bottle. |
| E118 | Waste bottle A is full. Please empty the bottle to continue. |
| E119 | Waste bottle B is full. Please empty the bottle to continue. |
| E120 | Remove plates redundant. |
| E121 | Error relate to the data folder. The path is not valid, or cannot be accessed. Or there are not enough disk space. |
| E122 | Error relate to the export folder. The path is not valid, or cannot be accessed. Or there are not enough disk space. |
| E123 | The read interval is too short. Please check settings in Read views. |
| E124 | Invalid aspirate parameters [in the protocol]. |
| E125 | Invalid dispense parameters [in the protocol]. |
| E126 | The wash tips parameters are invalid. |
| E127 | Invalid fluid bottle specified in wash tips process. |
| E128 | Invalid Read Mode configuration. You tried to take fluorescence (with excitation) and luminescence (without excitation) reading in one protocol. This is not allowed. |
| E129 | Invalid Read Mode configuration. Camera gain or gate time is out of range. Please check your protocol settings. |
| E130 | The protocol has transfer cells and/or wash Cell Reservoir operations even though the Cell Reservoir is not installed. |
| E131 | Cell Reservoir is not ready. |
| E132 | Cell flask spinning rate could not be set. Please check your instrument status. |

Instrument Errors

Instrument errors are errors that are generated by executing a command. These errors are detected during the execution of the command by the instrument. They are reported in the status response in the function line. These errors consist of a three-digit numeric code.

Instrument errors fall into two general categories, recoverable and fatal errors. When fatal errors occur, the instrument can not continue to do any operation. It aborts any pending operations. When recoverable errors occur with user interaction, the instrument can continue.

The **status** command returns `ERRORCODE` or `FATALCODE`.

- Error codes 100–199 are call Molecular Devices Tech support errors.
- Error codes 200–299 are instrument recoverable errors.
- Error codes 300–399 are instrument non-fatal errors.
- Error codes 400–499 are instrument fatal errors.
- Error codes 500–599 are instrument diagnostic errors.

A full list of these errors can be found in [Troubleshooting on page 209](#).

Remote Interface Revision History

V1.0

First release.

V1.1

Added the following new commands:

- `version`
- `loadtips`
- `unloadtips`
- `cyclecameratemp`
- `tempcontrolonoff`
- `washtips`

Error code response include text description after code, for example, `c45, folder does not exist`.

V1.2

Added Error Code:

`C01, Instrument is not in Remote Mode. Only status command can be sent while instrument is in Manual Mode.`

Bug Fix:

Commands made to the remote interface could sometimes take a long time and cause the remote interface to not respond to the client application within an acceptable timeframe. Client applications were able to work around this bug with massive TCP/IP timeouts (minutes) which made a true hang or network dropout undetectable. This workaround is no longer required.

The interface is unchanged, and if the client is correctly parsing the **status** response, no changes should be required in the client application, with the exception of reducing the TCP/IP timeout back to an acceptable value, for example, < 10 seconds. As always, rigorous regression testing of client code is encouraged.

V1.3

Added the following new commands to support new functionality in the Tetra Plus:

- `statusex`
- `configuration`
- `cellflaskcontrol`
- `washcellreservoir`



Note: Tips location has changed from the **Source 3** location to the **Source 1** location.

Barcode Specifications

About Bar Codes

The FLIPR Penta System bar code scanner is an OEM module with decode capability to most commonly used bar code symbols in industry worldwide. The list includes:

- WPC (UPC /EAN/JAN)
- IATA
- Code 39
- Interleaved 2 of 5
- Industrial 2 of 5
- Code 128
- Code 93
- MSI/Plessey
- Codabar (NW-7)

Bar Code Recommendations

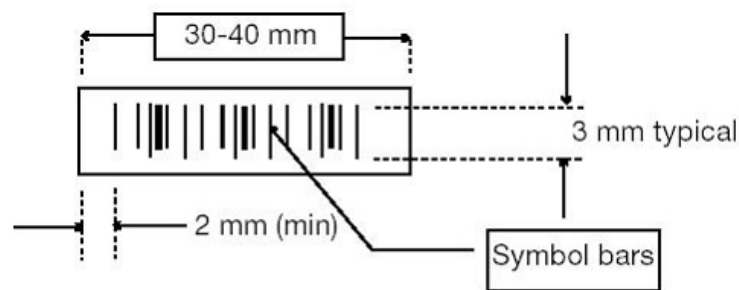
The bar code reader implementation allows a bar code read down to 7.5 mil (0.20 mm) on photographic-quality symbols and ambient light < 450 ft-candles.

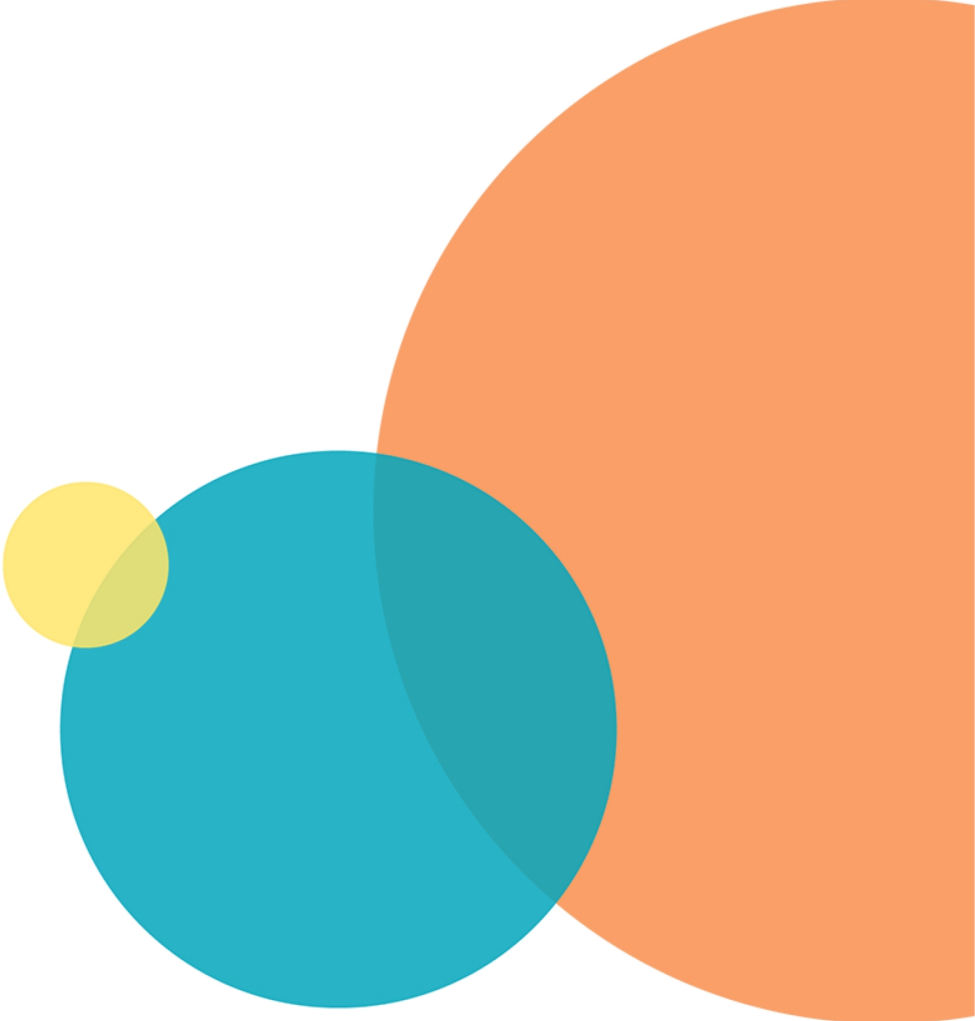
We recommend, however, the user stay above 7.5 mil (0.26 mm) since most bar code printers might not have the resolution needed to match the required aspect ratio for a minimum of 3 mm-high symbol lines. Before adjusting the bar code scanner to the bar code label location on the microplate, some general guidelines to the preparation and quality of the bar code label need to be followed for optimum performance:

- The minimum height or size of the symbol bars depends on the code symbol and the density used. Try to keep the height of the symbol bars to more than 0.12 in. (3.0 mm).
- Do not change the aspect ratio required by the label printer. Do not force the bar code symbol size to fit on your label space i.e., maintain the aspect ratio free of distortion.

- Align the bars horizontally (side-to-side) and vertically (top-to-bottom) if possible, to keep space available on the bar code label.
- Keep a few millimeters (2–4) of empty dead or silent zone on both ends of the code-symbol lines. Keep all symbol lines away from the label corners and lifted edges. Optimize your characters so there is no need to cram a lot of lines on the label.
- Keep the longitudinal dimension of the labels/symbols to less than 1.25–1.50 in. (30–40 mm).
- Choose the media (label substrate material or paper) print or transfer ink to give good contrast and optimal reflection. Metalized, very shiny or clear background labels may be less reliable. Test a sample if using a new color or surface material.

Bar Code Specifications





Appendix H: Electromagnetic Compatibility



H

Regulatory for Canada (ICES/NMB-001:2006)

This ISM device complies with Canadian ICES-001.

Cet appareil ISM est conforme à la norme NMB-001 du Canada.

ISM Equipment Classification (Group 1, Class A)

This equipment is designated as scientific equipment for laboratory use that intentionally generate and/or use conductively coupled radio-frequency energy for internal functioning, and are suitable for use in all establishments, other than domestic and those directly connected to a low voltage power supply network which supply buildings used for domestic purposes.

Information to the User (FCC Notice)

This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to part 18 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at their own expense. Changes or modifications made to this equipment not expressly approved by the party responsible for compliance may void the FCC authorization to operate this equipment.

Contact Us

Phone: [+1-800-635-5577](tel:+1-800-635-5577)
Web: moleculardevices.com
Email: info@moldev.com

Visit our website for a current listing of worldwide distributors.