



# ClonePix 2

Mammalian Colony Picker

Hardware and Software

## User Guide

## ClonePix 2 Mammalian Colony Picker 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.  
©2024 Molecular Devices, LLC.  
All rights reserved.



# Contents

|   |           |
|---|-----------|
| <b>Safety Information</b> .....                   | <b>6</b>  |
| Warnings, Cautions, Notes, and Tips .....         | 6         |
| Symbols on the Instrument .....                   | 7         |
| Electrical Safety .....                           | 9         |
| External or Implanted Medical Device Safety ..... | 9         |
| Heat and Burn Safety .....                        | 9         |
| Chemical and Biological Safety .....              | 10        |
| Moving Parts Safety .....                         | 10        |
| <b>Chapter 1: Introduction</b> .....              | <b>11</b> |
| Workflows .....                                   | 11        |
| Before You Start .....                            | 14        |
| Software Installation .....                       | 15        |
| Powering On the System .....                      | 19        |
| Powering Off the System and the Software .....    | 19        |
| <b>Chapter 2: Instrument Overview</b> .....       | <b>21</b> |
| Process Area .....                                | 22        |
| Computer Connections .....                        | 24        |
| Front Panel Controls and Display .....            | 25        |
| Electrical Connections .....                      | 25        |
| Bottles Connections .....                         | 26        |
| Compressed Air Supply .....                       | 26        |
| Microplate Stackers .....                         | 27        |
| Filters .....                                     | 29        |
| Using Barcodes .....                              | 29        |
| Camera Alignment .....                            | 30        |
| Safety Features .....                             | 30        |
| <b>Chapter 3: Software Overview</b> .....         | <b>31</b> |
| Main Navigation Screen .....                      | 31        |
| Menu Options .....                                | 33        |
| <b>Chapter 4: Picking Processes</b> .....         | <b>35</b> |
| Prepare for Pick Run .....                        | 35        |
| Pick Run Process .....                            | 44        |
| Imaging Run Process .....                         | 73        |

|  |            |
|--|------------|
| Review Results Process .....   | 75         |
| <b>Chapter 5: Utility Processes .....</b>                                | <b>79</b>  |
| Picking Head Management .....  | 79         |
| Utility Process .....  | 81         |
| <b>Chapter 6: Maintenance and Troubleshooting .....</b>                  | <b>83</b>  |
| Doing Preventive Maintenance .....                                       | 83         |
| Changing Pins .....  | 84         |
| Cleaning the Instrument .....  | 85         |
| Cleaning the System Fluid Supply .....                                   | 86         |
| Cleaning the Head and Pins .....   | 87         |
| Cleaning the Wash and Waste Bottles .....                                | 90         |
| Cleaning the Wash and Waste Tubing .....                                 | 90         |
| Cleaning the Picking Tubing .....  | 91         |
| Cleaning the Ethanol Wash Bath .....                                     | 91         |
| Cleaning the Source Plate Imaging Cover .....                            | 92         |
| Draining the Compressor .....  | 92         |
| Moving and Storing the Instrument .....                                  | 94         |
| Commissioning Process .....  | 94         |
| Troubleshooting .....  | 94         |
| Obtaining Support .....  | 96         |
| <b>Appendix A: Technical Specifications .....</b>                        | <b>99</b>  |
| Instrument Dimensions .....  | 101        |
| <b>Appendix B: Plate Compatibility .....</b>                             | <b>103</b> |
| Plate Holders .....  | 103        |
| Lid Lifter Assembly .....  | 104        |
| <b>Appendix C: Replacement Parts, Accessories, and Consumables .....</b> | <b>105</b> |
| Replacement Parts .....  | 105        |
| Accessories and Consumables .....  | 105        |
| Reagents and Media .....   | 105        |
| <b>Appendix D: Imaging Definitions .....</b>                             | <b>107</b> |
| <b>Appendix E: Guide to Fluorophore Configurations .....</b>             | <b>111</b> |
| Measuring Fluorescence .....   | 111        |
| Example of Fluorescent Statistic Generation .....                        | 112        |
| Interior Intensity Statistics .....                                      | 112        |
| Exterior Intensity Statistics Calculation .....                          | 113        |
| Exterior Intensity Statistics .....                                      | 115        |

|  |            |
|--|------------|
| <b>Appendix F: ClonePix 2 Data Viewer Software</b> ..... | <b>117</b> |
| Software Prerequisites .....                             | 117        |
| Installing ClonePix 2 Data Viewer Software .....         | 117        |
| Licensing ClonePix 2 Data Viewer Software .....          | 117        |
| Using ClonePix 2 Data Viewer Software .....              | 118        |

Information about the safe use of the instrument from Molecular Devices includes an understanding of the user-attention statements in this guide, the safety labels on the instrument, precautions to follow before you operate the instrument, and precautions to follow while you operate the instrument.

Make sure that everyone involved with the operation of the instrument has:

- Received instruction in general safety practices for laboratories.
- Received instruction in specific safety practices for the instrument.
- Read and understood all Safety Data Sheets (SDS) for all materials being used.

Read and observe all warnings, cautions, and instructions. 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.

---










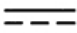

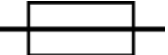














**Tip:** A tip provides useful information or a shortcut but is not essential to the completion of a procedure.

---

## Symbols on the Instrument

Each safety label on the instrument contains an alert symbol that indicates a type of potential safety hazard.

| Symbol  | Indication   |
|---|--|
|    | 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.  |
|    | Potential electrical-shock hazard from a high-voltage source. All safety instructions must be read and understood before proceeding with the installation, maintenance, and servicing of all modules. 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. |
|    | Potential lifting hazard. For information about the weight of the instrument,  |
|    | Potential pinch hazard.  |
|   | Potential heat hazard.   |
|  | Rotating parts hazard.   |
|  | Power on. See <a href="#">Powering On the System on page 19</a> .  |
|  | Power off. See <a href="#">Powering Off the System and the Software on page 19</a> .   |
|  | Alternating current.   |
|  | Direct current.  |
|  | Location of the protective earth (ground) terminal.  |
|  | Location of a fuse.  |

| Symbol  | Indication   |
|---|--|
|    | Instrument serial number.  |
|    | Instrument manufacture date.   |
|    | CSA certification.   |
|    | European technology conformity.  |
|    | United Kingdom technology conformity.  |
|    | South Korea technology conformity.   |
|    | Compliance with Australian radio communication requirements.   |
|   | Compliance with Chinese RoHS Pollution Control Requirements.   |
|    | <p>Compliance with the Waste Electrical and Electronic Equipment (WEEE) Directive of the European Union. You must not discard this electrical or electronic product or its components in domestic household waste or in the municipal waste collection system.</p> <p>For products under the requirement of the WEEE directive, contact your dealer or local Molecular Devices office for the procedures to facilitate the proper collection, treatment, recovery, recycling, and safe disposal of the device.</p> |
|    | There is an authorized representative in the European community.   |
|    | Instrument manufacturer.   |
| <p>Info for USA only: California Proposition 65</p>  <p><b>WARNING</b><br/>Cancer &amp; Reproductive Harm<br/><a href="http://www.P65Warnings.ca.gov">www.P65Warnings.ca.gov</a></p> <p style="font-size: small;">CH1063AA</p> | Compliance with California Proposition 65, which requires businesses to warn Californians about significant exposures to chemicals that cause cancer, birth defects or other reproductive harm.  |



## 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.

## External or Implanted Medical Device Safety

Persons with external or implanted medical devices must evaluate the risks related to these devices before entering an area where the instrument is in use. Keep magnetic storage devices or strips, such as hard drives and credit cards, away from the instrument.

## Heat and Burn Safety

The instrument is fitted with a high-temperature halogen dryer. The casing can become hot during the drying cycle.



**WARNING!** The casing of the halogen dryer can become hot during the drying cycle. Before touching this area, make sure that the dryer has had time to properly cool to a safe temperature.

---

## 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 door automatically 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.

In an emergency, press the Emergency Stop button on the front of the instrument to immediately stop all motion and turn off the instrument. Before you can restart the instrument, you must pull out the Emergency Stop button and then press the Start button.



**WARNING! Do not obstruct or otherwise prevent access to the Emergency Stop button.**

---

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.

ClonePix™ 2 software controls the process of selective mammalian cell picking by the ClonePix 2 system. The software is designed to simplify the day-to-day interaction with the instrument. It permits multiple users to select and collect colonies of suspension cells from semi-solid medium or adherent colonies from liquid medium. The colony detection feature contains powerful algorithms for selecting cell colonies based on physical characteristics and proprietary multi-channel fluorescence technology for detecting protein secretion or production.

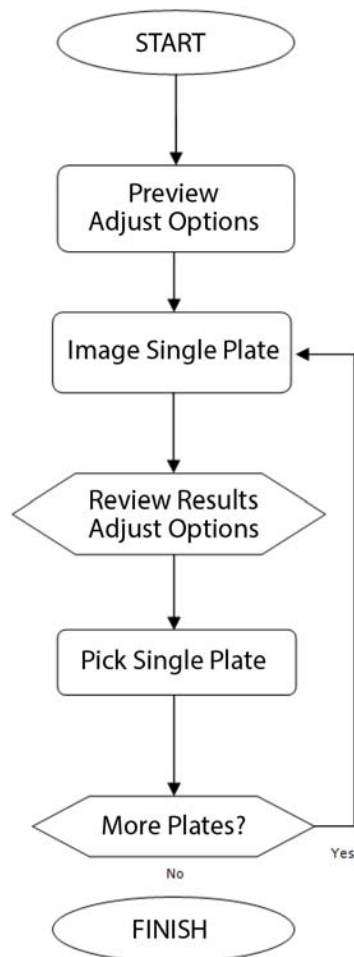
There are two software packages. ClonePix 2 software is required to run the instrument and ClonePix 2 Remote Data Viewer is available for purchase to access results remotely. For more information, contact Molecular Devices Technical Support. See [Obtaining Support on page 96](#).

## Workflows

ClonePix 2 software uses three different workflow processes: Pick Run, Batch Run, and Imaging Run.

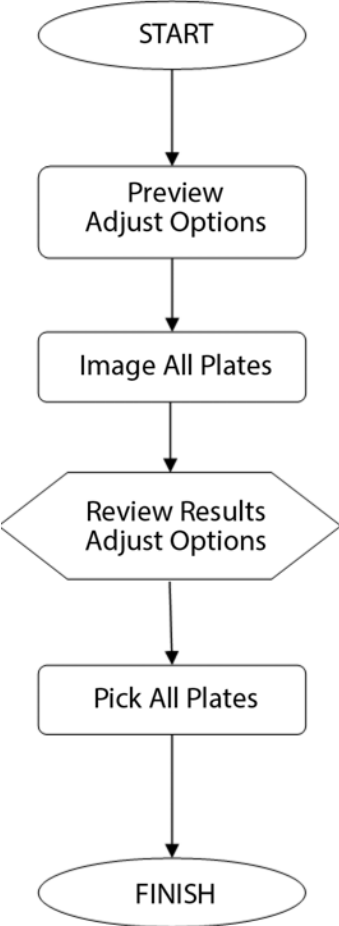
### Pick Run

The Pick Run processes images and picks the first plate, then the second plate, and so on, from however many plates you choose.



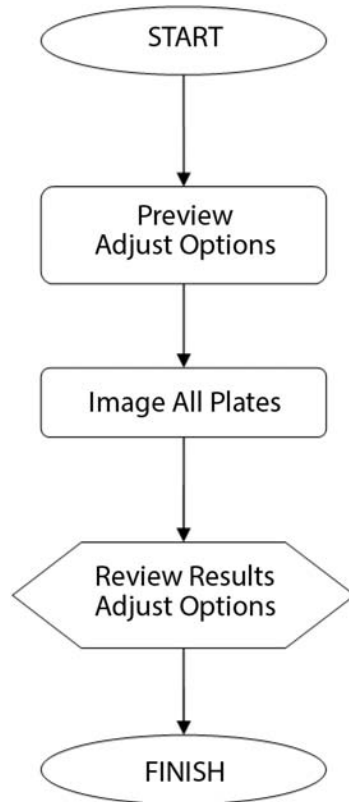
### Batch Run

The Batch Run process is set up within the Pick Run process. This process images all plates in the stacker cassette and then aggregates statistics for all plates before the picking of the selected clones from the entire batch.



## Imaging Run

Use the Imaging Run process to observe the status of plates during incubation or to monitor colony growth and secretion over time. The Imaging Run process images all plates without picking. After the plates are imaged, you can review the results in ClonePix 2 software or on another computer using the Remote Data Viewer Software. You can then re-analyze the images with different statistical criteria.



## Before You Start

Before you power on the ClonePix™ 2 Mammalian Colony Picker and start the software, do the following:

- Verify that the instrument and compressor are plugged in.
- Verify that the correct picking pins are installed in the picking head for the type of cells to be picked.



**Note:** Before a Pick Run, picking pins should be cleaned by sonication in a 2% solution of QuClean (K2505) and autoclaved. Also autoclave the Picking Pin Removal Key (X4948) at the same time. See [Sonicating the Pins on page 90](#). If the picking pins need to be changed, remove the Picking Head and swap the pins. See [Picking Head Management on page 79](#).



**Tip:** The picking pin type is engraved at the base of each pin. You might need to remove the pin from the head to confirm the pin type.

Picking pin types include:

- **F1 Picking Pins** (400 µm internal diameter; X4961) for suspension cell picking from semi-solid medium.



**Note:** These are the standard pins that come with the instrument.

- **F2 Picking Pins** (700 µm internal diameter; X4962) for adherent cell picking from liquid medium.
- Verify that instrument bed is clear of obstructions and loose items. See [Process Area on page 22](#).
- Verify that all motor tracks are free of obstruction.
- Verify that there are no obstructions to movement of the picking head.
- Verify that the main instrument axes, XY, are roughly in the center of their respective travel positions. The axes need to be away from their respective end-stops on startup so that they can reference themselves correctly to their home positions.
- Wipe the instrument bed with 70% ethanol using a lint-free cloth.
- Verify that no plates or cassettes are in the stackers.
- Verify that the Emergency Stop button on the front panel of the instrument is pulled out. See [Front Panel Controls and Display on page 25](#).
- Fill the ethanol feed bottle with 70% ethanol and empty the ethanol waste bottle, as needed.
- Separately autoclave the picking feed bottle and cap assembly, then fill the Picking Feed bottle with sterile deionized water and cap the bottle in a sterile tissue culture hood.
- Ensure that your picking waste bottle is empty before continuing.
- Verify that the enclosure front door is closed.
- Wait for approximately 2 minutes after the instrument is powered on before opening the software to ensure that all drives 'home' correctly. See [Powering On the System on page 19](#).



**WARNING!** For safety reasons the front door interlock should never be interfered with or overridden. If the door is opened at any stage, this triggers the actuator head or germicidal lamp to stop until the door is closed.

## Software Installation



**Note:** Molecular Devices installs the ClonePix 2 software and the software updates.

The ClonePix 2 software controls the ClonePix 2 Mammalian Colony Picker instrument and is based around the concept of the instrument running processes.

### Computer System Requirements

The ClonePix 2 software requires the following computer specifications:

#### Computer System Requirements

| Item              | Description   |
|-------------------|---|
| Operating system  | Windows 11<br>Windows 10, 64-bit<br>Windows 7, 32-bit |
| Memory            | 8 GB RAM or more                                      |
| Data Connection   | 10/100 Ethernet port                                  |
| Camera Connection | SBIG: USB 2.0 port<br>Basler Pylon: USB 3.0 port      |



**CAUTION!** Do Not replace the computer operating system with one of your own computers. The computer supplied with your system includes hardware components specifically configured to control your instrument.

### Backing Up the Previous Version Configuration

Before making any major computer changes, such as installing drivers or software, back up the current configuration.

#### Add a Windows System Restore Point

If you are using the Windows 7 operating system, do the following:

1. Open the computer Control Panel, go to **System and Security**, and then click the **System** link.
2. In the left panel, click the **System Protection** link.
3. In the System Properties dialog, click the **System Protection** tab and then click **Create**.
4. In the Create a Restore Point dialog, choose a name that you will remember, such as the current date or the name of a program you are about to install.
5. When *The restore point was created successfully* message displays, click **Close**.
6. In any other dialogs, click **Close** to finish.

If you use the Windows 10 or 11 operating system, do the following:

1. Open the computer search field, enter **CREATE A RESTORE POINT**, and select **Control panel**.
2. In the System Properties dialog, on the System Protection tab, select the main system drive, and click **Create**.
3. Enter a description to identify the restore point.
4. Click **Create**.
5. When the message *The restore point was created successfully* displays, click **Close**.
6. In any other dialogs, click **Close** to finish.

## Back up ClonePix 2 Software Configuration File

To back up your current configuration:

1. Click **Start > All Programs > Molecular Devices > ClonePix 2 software > Configuration Manager**.
2. In the Fusion Configuration Manager, click **Backup To File**.
3. In the Backup Fusion Configuration dialog, enter the version number of the currently installed software and the date and then click **OK**.
4. In the Save Configuration File dialog, locate and select the folder where you want to save the backup file, enter a file name, and then click **Save**.

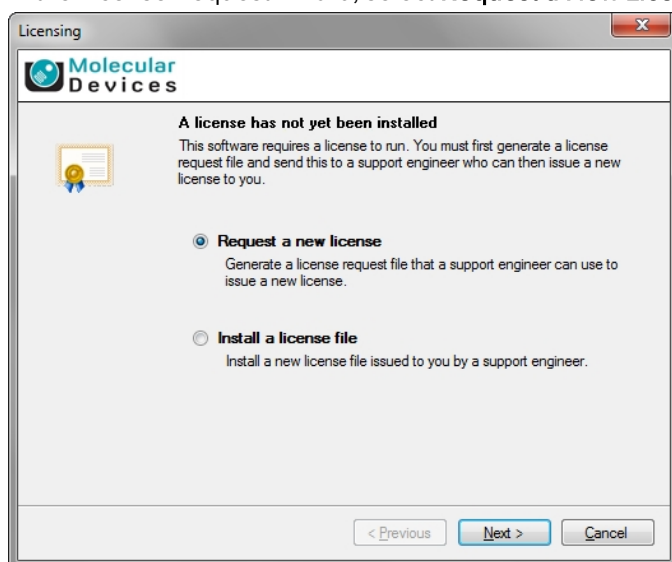
## Obtaining a License for the Software

The first time that you start the ClonePix 2 software after completing the installation, the software prompts you for a license.

### Requesting a Software License

To request a new software license:

1. In the License Message dialog, click **Yes**.
2. In the License Request wizard, select **Request a New License**, and then click **Next**.





3. Enter the requested information and click **Next**.

The screenshot shows a window titled "Licensing" with the Molecular Devices logo. The text reads: "Please provide the following details. This will help our support engineer to create a license for you." Below this are three input fields: "Registered User Name\*" with the value "User", "Company/Institute Name\*" with the value "Company", and "Instrument Serial Number (optional)". There is also a "Notes/Comments:" text area. At the bottom, it says "\*These fields are required" and has buttons for "< Previous", "Next >", and "Cancel".

4. To save the request file to a location on the computer, click **Save**.

The screenshot shows the same "Licensing" window. The text now reads: "Click the 'save' button to store the license request into a file. Take note of where you save the file to so that you can locate it later." A "Save..." button is centered on the screen. The "< Previous", "Next >", and "Cancel" buttons remain at the bottom.

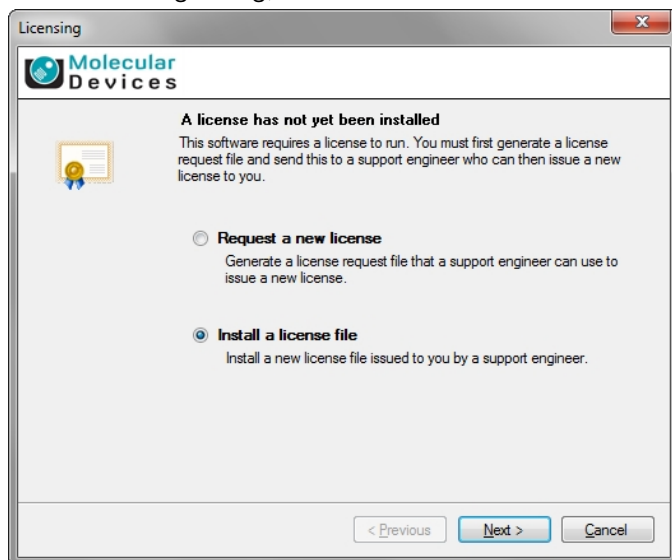
5. In the Save the Request to a File dialog, save the file in a location where you can find it.
6. To close the wizard, click **Finish**.
7. Change the **LicenseRequest.req** file name to a **LicenseRequest.txt** file, and then email it to Molecular Devices Technical Support. See [support.moleculardevices.com/](http://support.moleculardevices.com/).
8. After you receive the license file, save it on the computer where it is accessible, and then install the license.

## Installing a Software License

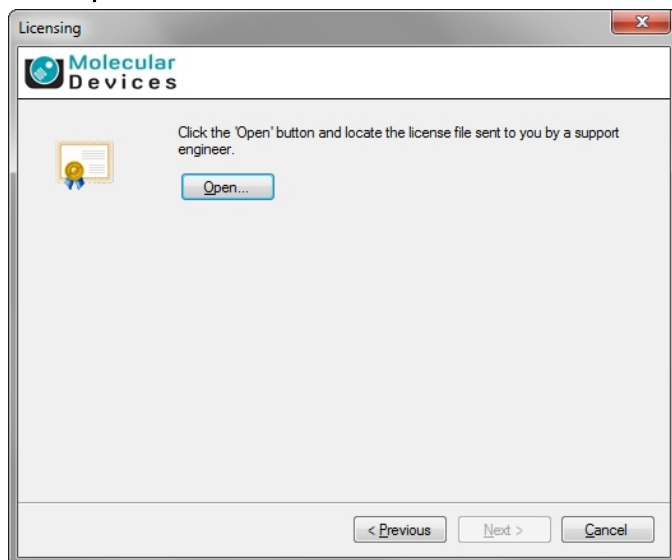
After you receive the license from Molecular Devices Technical Support, do the following:

To install a software license:

1. From the computer desktop, double-click the ClonePix 2 icon to start the software.
2. In the Licensing dialog, select **Install a License File** and click **Next**.



3. Click **Open**.



4. In the Select the License File to Install dialog, select the license file you saved and click **Open**.
5. After the license file installs, click **Finish**.

## Powering On the System

Before continuing, do as instructed in [Before You Start on page 14](#).

Every time you use the instrument the three axes sequentially run through their “Initialize drives” routine. This enables the drives to find their respective home positions. The system must complete this routine without interference to ensure that there is no damage to the instrument or its auxiliary equipment.

To power on the system:

1. Power on the compressor and ensure compressed air gauge is set to 80 psi (5.5 bar).
2. On the front panel of the instrument, push the **Start** button.

The Power On light illuminates.



**Tip:** If the power to the system does not turn on, make sure that the door is closed and the Emergency Stop button is pulled out.

---

The instrument cycles through various start-up processes indicated on the front indicator panel.

3. Power on the computer and wait for it to finish initializing.
4. From the computer desktop, double-click on the ClonePix 2 icon.

## Powering Off the System and the Software

To power off the ClonePix 2 Mammalian Colony Picker:

1. To exit the ClonePix 2 software, on the Main Navigation screen, select **File > Exit**.
2. To shut down the computer, click **Start > Shut down**, and then wait for computer to power off completely.
3. To power off the instrument, on the front of the instrument, press the **Stop** button.
4. Turn the power off at the mains.

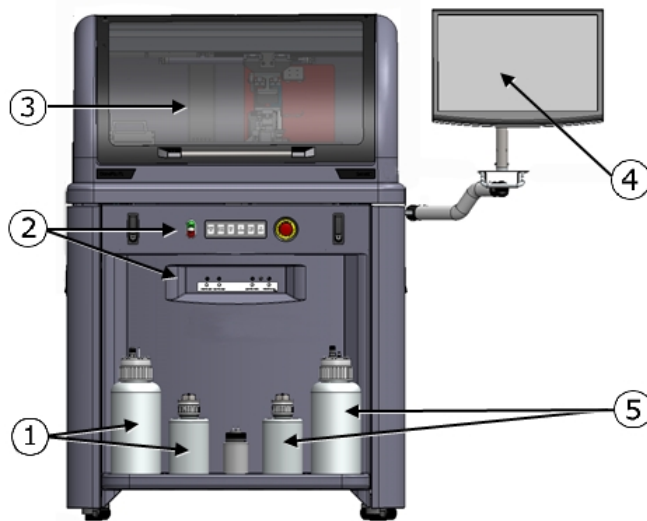


The ClonePix 2 Mammalian Colony Picker is installed at your site by Molecular Devices.

Before you operate the instrument or do maintenance operations, make sure that you are familiar with the safety information in this guide. See [Safety Information on page 6](#).

The ClonePix 2 system provides a controlled environment for sample protection conducive to mammalian cell picking. Protection within the enclosed process bed is provided by a rear mounted HEPA filtered air supply and a front mounted exhaust fan, which creates a horizontal laminar flow of sterile air across the picking area of the bed when the front door is closed.

The process area of the instrument contains source and destination stacker lanes, and wash baths that clean the pins before and during a process. See [Process Area on page 22](#).



**ClonePix 2 Mammalian Colony Picker Overview**

| Item | Description          |
|------|----------------------|
| 1    | Feed Bottles         |
| 2    | Front Control Panel  |
| 3    | Enclosure Front Door |
| 4    | Computer Monitor     |
| 5    | Waste Bottles        |

For information on cleaning the instrument, see [Cleaning the Instrument on page 85](#).

For more information on regular instrument maintenance, see [Maintenance and Troubleshooting on page 83](#).

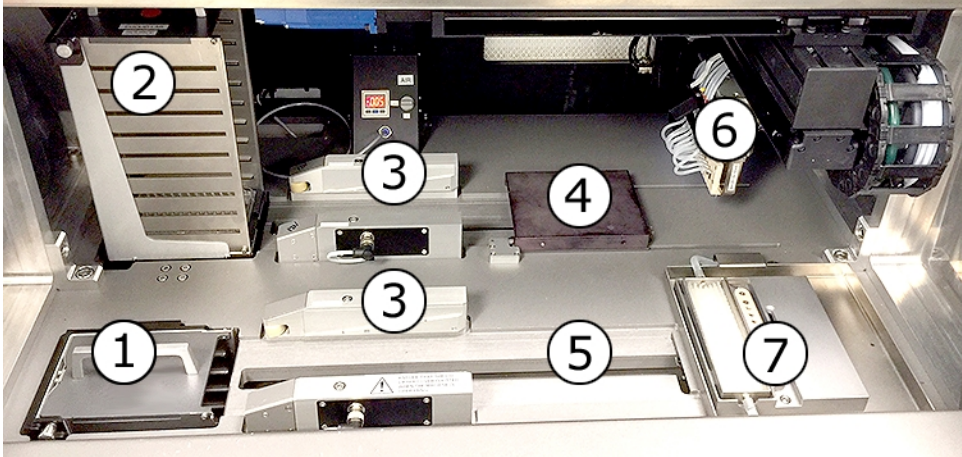
For technical specifications, see [Technical Specifications on page 99](#).

## Process Area



**ClonePix 2 Mammalian Colony Picker Process Area**

| Item | Description                    |
|------|--------------------------------|
| 1    | Destination Stacker            |
| 2    | Source Stacker                 |
| 3    | Microplate Lid Lifters         |
| 4    | Enclosure Front Door           |
| 5    | Picking Bed and Source Imaging |
| 6    | Destination Dispensing Bed     |
| 7    | Picking Head                   |
| 8    | Picking Pins Wash Station      |

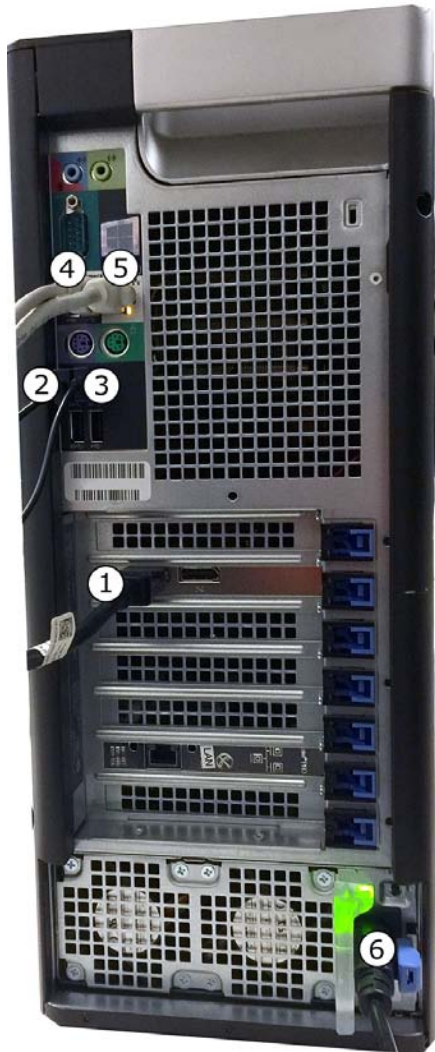


ClonePix 2 Mammalian Colony Picker Process Area Close-Up

| Item | Description                    |
|------|--------------------------------|
| 1    | Destination Stacker            |
| 2    | Source Stacker                 |
| 3    | Microplate Lid Lifters         |
| 4    | Picking Bed and Source Imaging |
| 5    | Destination Dispensing Bed     |
| 6    | Picking Head                   |
| 7    | Picking Pins Wash Station      |

## Computer Connections

The ClonePix 2 system uses the following computer cable connections:



ClonePix 2 Mammalian Colony Picker Computer Connections

| Item | Description |
|------|-------------|
| 1    | Monitor     |
| 2    | Mouse       |
| 3    | Keyboard    |
| 4    | Camera      |
| 5    | Ethernet    |
| 6    | Power cable |

For computer details, see [Computer System Requirements on page 15](#).



Front Panel Controls and Display

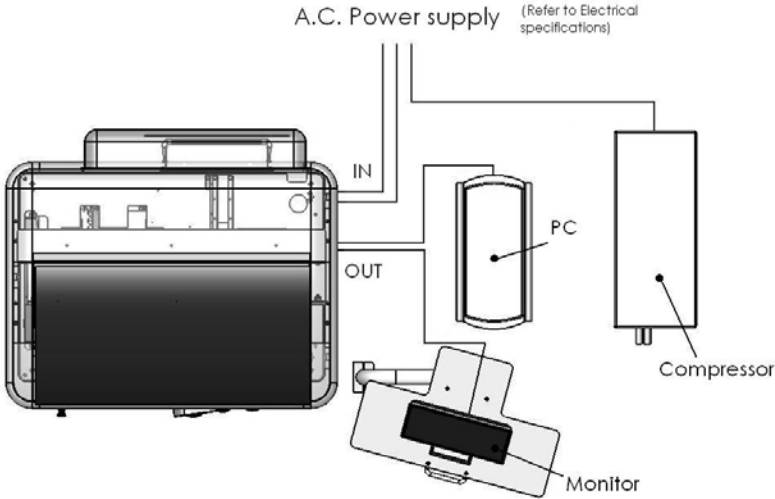


ClonePix 2 Mammalian Colony Picker Front Panel Controls and Display

| Item | Description                 |
|------|-----------------------------|
| 1    | Feed Tube Connection Ports  |
| 2    | Stop Button                 |
| 3    | Start Button                |
| 4    | Activity Status Lights      |
| 5    | Emergency Stop Button       |
| 6    | Waste Tube Connection Ports |

Electrical Connections

The ClonePix 2 system uses the following power cable connections:



## Bottles Connections

The ClonePix 2 system uses the following bottle-contained fluid carrying tubing system:



The Picking Feed bottle connection carries sterile water to the Wash Station and back to the Picking Waste bottle. The Ethanol Feed bottle connection carries ethanol to the Ethanol Wash Station and back to the Ethanol Waste bottle. The outlets on the instrument are in the same order as the bottles.

Make sure that the bottles are marked so that the tubes can be re-connected correctly when you need to refill and empty.

## Compressed Air Supply

Compressed air is required for the picking movement of the picking pins and the movement of the imaging cover. Laboratories with built-in compressed air systems can connect directly to the filter-regulator inlet on the side of the ClonePix 2 instrument. Alternatively, an optional oil-free compressed air unit can be connected that draws air from the local environment and delivers it to the instrument through the filter-regulator inlet and filters it to 0.02 microns using a 3-stage filtration process.



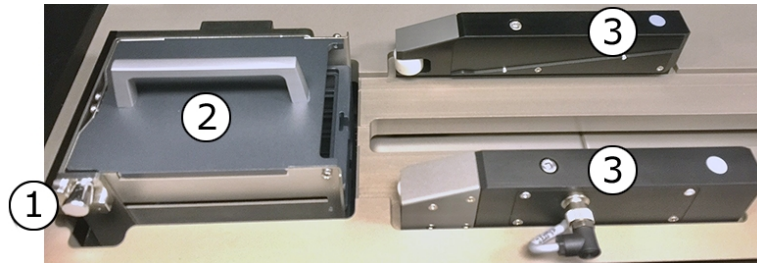
### ClonePix 2 Mammalian Colony Picker Compressed Air

| Item | Description                                     |
|------|---|
| 1    | Air inlet from house air or optional compressor |
| 2    | Air outlet to instrument                        |

For the inlet air-pressure requirements, see [Technical Specifications on page 99](#).

## Microplate Stackers

The stacker lanes and cassettes enable the ClonePix 2 Mammalian Colony Picker to run processes with multiple microplates. The stacker can be used in every process. The stacker cassettes are tall metal cases designed to hold up to 10 microplates and lid pairs. The source stacker cassette and destination stacker cassette must always be used together in a lane when processing microplates and lid pairs. The source stacker cassette is used for automating the loading of the microplates onto the lane so that they can be automatically moved into the instrument for colony inoculation. The microplates are automatically collected from the instrument in the destination stacker cassette after colony inoculation. The lid is removed as the microplate is pushed through the lid lift mechanism. The lid is replaced as the microplate returns through to the stacker cassette.



**ClonePix 2 Mammalian Colony Picker Microplate Stacker**

| Item | Description           |
|------|-----------------------|
| 1    | Lock pin              |
| 2    | Stacker cassette      |
| 3    | Microplate Lid Lifter |

### Removing a Stacker Cassette

To remove the stacker cassettes:

- Pull the lock pin outwards and lift the cassette out of the stacker hole.



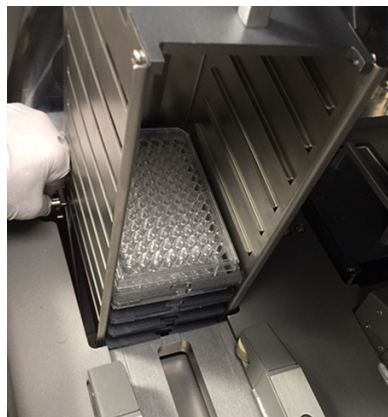
## Loading and Installing a Stacker Cassette

To load and install a stacker cassette:

1. With the cassette removed from the instrument, load the microplates with lids on and the A1 well at the front of the slot.



2. Pull the lock pin outwards and lower the cassette into the stacker hole until the lock pin clicks into place.



## Filters

The ClonePix 2 Mammalian Colony Picker supports the following filter options:

| Description   | Excitation Wavelength (nm) | Filter Position | Emission Wavelength (nm) | Filter Position |
|---|----------------------------|-----------------|--------------------------|-----------------|
| CFP AF430   | 440 (Royal Blue)           | 4               | 505 (Cyan)               | 4               |
| CloneDetect FITC* Fluorescein<br>GFP AF488          | 470 (Blue)                 | 3               | 535 (Green)              | 3               |
| YFP   | 500 (Cyan)                 | 2               | 550 (Yellow)             | 2               |
| CloneDetect 549* R-phycoerythrin<br>Rhodamine AF546 | 530 (Green)                | 1               | 590 (Red)                | 1               |
| CloneDetect 649* Cy5 FMAT Blue<br>AF647             | 622 (Red)                  | 6               | 700 (Far Red)            | 6               |



**Note:** White light imaging uses excitation filter position 7 and emission filter position 5.

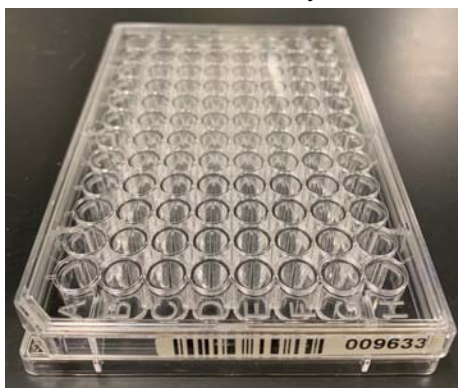
## Using Barcodes

If you use barcodes, the following barcode parameters are required:

- Use a legible barcode of the following types: 1D (linear) barcodes with code 11, 39, 93 and 128.



- Do not use special characters, such as hyphens, in the barcode. Special characters can cause missed reads and other errors downstream.
- Place the barcode centrally on one of the short sides of the plate.



**Plate With Barcode**

For troubleshooting, see [Troubleshooting Barcodes on page 94](#).

## Camera Alignment

For pin-to-spot accurate picking, you must correctly calibrate and align the camera. The ClonePix 2 Mammalian Colony Picker camera calibration and alignment is done by Molecular Devices and verified during on-site installation.

Knocking the actuator can impact picking accuracy. If you need to re-calibrate the camera, contact Molecular Devices Technical Support [support.moleculardevices.com/](https://support.moleculardevices.com/).

## Safety Features

The ClonePix 2 system uses the following safety features:

### Enclosure Front Door

The enclosure front door uses an electromagnetic switch that prevents the instrument from running when the door is open. The door is made of acrylic, which traps UV light in the instrument.

The electromagnetic switch protects the following:

- It prevents the motors from running, which protects against any physical damage.
- It disables the UV light, which protects against the risk of damage from UV radiation.



**WARNING!** Never tamper with the electromagnetic door switch.

---

### UV Light

Inside the process area is a UV germicidal lamp with timer. It is a 30W linear discharge lamp with a defined output at 253.7 nm to make the lamp an effective source of germicidal radiation.



**WARNING!** Dispose according to your local regulations.

---

### Emergency Stop Button

Pressing the Emergency Stop button immediately stops the instrument and cuts power from the drives. The button must be pulled out before the machine can be restarted.



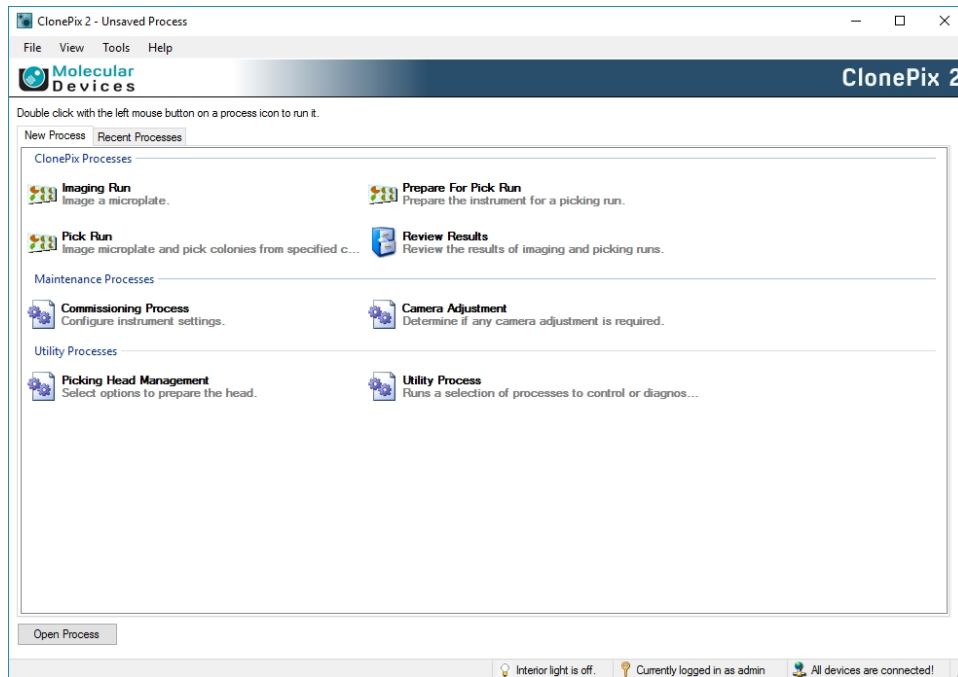
**CAUTION!** If the computer is powered through the instrument, power to the computer is also be cut and your current data is not saved.

---

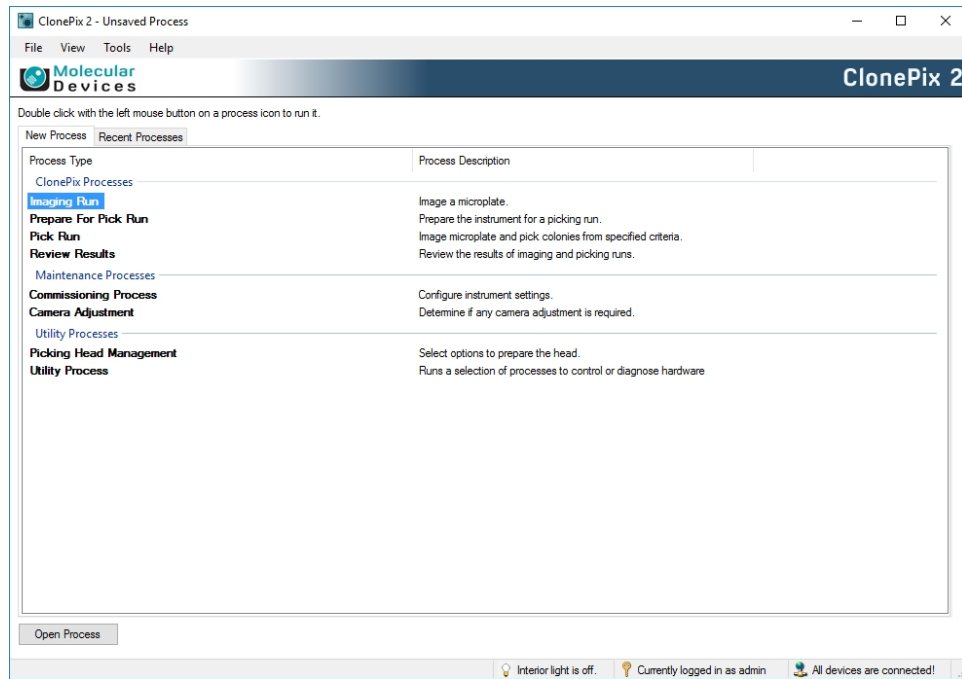
The ClonePix 2 software controls the ClonePix2 instrument and is based around the concept of the instrument running processes. A Process is a standard program for carrying out a task. Using the default set of processes, you can modify and save them. Processes are managed from the Main Navigation screen, that displays when the software starts.

## Main Navigation Screen

The Main Navigation screen displays when the ClonePix 2 software starts. From this screen, you select to run processes.



Use the right-click option menu to view the process options as icons or as a list.



The Navigation screen includes the following functionality:

- **New Process Tab** displays the available processes, which are grouped into the following categories:
  - **ClonePix 2 Processes**
  - **Maintenance Processes**
  - **Utility Processes**
- **Recent Processes Tab** displays a list of recently opened processes. This is the same list of processes available from the **File > Recent Processes** menu option.
- **Interior Light** in the status bar at the bottom of the screen, the interior light can be activated or deactivated at any time by double-clicking the interior light icon.
- **Open Process** button displays the Open dialog where you can select and open a saved process file. Can also be opened using the **File > Open Process** menu option.
- **Login Credentials** in the status bar at the bottom of the screen, displays which user-login is connected, if any.
- **Connection Status** appears in the bottom right corner of the Navigation screen and displays the status of the Ethernet connection between the computer and the ClonePix 2 system.



## Menu Options

The available options in the menus at the top of the Main Navigation screen change depending on the screen view and selections.

### File Menu

- **Open Process** allows you to open a previously saved process.
- **Save Process** saves the current process.
- **Save Process As** allows you to save the current process with a different name or location.
- **Close Process** closes the current process and returns you to the Process Selection Window.
- **Save As Template** saves the current process as a template that can be used for creating new processes.
- **Recent Processes** allows you to select and open a recent process.
- **Switch User** can be used in a multiple user environment to log off the current user so that a different user can log on.
- **Exit** closes the Navigation window and shuts down the software.

### View Menu

- **Properties** displays the properties of the process that is being set before starting the process.
- **Progress** displays the progress of a running process.
- **Administrate Properties** lets you change the Properties view and default values.

### Tools Menu

- **Configuration** displays the Edit Configuration dialog.



**CAUTION!** We recommend that only trained personnel configure these settings.

---

- **Prepare Error Report** starts the Error Reporting Wizard to create a data file containing the configuration and recent log files to help troubleshoot your problem.

### Help

- **About** displays the version numbers of the software modules.
- **Online Support** displays the support web page. This option requires an active internet connection.

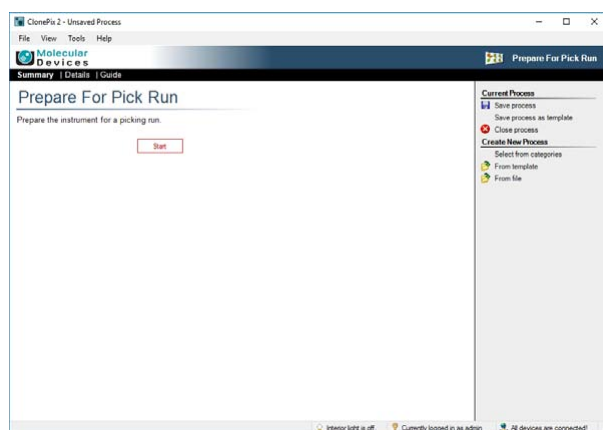


### Prepare for Pick Run

Run the Prepare for Pick Run process before you run a Pick Run process to ensure that the system is ready for picking. Any section of the following process can be bypassed by clicking **Next**.

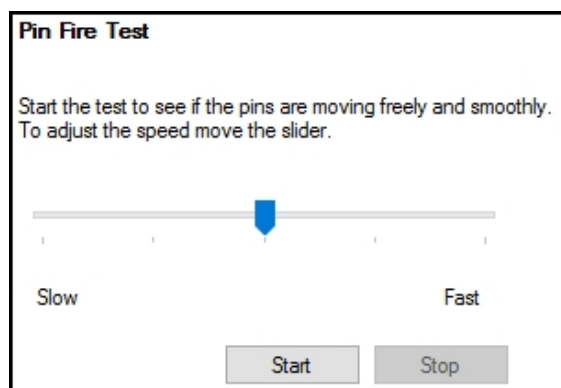
**\* Tip:** This process is typically run only once a day, before the first pick run. When multiple pick runs are planned, Prepare for Pick Run is not required before each picking run.

Click **Start** to begin the process.



### Pin Fire Test

Run the Pin Fire Test to confirm that the pins are firing correctly.



To use Pin Fire Test:

1. Set the speed to fast and click **Start**.
2. Observe that the pins move freely.
3. Click **Stop** to end the test.
4. Click **Next** to move to [Alignment on page 36](#).



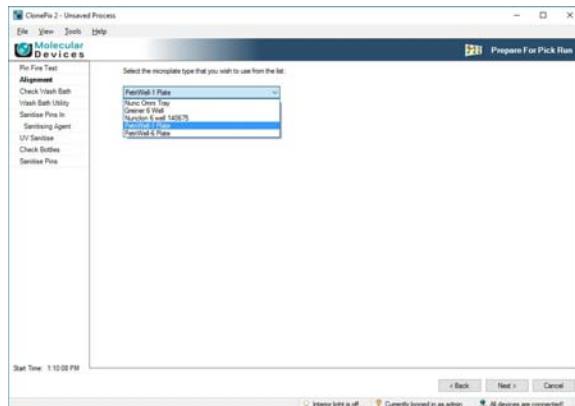
**CAUTION!** If any pin is not firing and retracting freely, do not continue. Run the Pin Fire Test continually for up to one hour. If this does not free up the pin action, do not continue. Remove the Picking Head and contact Molecular Devices Technical Support to arrange for reconditioning. The Picking Head cleaning procedures described in the General Maintenance section of the Robot User Guide are for sterilization purposes only and are unlikely to resolve any pin movement issue.

## Alignment

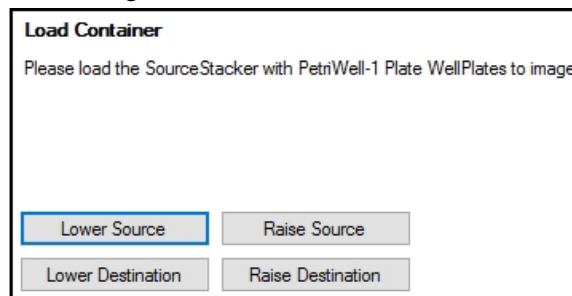
Run the Alignment step to align the source plate and the primary pin to the camera.

To use Alignment:

1. Select the source plate type from the drop-down list and click **Next**.



2. If needed, to raise the source stacker click **Raise Source**. If the destination stacker is obstructing access to the source stacker, to lower it, click **Lower Destination**.



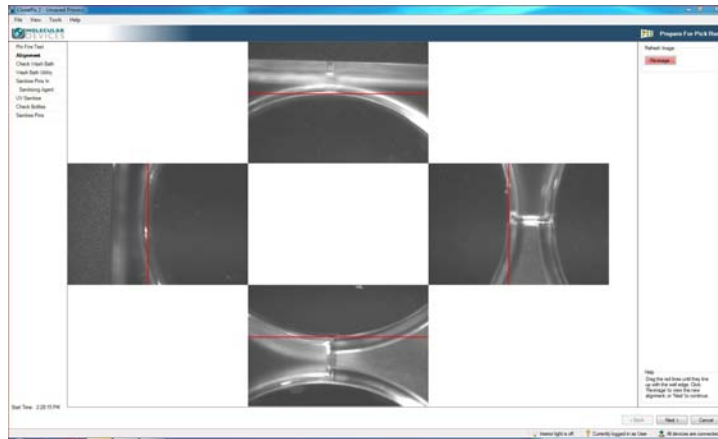
3. Load an empty source plate into the source stacker cassette.



**Tip:** The empty source plate must be cell-free and media-free. Verify that your source plate is loaded into the appropriate color-coded plate holder before inserting into the cassette.

4. Click **Next** to initiate imaging of the edges of well A1.

- If alignment is needed, drag the red lines to line up the well and then click the **Re-image** button.



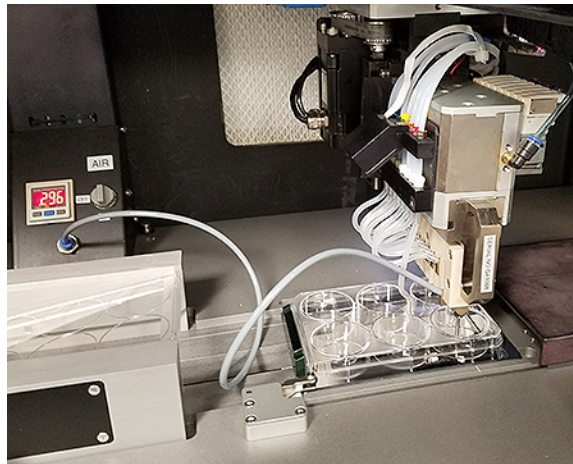
- Click **Next** to move to the alignment of the primary pin Z height.

**\* Tip:** In normal usage, the Z height of the primary pin should not require adjustment. Moving onto the next stage without modification is recommended.

If pin height adjustment is required:

- Click **Retract Pin**.
- Open the instrument door and remove the picking tube from the pin 1.
- Attach the tube from the air pressure gauge to the pin 1.

The air pressure gauge displays a value in red.



- Turn the pressure gauge to the **ON** position by twisting the gray knob clockwise so that the handle is vertical.

- Click **Go To Source Stacker**.



**Note:** The datum point can be set only with the pin extended.

Visit Position

Stop Short Above

Set Position

Click the buttons to jog the specified distance.

Vertical

Lateral

Move Size:

Vertical Position

The datum can only be changed with the pin extended. Lower the pin until it just touches the bottom of the well. You can use the Air Pressure Meter to find the bottom of the well. Make sure you attach this at the beginning of the datuming process because the process of attaching it can change you datum point ever so slightly. Press 'Set' to set the tip of the extended picking pin as the new Datum point. When this is complete press the 'Next' button.

- Use the Vertical arrow buttons to lower the pin in small increments until the value on the pressure gauge turns from red to green.



**Tip:** To change the move increment, select a new value from the Move Size list. When the pressure gauge value starts to increase, reduce the move increment. Record the previous Z datum value for guidance and comparison.

- Before you set the datum, adjust the Z position so that the pin is just touching the well bottom using the smallest move increment 0.01 mm, and then back off by 0.01 mm.
- To set this datum point, click **Set**, and then when prompted, to save the datum, click **Yes**.
- Before you continue, click **Retract Pin**, turn off the pressure gauge, remove the air pressure tube from pin 1, and then reconnect the picking tube.
- To move to alignment of the primary pin X & Y coordinates, click **Next**.

To align the pin correctly:

1. Click **Go To Source Stacker**.

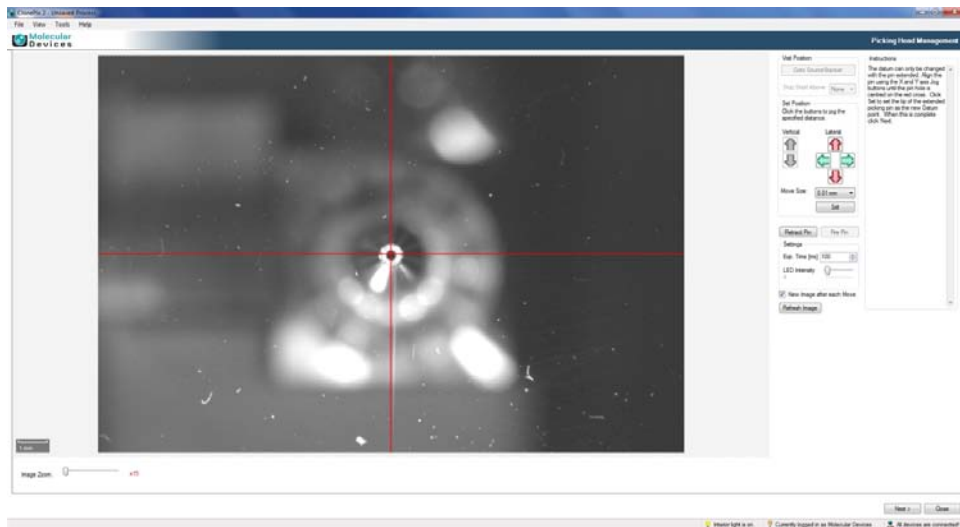


**Note:** There is no need to choose Stop Short Above as the pin already stops short by 0.5 mm.

2. Use the **Lateral jog** buttons to move the pin until the pin hole is centered on the red cross.



**Tip:** Use the **Image Zoom** slider to zoom in on the image for accurate alignment.



3. When the pin is in the required position, click **Set**.



**Note:** If you do not click Set, the new datum point will not be fixed.

4. If the image is not bright enough, increase the **Exposure Time** or **LED Intensity**.
5. Click **Next** to set up for sterilization of the fluid system.

## Check Wash Bath

To check the wash bath:

1. Check to see if there is any 70% ethanol in the wash bath.
  - If the system has not been used, the bath might be empty, fill it in the next stage of the process.
  - If the system has been used already, the bath might be filled enough.
2. Click **Next**.

## Wash Bath Utility

The Wash Bath Utility step allows the Wash Bath to be filled and the Auto Replenish to be set as required.


To use the Wash Bath Utility:

1. Click **Start** to fill the Wash Bath.
2. Click **Stop** when the Wash Bath is filled.
3. Select **Auto Replenish Wash Bath** to turn on the Auto Replenish function. This keeps the wash bath filled while the software is running. Change the default timings, if necessary.
4. Click **Next**.

## Sanitize Pins in Sterilant

The Sanitize Pins in sterilant step sanitizes the pump system between the sterile wash bottle and the picking pins.


---

 **Tip:** Sanitize the pins at the start of each day. We recommend using Spor-Klenz® Ready-To-Use Cold Sterilant.

---

As an alternative, you can use 70% ethanol to sanitize the pins and feed system. See [Alternative: Sanitize Pins in 70% Ethanol on page 41](#).

---

 **CAUTION!** Do not leave sterilant in the system for more than an hour.

---


To use Sanitize Pins with sterilant:

1. Connect a wash bottle containing the freshly prepared sterilant to the wash supply line. For preparation instructions, see the documentation that came with the sterilant.
2. To pump the sterilant through the entire system, click **Start**. This takes several minutes. The sterilant must be left for 10 minutes. This can be done at the same time as UV sterilization. See [Ultra Violet Sanitize on page 41](#).
3. Click **Next**.



### Alternative: Sanitize Pins in 70% Ethanol

This alternative sanitize step uses 70% ethanol to sanitize the pump system between the sterile wash bottle and the picking pins.

 **Tip:** Sanitize the pins at the start of each day.



**CAUTION!** Do not leave ethanol in the system for more than an hour.

**Sanitise Pins**

Purge the pump system with sanitising agent

Sanitise Options

|   |    |
|---|----|
| Number of cycles for the head to purge                      | 30 |
| Number of cycles the pins will be scrubbed in the Wash Bath | 5  |
| Number of seconds the Dryer will be on for                  | 10 |

To use Sanitize Pins with 70% ethanol:

1. Prepare a solution (about 1 liter) of 70% ethanol (70% ethanol: 30% deionized water) in a ClonePix feed bottle.
2. Connect the bottle containing 70% ethanol. Run purge to fill the system. Leave the solution in the system to disinfect for 20 minutes.
3. Replace the ethanol feed bottle with a feed bottle containing high-quality sterile water, connect the sterile water bottle and purge to rinse out the ethanol.
4. Click **Next**.

### Ultra Violet Sanitize

The Ultra Violet Sanitize process activates the interior germicidal lamp.

While the UV sanitise is in progress you can check the fluid level in the waste bottles and make sure the feed bottles are full.

Ultra Violet Sanitise

Duration of Ultra Violet exposure:  seconds

To use Ultra Violet Sanitize:

1. Set the required duration. 10 minutes (600 seconds) is recommended.
2. Click **Begin**. The lamp turns off after the set time. If the protective door is opened, the lamp switches off until the door is closed. After the door closes, the duration time resets to 10 minutes and can be activated again.
3. Click **Next**.

## Check Bottles

The feed system must be purged with sterile water.

To check bottles:

1. Disconnect the fluid supply line from the bottle that contains sanitizing solution (Sterilant or 70% ethanol) and connect to an autoclaved bottle of high-quality sterile water.
2. Empty the waste bottle.
3. To set up purge conditions, click **Next**.



**Tip:** The connectors seal automatically when disconnected. As a sanitation precaution, liberally spray both ends with 70% ethanol before you connect the wash bottle.

## Sanitize Pins

**Sanitize Pins**

Purge the pump system with sanitising agent

Sanitise Options

|   |    |
|---|----|
| Number of cycles for the head to purge                      | 30 |
| Number of cycles the pins will be scrubbed in the Wash Bath | 5  |
| Number of seconds the Dryer will be on for                  | 10 |

Start

To start the Sanitize Pins purge:

1. Click **Start** to pump sterile water through the entire system. This takes several minutes.

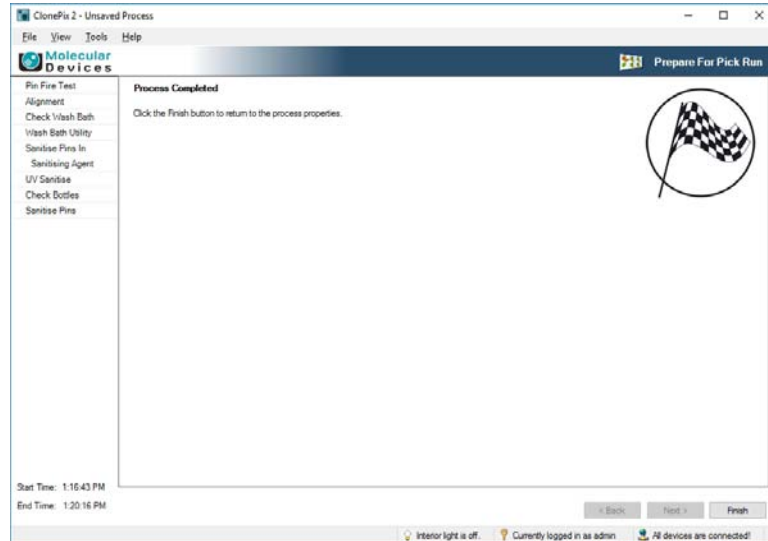


**CAUTION!** Never bypass sanitize steps. If the system does not need to be sterilized, it is still essential to remove any air bubbles that have appeared if the system has been idle for any time, especially overnight. Air in the line can seriously affect picking success rates.

2. Click **Next**.

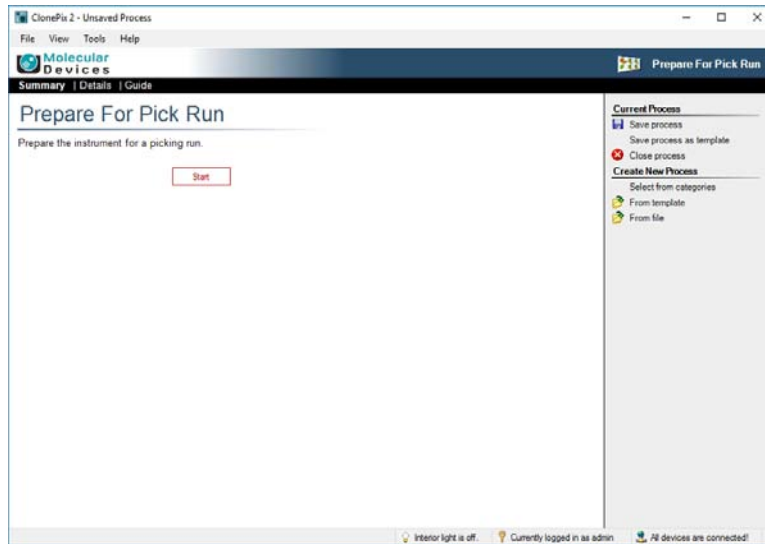
The system parks the head safely, returns the source plate, and closes the illumination cover.

## Finish



To finish:

1. Click **Finish** to return to the Prepare for Pick Run process Start page.



2. Click **Close Process** to return to the Main Navigation screen. See [Main Navigation Screen](#) on page 31.

## Pick Run Process

The Pick Run process includes default parameters required for a full imaging and picking run, which involves image acquisition, colony detection, group selection, and automated picking. Use the Pick Run process to create multiple user-defined Pick Run or Batch Pick Run processes that let different users reconfigure the robot to their own preferences.



**Note:** When a Pick Run finishes, the data is accessible in Review Results, but the data is locked and cannot be modified. See [Review Results Process on page 75](#).

ClonePix 2 - Unsaved Process

File View Tools Help

Molecular Devices

Summary | Details | Guide

### Pick Run

Image microplate and pick colonies from specified criteria.

#### Imaging Settings

Run Annotation: PetriWell-1 Plate  
 Source Microplates: Read Barcode: True  
 Barcode Options: Auto-assign barcode in case of failure: False  
 Batch plates: False  
 Source Plate Options: Prompt for more plates when cassette is empty  
 Acquisition Options: Undefined  
 Review Colony Selection: Prime: Undefined  
 First Cycle

#### Picking Settings

Destination Microplates: PetriWell-96 Plate  
 Destination Wells: All wells will be available  
 Barcode Options: Read Barcode: True  
 Auto-assign barcode in case of failure: False  
 Deposit Options: Fill Destination Plates.  
 Dest. Plate Options: Prompt for more plates when cassette is empty  
 Pick Number Options: Organise By: Plate  
 Limit the number of colonies picked: false  
 Pin Options: Aspirate Volume: 5  
 Dispense Volume: 7  
 Adherent Colonies: True  
 Suspension Colonies: False  
 Picking Height Adjustment: -0.1  
 Audit Options: Allow Target and Aspirate images to be acquired but not saved.  
 Dispersal Options: Use Dispersal: False  
 Dispersal Volume: 20  
 Dispersal Cycles: 10

#### Sanitise Pin Options

Sanitise Pin Options: Purge Cycles: 3  
 Bath Cycles: 3  
 Dry Time: 10

Start

Current Process

- Save process
- Save process as template
- Close process

Create New Process

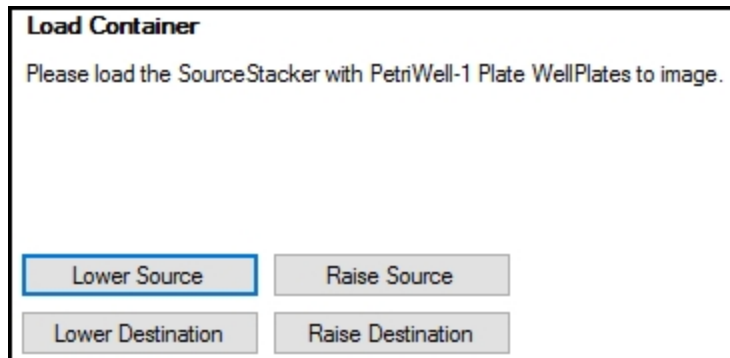
- Select from categories
- From template
- From file

Interior light is off. Currently logged in as admin All devices are connected!

## Start Pick Run

To start Pick Run:

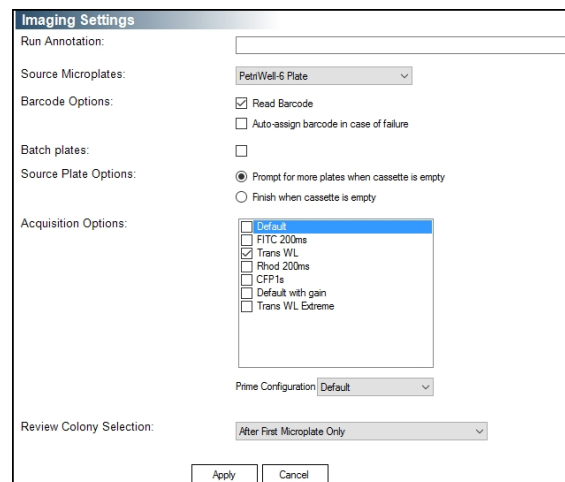
1. Before you continue, make setting changes. On the Pick Run > Start screen:
  - [Change Imaging Settings on page 45](#)
  - [Change Picking Settings on page 47](#)
  - [Change Sanitize Pin Options on page 49](#)
2. Click **Start** to start the run process .
3. When the Load Container dialog displays, load the source plates into the source stacker.



4. If needed, raise or lower the source stacker.
5. Click **Next**.

## Change Imaging Settings

To edit Imaging Settings, click the **Imaging Settings** title, or alternatively, the **Details** tab or **Guide** tab.



To change the following:

- **Run Annotation** - Enter a name to identify the run.
- **Source Microplates** - Select the type of source well plates.

- **Barcode Options** - Select one of the following options:
  - **Read Barcode** - Activates the source plate barcode reader. Must be activated for a Batch Pick Run.



**Tip:** Using barcodes is recommended. Barcodes must be placed on the front short side of the microplate (A1 end). Also see [Using Barcodes on page 29](#).

---

- **Auto Assign Barcode In Case Of Failure** - Automatically generates a unique plate identifier if the barcode fails to read.  
Clear this checkbox to have the system request an identifier to be entered manually.



**Note:** Compatible barcodes for the system are 39, 93, and 128. Also see [Using Barcodes on page 29](#).

---

- **Batch Plates** - Selects the **Batch Pick Run** option where all plates in the cassette are imaged and statistics are generated as an aggregate for all plates. The colony selection for the whole series of plates can be reviewed and adjusted before picking.
- **Source Plate Options** - Select one of the following options:
  - **Prompt For More Plates When Cassette Is Empty** - Prompts you to add more plates to the cassette after all the plates in the cassette are imaged.
  - **Finish When Cassette Is Empty** - Ends the imaging process after all the plates in the cassette are imaged.
- **Acquisition Options** - Select from the list of acquisition options that have already been created. You can modify this in Preview.
- **Prime Configuration** - Use these settings for detection of all colonies. Generally, this is the white light acquisition option, although a fluorescent option designed for colony detection could be used, for example, viability assay.
- **Review Colony Selection** - Select one of the following options:
  - **After First Microplate Only** - Displays the Results screen after only the first microplate is imaged to allow you to adjust the Groups before you commit to picking. These settings apply to any other plates in the cassette.
  - **After Every Microplate** - Displays the Results screen after each microplate is imaged to allow you to adjust the Groups before you commit to picking.
  - **Batch – Review All** - Must be selected when you use the Batch Pick Run. All source plates are imaged to allow the groups to be adjusted for the whole series of plates before you commit to picking.

## Change Picking Settings

To edit Picking Settings, click the **Picking Settings** title, or alternatively, the **Details** tab or **Guide** tab.

### Picking Settings

Destination Microplates: PetriWell-96 Plate ▼

Destination Wells:

|   |   |   |   |   |   |   |   |   |   |    |    |    |
|---|---|---|---|---|---|---|---|---|---|----|----|----|
|   | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
| A |   |   |   |   |   |   |   |   |   |    |    |    |
| B |   |   |   |   |   |   |   |   |   |    |    |    |
| C |   |   |   |   |   |   |   |   |   |    |    |    |
| D |   |   |   |   |   |   |   |   |   |    |    |    |
| E |   |   |   |   |   |   |   |   |   |    |    |    |
| F |   |   |   |   |   |   |   |   |   |    |    |    |
| G |   |   |   |   |   |   |   |   |   |    |    |    |
| H |   |   |   |   |   |   |   |   |   |    |    |    |

Barcode Options:  Read Barcode  
 Auto-assign barcode in case of failure

Deposit Options:  Match Destination plate to Source plate

Dest. Plate Options:  Prompt for more plates when cassette is empty  
 Finish when cassette is empty

Pick Number Options:  Collate by Well  
 Limit Colonies  
0 Number of Colonies From Each Plate

Pin Options:  Adherent  
 Suspension 500 Semi-Solid Media Delay (ms)  
Pick Height Adjustment -0.10 Below Well Bottom  
Aspirate Volume (µl) 5  
Dispense Volume (µl) 7

Audit Options:  Save Target and Aspirate Images

Dispersal Options:  Use Dispersal  
Dispersal Cycles 10  
Dispersal Volume 20

You can change the following:

- **Destination Microplates** - Select the type of destination well plates.
- **Destination Wells** - Specify the deposit wells to use. Right-click to select wells and left-click to deselect wells. All destination plates are filled using this template.

- **Barcode Options** - Select one of the following options:
  - **Read Barcode** - Activates the source plate barcode reader. Must be activated for a Batch Pick Run.
  - **Auto Assign Barcode In Case Of Failure** - Automatically generates a unique plate identifier if the barcode fails to read.  
Clear this checkbox to have the system requests an identifier to be entered manually.



**Note:** Compatible barcodes for the system are 39, 93, and 128.

- **Deposit Options** - Select to have the destination plate matched to the source plate such that the completion of picking from each source plate prompts the return of the current destination plate. When not selected (default), all destination plate wells are sequentially filled.
- **Destination Plate Options** - Select one of the following options:
  - **Prompt For More Plates When Cassette Is Empty** - Prompts you to add more plates to the cassette after all the plates in the cassette are imaged.
  - **Finish when cassette is empty** - Ends the imaging process after all the plates in the cassette are imaged.
- **Pick Number Options** - Colonies can be picked by source plate or by source well. By default picks as if all wells of the current plate contain the same sample and are organized by source plate. Select one of the following options:
  - **Collate By Well** - When selected the system picks all specified colonies from well A1 first and then proceeds to the next well. Select this when, for example, wells of a multi-well source plate contain different samples.
  - **Limit Colonies** - Limits the number of colonies picked from each plate or well.
- **Pin Options** - Select one of the following options:
  - **Adherent or Suspension** - Select the **Adherent** option to collect colony monolayers and select the **Suspension** option to collect colonies from semi-solid medium. By default, there is a 500ms delay for Suspension picking to permit the colonies to be fully collected from the semi-solid medium.
  - **Pick Height Adjustment** - The pick height is automatically adjusted for optimal Adherent or Suspension picking (by default these are 0.1 mm below and 0.4 mm above well bottom, respectively).
  - **Aspirate Volume** - General recommendation is no less than 10  $\mu$ L.
  - **Dispense Volume** - General recommendation is no less than 12  $\mu$ L.
- **Audit Options** - Saves the target and aspirate images before and after picking a colony. If you select this option, all images are saved. If you do not select this option, you can toggle the images on and off while picking but the images are not saved.



**Note:** This option slows picking dramatically.


- **Dispersal Options** - Selecting dispersal separates the cells of a picked colony by aspirating and dispensing in the destination plate. If dispersal is required, you can set the following parameters:
  - **Dispersal Cycles** - 3 to 6 for CHO cells and 6 to 10 for hybridomas. Small colonies in chemically defined media can be left undispersed.
  - **Dispersal Volume** - Maximum volume that can be dispersed is 20  $\mu$ L.

Click **Apply** To save the settings.



## Change Sanitize Pin Options

To edit Sanitize Pin Options, click the **Sanitize Pin Options** title, or alternatively, the **Details** tab or **Guide** tab.

 **Tip:** The default settings are appropriate for most situations.

### Sanitize Pin Options

Sanitize Pin Options:

|   |    |
|---|----|
| Number of cycles for the head to purge                      | 3  |
| Number of cycles the pins will be scrubbed in the Wash Bath | 3  |
| Number of seconds the Dryer will be on for                  | 10 |

Apply
Cancel

To edit Sanitize Pin Options:

1. Specify the number of purge cycles.
2. Specify the number of cycles the picking pins are scrubbed in the ethanol bath.
3. Specify the number of seconds the halogen dryer is on.

If you create new settings, click **Save Process**.

## Preview

The Preview screen starts the image acquisition and colony detection set up.

The first tab that displays in Preview mode is the Images tab. This initially displays as a gray until you select an image in the well map to acquire the images.

The screen is divided into two panels.

- The panel on the left displays the acquired images, with composite, white light and fluorescent channel thumbnails. To display these thumbnails in the central image frame, left-click. You can only capture and view one image area of the well at a time, represented by a light blue rectangle in the Well Map.
- The panel on the right, displays the Well Map and several tabs at the bottom that enable you to adjust imaging settings. Tabs include:
  - **Acquisition** - See [Image Acquisition Settings on page 50](#)
  - **Detection** - See [Image Detection Settings on page 51](#).
  - **Non-cellular** - See [Non-Cellular Tab on page 65](#).
  - **Groups** - See [Groups Tab on page 66](#).
  - **Statistics** - See [Statistics Tab on page 69](#).

**\* Tip:** The most important settings to set up correctly are the image Acquisitions and the Detection of the colonies. Altering these settings later in the process can take a long time.

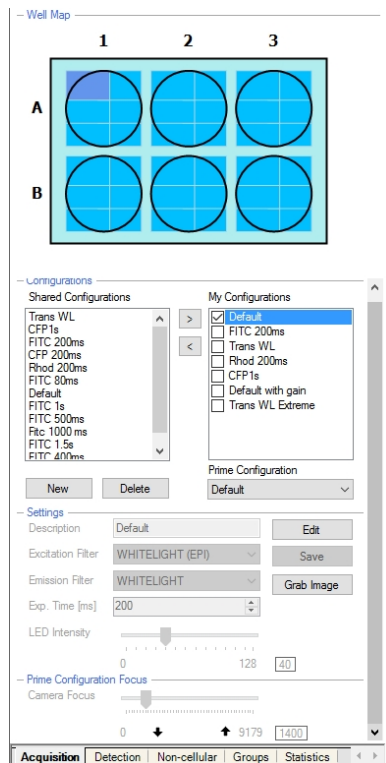
### Image Acquisition Settings

Use the **Preview > Image > Acquisition** tab to define settings for image capture. Make sure that the images are bright but not over exposed to the white or fluorescent light.

Image overexposure displays as red pixels on the image. Where these are seen, reduce the exposure time.

To select Configurations for image acquisition, double-click each configuration name.

To preview the chosen images in the current image area, click **Grab Image**.

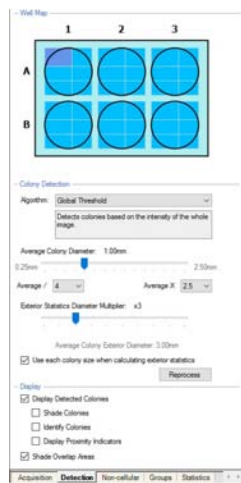


Settings include the following:

- **Well Map** - A map of the plate and the selected well displays above each of the settings tabs. Click any of the wells imaged (displayed in light blue) to display the images for that well.
- **Configurations** - Defined acquisition options are shown in the My Configurations box after they are created. To share personal configurations with other users, click on the appropriate option in My Configurations to highlight the text in blue and then click < to transfer to the Shared Configurations list.  
**Prime Configuration** - Use for colony detection. Generally, this is the white light acquisition option or a fluorescent option designed for colony detection, for example, viability assay such as Live Detect.
- **Settings** - Use to set up and modify acquisition options.
  - To modify a configuration, click to highlight the text in blue and then click **Edit**.
  - Select appropriate settings and then click **Save**.
  - To create a new configuration click **New**, then click the new configuration to highlight the text in blue, and then click **Edit**.
  - Select settings and then click **Save**.
  - For white light imaging, use the **Trans** (transillumination) option. For fluorescent imaging, use the maximum LED intensity.
- **Prime Configuration Focus** - This is the focus value that the camera uses when acquiring the images for each configuration.

### Image Detection Settings

Use the **Preview > Image > Detection** tab to define optical settings for detecting colonies within the prime configuration (white light). It is important at this stage to make sure the all the desired colonies are detected. Detected colonies are displayed with a colored circle around the perimeter of the colonies (the color of the circle corresponds to the groups it is automatically assigned to by default).



Settings include the following:

- **Well Map** - A map of the plate and the well you select displays above each of the settings tabs. Click any of the wells imaged (displayed in light blue) to display the images for that well.
- **Colony Detection** - Select one of the following Algorithms:
  - **Global Threshold** - Detection based on single background threshold intensity of the whole image.
  - **Local Threshold** - Detection based on background intensity around the vicinity of the colony.
  - **Flatten Detection** - Locates features by first removing background variations.
  - **Edge Detect** - Locates colonies by detecting the transition at the edge of each colony. This is better suited to detect adherent colonies.

**Average Colony Diameter** - Use the scale bar on the image to set the most appropriate average colony diameter. The software searches for colonies within a ten fold range of colony diameters therefore excluding any large or small features that are not colonies. Minimum and maximum diameters are shown for the diameter you specify. It is possible to extend this range by changing the upper and lower limits displayed as drop-down lists. The default values are the average colony diameter divided by 4 for the lower limit and the average colony diameter multiplied by 2.5 for the upper limit.

**Exterior Statistics Diameter Multiplier** - Move the cursor over any colony on the image to display a red circle around the colony. This defines the area that is used to calculate Exterior statistics. Use the Diameter Multiplier slider to set the required diameter to the red circle as multiples of the Average Colony Diameter.

**Use Each Colony Size When Calculating Exterior Statistics** - Select to use the actual size of each colony, rather than the Average Colony Diameter so that the larger colonies have a larger exterior red circle.

To view the effect of any change, click **Reprocess**.

- **Display** - Use the following options to display various aspects of the images:
  - **Display Detect Colonies** - Clear this checkbox to hide colony detection from view to aid visualization of the underlying raw image.



**Note:** Be sure to re-select the display, otherwise subsequent processing is hidden.

---

- **Shade Colonies** - Select this option to fill the colony outline with the corresponding color for better visualization.
- **Identify Colonies** - Superimposes a feature ID for each colony.
- **Display Proximity Indicators** - Select to display the nearest adjacent colony for every detected colony in the image.



**Note:** This is not the same as the Group Proximity (colony exclusion).

---

- **Shade Exclusion Zone** - Selected to display the exclusion areas on the image as hatched red lines. Clear to remove the exclusion areas around the image.

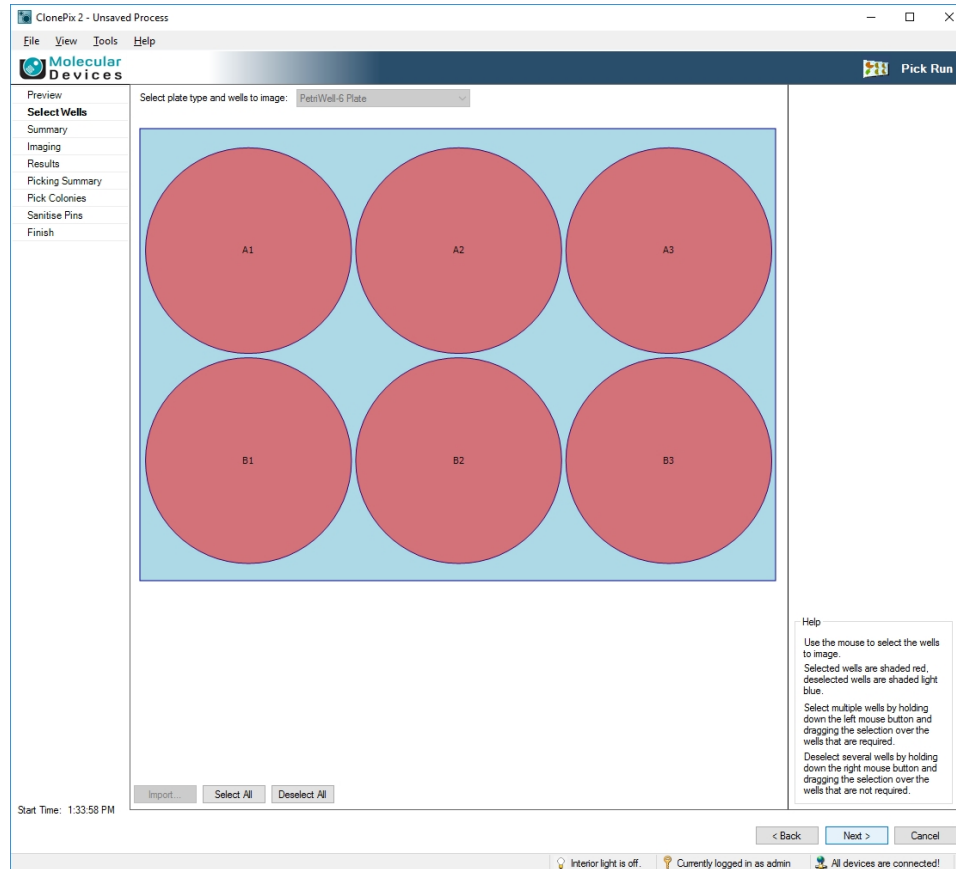
## Select Wells

Use the Select Wells screen to choose which wells to image.

To deselect the wells (showing them as white), right-click on a well .

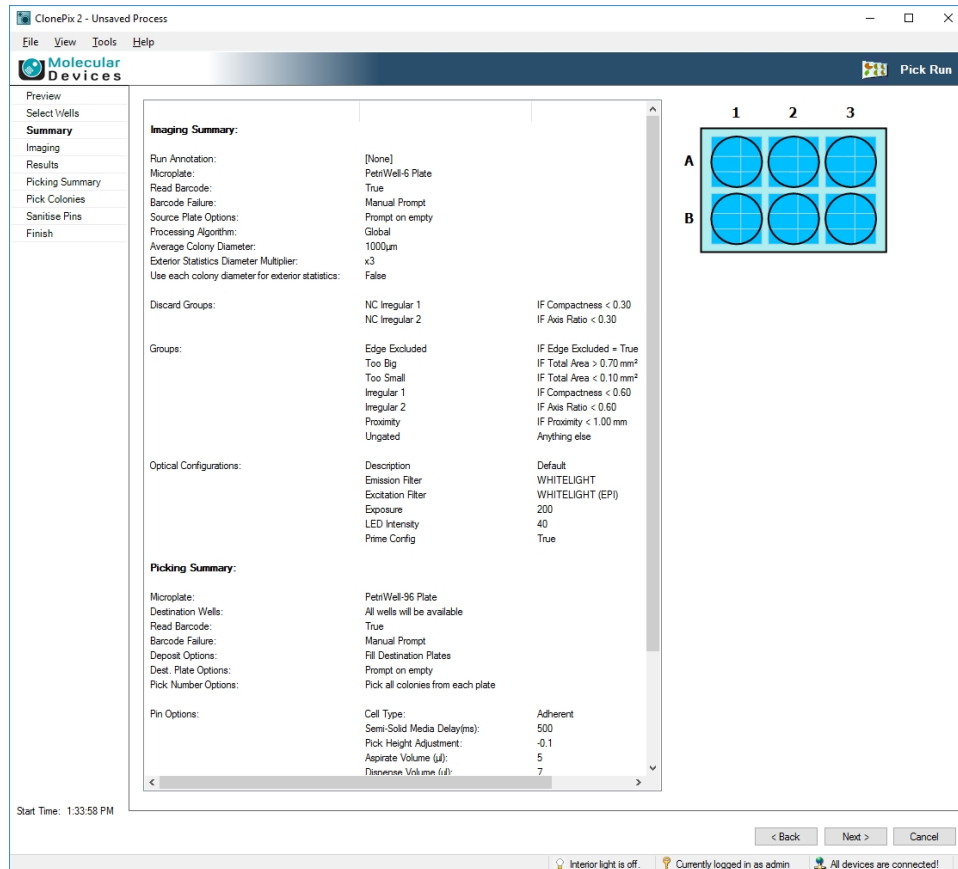
To select the well (showing them as pink), left-click on a well .

Alternatively, click **Select All** or **Deselect All** to add or removed wells to be imaged.



## Summary

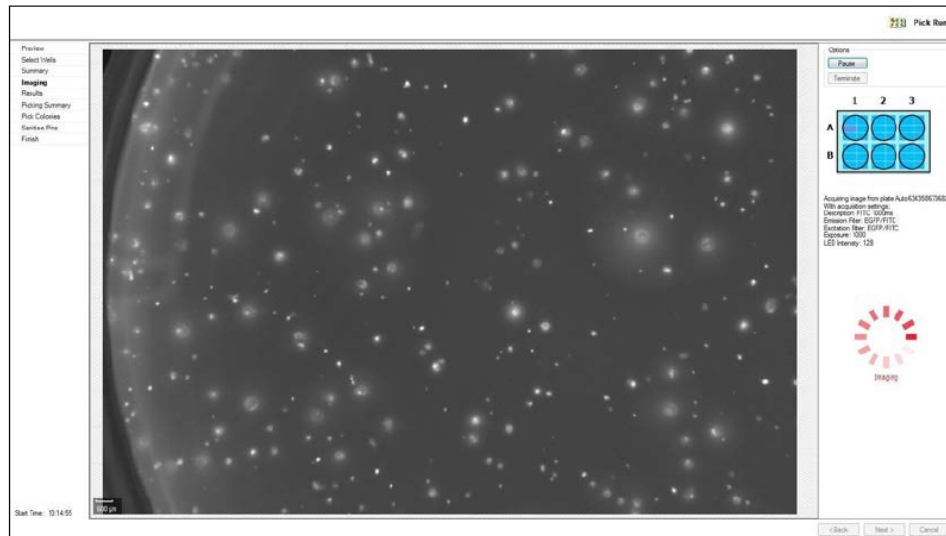
Use the Summary screen to confirm that all settings are correct before you proceed to imaging. Both imaging and picking summaries appear. A well map to the right of the summary identifies the wells that are imaged.



## Imaging

The Imaging screen displays the images that are being captured as imaging progresses. The selected wells image in sequence with the acquisition options defined. The images are then processed together for colony detection according to the defined settings.

A well map displays the wells being imaged (displayed in light blue) and the current image being taken (displayed in dark blue). A small summary appears below the well map showing the acquisition details of the image being captured.



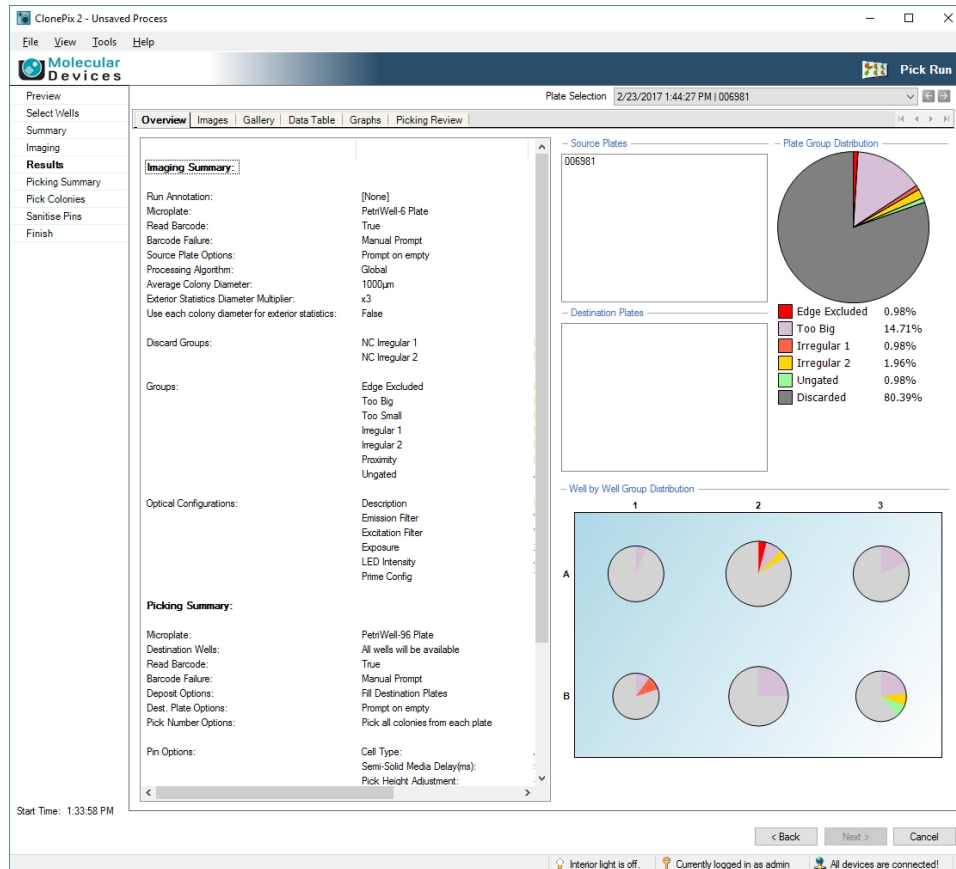
To cancel during imaging, click **Pause**, and then click **Terminate**.

## Results

The Results screen displays when imaging ends to show you the data in several tabs.

This screen has a similar format to the Preview screen except:

- You can use the Well Map to navigate all image areas.
- The Acquisition tab displays only the configurations used during imaging. The only settings you can alter are selection of Prime Configuration.



Use the tabs on the Results screen to discard debris or unwanted colonies and then create selection groups for picking in the next stage.

To aid with viewing data between tabs, you can select a population by clicking and dragging on the Scatter / Ranking graphs, or by clicking while holding the **CTRL** key on the Image tab or Statistics tab. Colonies you select are visualized as follows:

- In the Scatter graph or Ranking plot on the Graphs tab, the data points are colored dark blue.
- Within images, the colonies are highlighted with a blue/white circle.
- On the Data Table tab, the data points are highlighted with a blue line above and below the row.

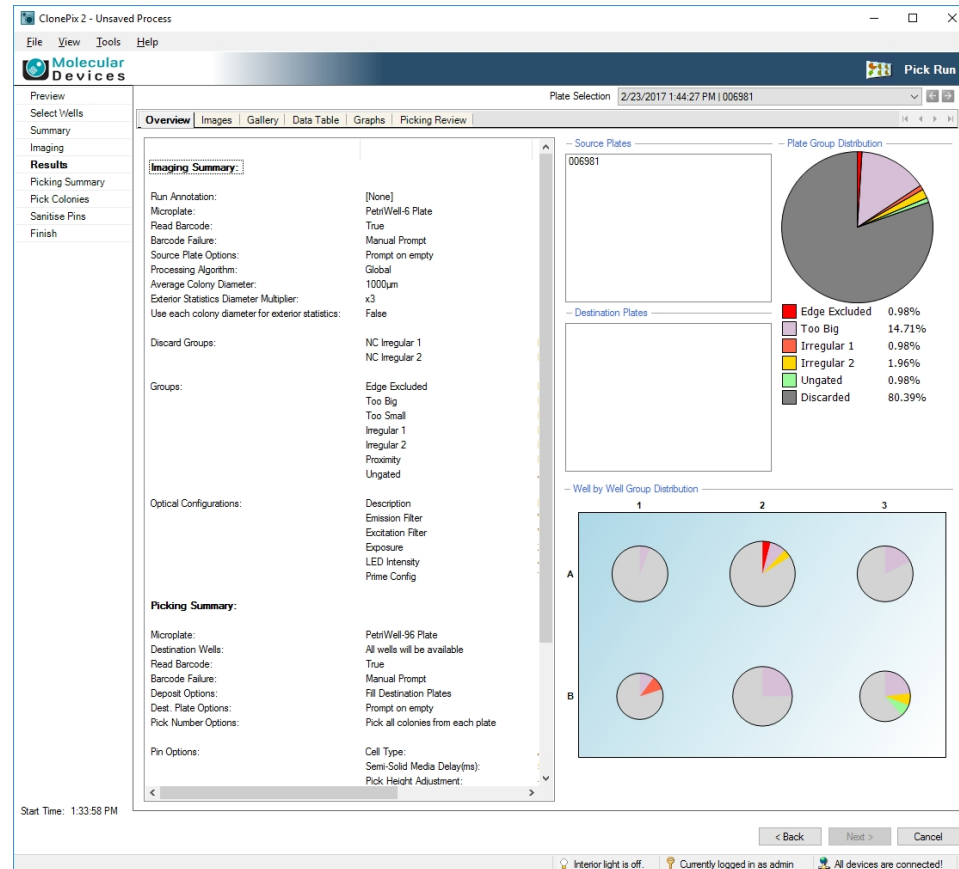


**Note:** If you select Review Colony Selection > After First Microplate Only, the Results screen is bypassed for subsequent plates. See [Change Imaging Settings on page 45](#).



## Overview Tab

The Overview tab displays a summary of the imaging and picking settings.

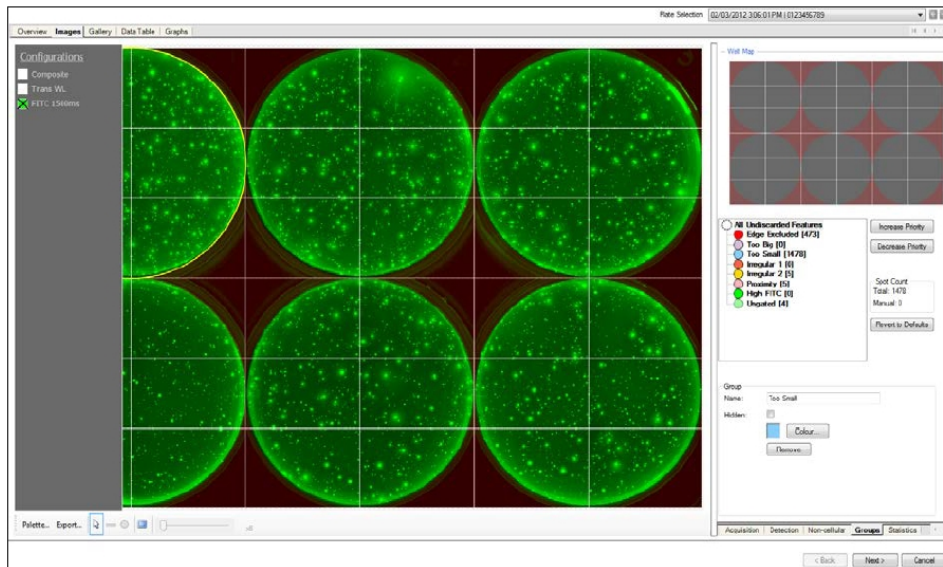


A summary of the plates imaged displays on the right, including the following:

- **Source Plates** - The source plates imaged are listed here. When clicked, highlighted in blue, the well map appears below in the Well by Well Group Distribution section.
- **Destination Plates** - The destination plates used in the picking process appear here. At this stage in the process, no plates are visible because no colonies have been picked.
- **Well by Well Group Distribution** - A selected plate appears here. Each imaged well appears as a pie chart with each slice of the pie representing a group. The pie size is also proportional to the number of colonies detected in the well.
- **Plate Group Distribution** - The group distribution of all the wells in all the plates imaged appears in a pie chart. The percentage of each group appear below the pie chart. When you move the cursor over this key, the percentages display as the actual number of colonies in each group. Any discard groups created on the Non-Cellular tab (see [Non-Cellular Tab on page 65](#)) appear together as the Discarded proportion of the pie chart in dark gray.

## Images Tab

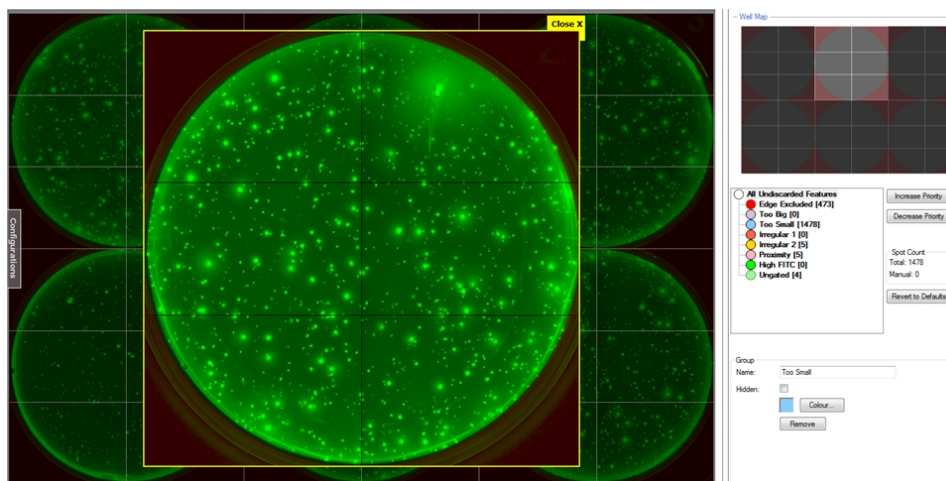
The Images tab displays every image frame from all the plates imaged.



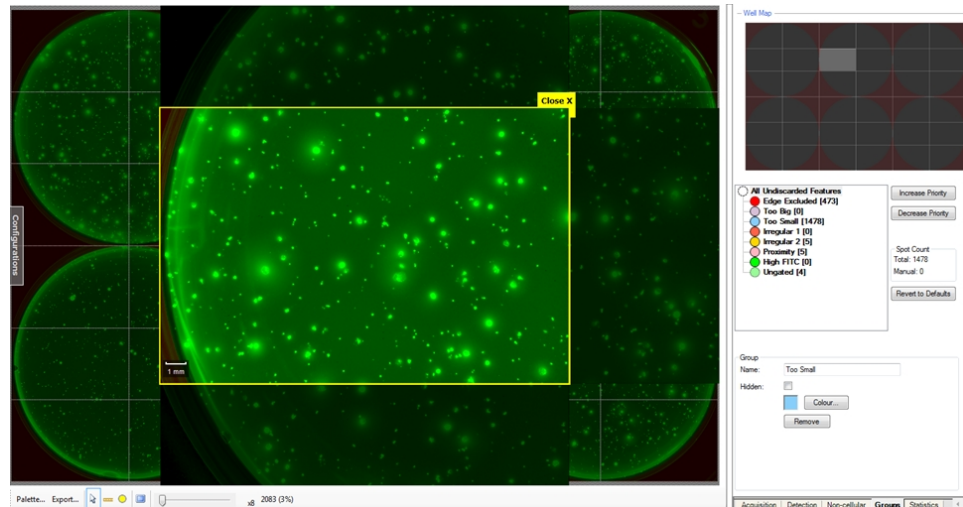
To view a specific image:

1. In the **Well Maps**, hover the cursor over the frame and click it.
2. Select the **Configurations** tab, to open the Configurations menu.
3. In the **Configurations** menu list of the different filters used to image the plate, select the checkbox for each filter name to view the image.

Hover over the whole plate image to cause either wells (in the case of plates containing multiple wells) or frames to be highlighted with a surrounding yellow box. When you select a well, a larger image of the well displays on top of the whole plate image, the Well Map updates to reflect this selection.



Hover over the whole well image to cause frames to be highlighted with a surrounding yellow box. When you select a frame, a larger image of the frame displays on top of the whole plate image and the Well Map updates to reflect this selection. This is Single Image mode.



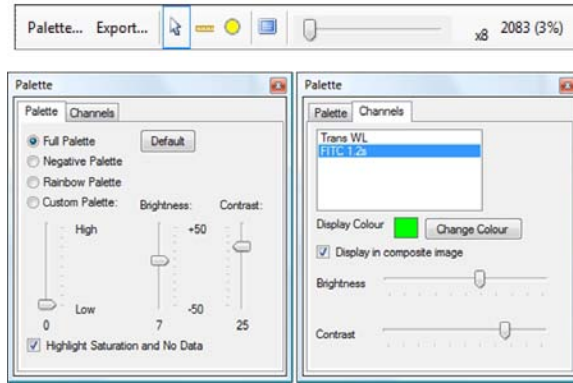
While in Single Image mode, frames next to the selected image also appear but display darker. Click these frames to slide the adjacent frame into view and load the newly selected image. The Well Map is also functional in Single Image mode and enables quick cross plate navigation.

Below the Well Map are the Acquisition, Detection, Non-cellular, Group, and Statistics tabs. Select the **Acquisition** tab or **Detection** tab to display settings that you selected during the Preview. See [Preview on page 49](#). You can change these settings but when you use Reprocess, all the captured images are reprocessed, which can take some time, especially if you use batch mode and ten plates were imaged. See [Non-Cellular Tab on page 65](#), [Groups Tab on page 66](#), and [Statistics Tab on page 69](#).

Controls include the following:

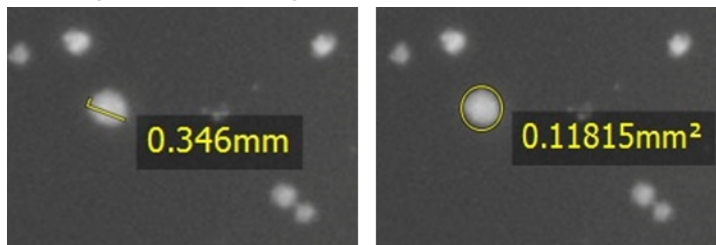
- **Composite** - Fluorescent (non-prime) images can be superimposed onto prime images in the Composite image view. It is possible to change the brightness and so on, of the image by selecting the checkbox and then using Palette.
- **Trans WL** - The destination plates used in the picking process appear here. At this stage in the process, no plates are visible because no colonies have been picked.
- **FITC 1.5 s** - Fluorescent checkboxes display below the white light checkbox. If there are several fluorescent channels used (for example, FITC, RHO, CY5) they are all displayed individually below the white light checkbox. It is possible to change the brightness and so on, of the image by selecting the checkbox and then using the Palette.

- Palette** - Palette button allows the images on-screen to be enhanced or modified. There is a palette available for each filter (Composite, White Light, and Fluorescent channels). To change any of the settings, select the checkbox and click palette. There are various palettes to choose from for improving the brightness and contrast of the image. The Highlight Saturation and No Data option (selected as default) displays the red (saturated) and blue (No Data) pixels on the image. This can be removed by clearing this checkbox.



You can also apply colors to the images on the Channels tab by selecting the channel of interest and giving it a color using the **Change Color** option. You can display the fluorescent channels in the composite image by selecting **Display** in composite image box. You can alter the brightness and contrast of the composite image using the sliders in this section of the palette.

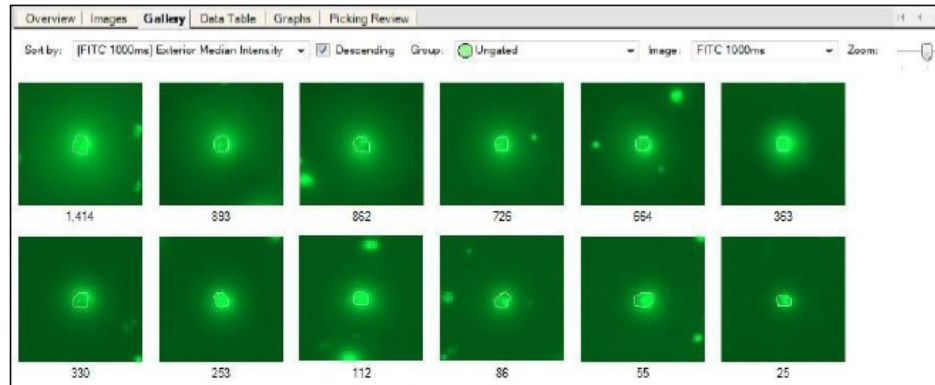
- Export** - Export Image permits the current image to be exported. The exported image can be generated with or without detection (user defined) and is shown in the color specified in the acquisition option. The image can be exported as a .bmp, .jpg, or a .png file.
- Select, Ruler, and Area**
  - Select** - The default tool, when the cursor is over a colony in Single Image mode, the properties of that colony appear on the Statistics tab. This tool also allows panning and zooming around the image.



- Ruler** - Measures diameter of a colony. Click **Ruler**, and then drag the cursor across a colony. Click **Select** after measurement.
  - Area** - Measures the area of a colony. Click **Area**, and then click the center of the colony and drag out to the perimeter. Click **Select** after measurement.
- Full Screen** - Enlarges the image to fill the full screen. Press **ESC** to return to regular view.
- Zoom Control** - Digitally zoom images. The magnification displays to the right of the Image Zoom bar, which turns red when the actual image size is exceeded. Beyond this point, the image becomes pixelated. The scale bar in the bottom left corner of the image pane adjusts with the Zoom Control.

## Gallery Tab

Use the Gallery tab to define the visualization order of the colonies to be picked. It is possible to order them according to a fluorescent or morphological statistic, group, and the acquisition option used (White light, FITC).



The colonies can display using the following settings:

- **Sort By** - Sorts the colonies according to the statistic chosen from the drop-down list. This can be a fluorescent statistic or a morphological statistic.
- **Descending** - Select the checkbox to arrange the colonies in descending order from left to right. Clear the checkbox to arrange the colonies in ascending order from left to right.
- **Group** - Enables all the features or any of the groups to be displayed.
- **Image** - Enables the white light or fluorescent image of the colonies to be displayed.
- **Zoom** - Magnifies into and out of the image by moving the slider from left to right.
- **Viewing Colonies** - Shows the number of colonies being viewed.

### Data Table Tab

The Data Table tab displays the statistics for colonies detected in the image. Each line represents a detected colony colored according to its Group. Fluorescent image parameters denoted by the prefix [acquisition option] display to the left and Prime (white light) image parameters display to the right. By default, parameters are listed in alphabetical order. Click a column header to sort data by the other parameters. Click a second time to invert the order.

The data table is interactive. Double-click on a row of colony data to link to the image of the colony (identified with a blue circle). Alternatively, right-click a row to display a list of link options.

The right-click menu also allows you to manually link colonies to a Group. Any modifications to automatic grouping are logged in the statistics column Manual Group.



Select a row to display the statistics for that colony on the Statistics tab to the right of the data table.

Statistics are ordered for easy viewing:

- **Colony Images** - When you select a colony in the data table, the corresponding image displays to the left of the table. you can change the size of the colonies in the thumbnails by using the Zoom slider at the bottom.
- **Select Statistics** - There are various statistics that appear as default, however you can display only the preferred statistics. To do this, open the select statistics window and select the desired statistics. After you click **Close**, only the chosen statistics display and the drop-down lists on the Graphs tab.



**Note:** Deselecting statistics hides only the data from view and can be reselected at any time.

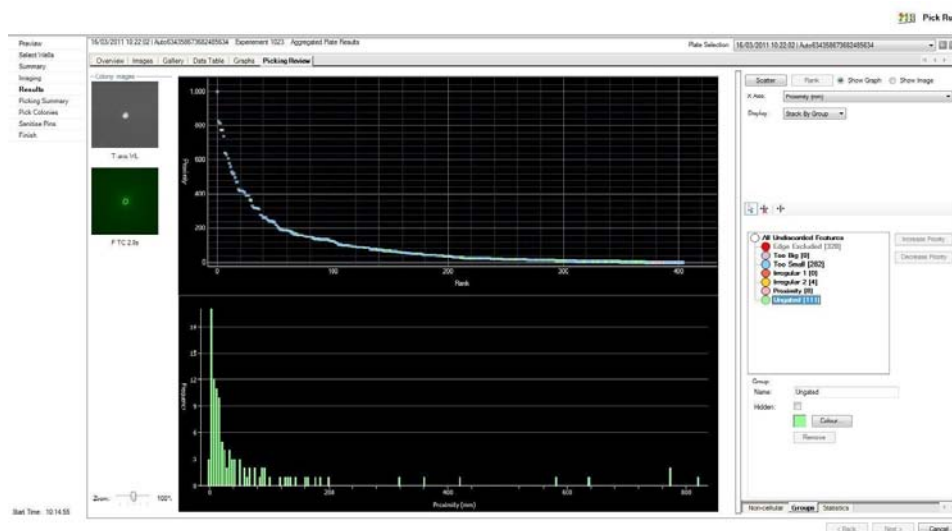
- **Export Statistics** - Select this to open a window to export data. Only statistics for picked colonies or all detected colonies can be exported. Data can be exported as:
  - CSV format for opening as a single sheet in Excel and other statistics software.
  - Excel-XML format for opening in multiple sheets in Excel and other statistics software.
  - XML format for compatibility with LIMS and other bespoke software. This is not functional in Preview.
- **Generate Report** - Selected to run a wizard to generate an XML-based summary of a run.

You can export the following information into the Report for each plate processed:

- **Imaging Settings** - All settings applied during set up.
- **Picking Settings** - All settings applied during set up. Only available for a completed Pick Run.
- **Statistics** - All the statistics you select in the data table are entered into the report. It is recommended that these are excluded from the report as the data table is extensive and this makes the report large. It is recommended that statistics are exported through Export Statistics.
- **Scatter Plot** - The last Scatter plot to be created on the Graphs tab is entered into the report.
- **Rank Plot** - The last Rank plot and Histogram created on the Graphs tab is entered into the report.
- **Pick Summary** - All the details from the Picking Review tab are entered into the report. As well as giving the report a name, you can enter the following information into the wizard; Title, Cell Type, Project Details, Plating Efficiency, Incubation Time, Summary.

## Graphs Tab

Use the Graphs tab to create groups. Each point that displays within the graphs is a detected colony. The color of each point indicates the group that the colony has been assigned to. Hover over each point on the graphs to display the corresponding colony in the image thumbnails to the left of the graphs.

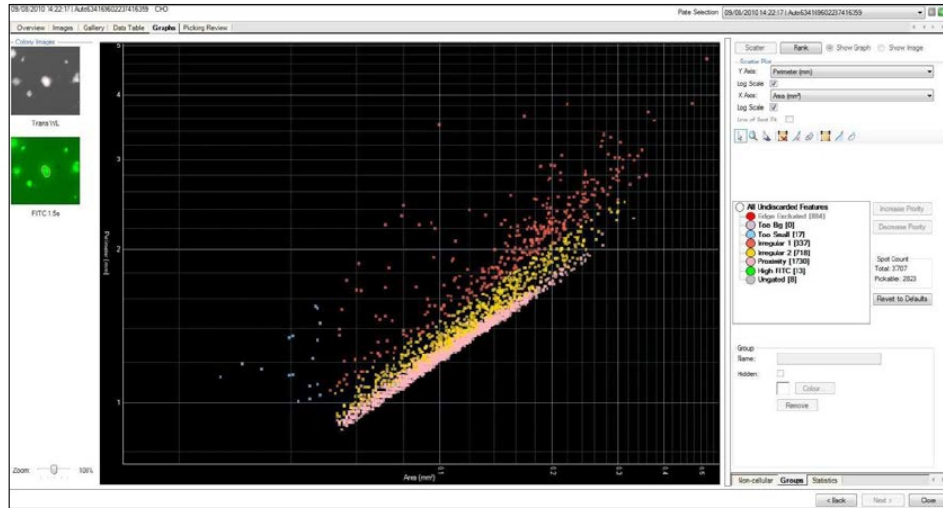


Graphing options include the following:

- **Scatter** - When you select a colony in the data table, the corresponding image displays to the left of the table. Use the Zoom slider at the bottom to change the size of the colonies in the thumbnails.

- **Rank** - There are various options when looking at results and creating groups in the ranking plot. When you select Rank, the Rank Plot displays at the top of the screen and a Histogram displays beneath the rank plot.

**Rank Plot** - Display either the rank plot or a well image. When you select Show Graph, the ranking plot displays above the histogram. The Rank Plot and Histogram statistics are linked. When you select a statistic in one graph, it is also displays on the other. You can select to Log Scale the Y Axis in the Rank Plot and also zoom into an area on the graph. When you select Show Image, the image of the well you select displays above the histogram. As you create or edit groups, the color change of the colonies update on the image.

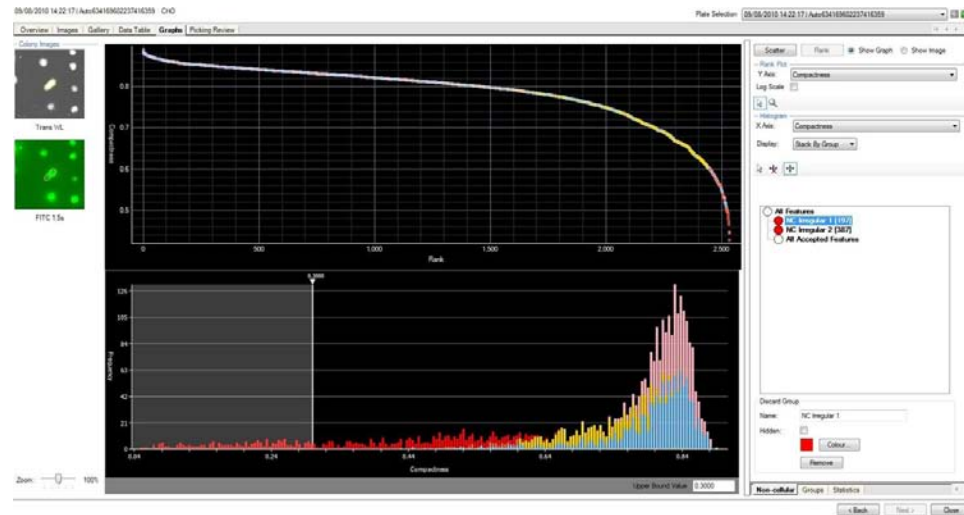


- Rank Plot** is interactive. Double-click a data point to link to the image of the colony (identified with a blue circle). Alternatively, right-click on a data point to displays a list of link options. Right-click on the background to copy the current graphic image.
- **Histogram** - Create groups using this histogram by selecting a statistic from the drop-down list and drawing a line gate where required. You can also change group default values here. **Display** allows you to visualize the grouping histogram by Stack By Group or by Totals Only.



## Non-Cellular Tab

Use the Non-Cellular panel to remove any kind of debris from the data set, for example, hairs, small scratches, and small bubbles, that are not colonies before you decide which colonies to pick. By removing debris from the data set before you create groups, the debris is not considered as being part of the data. This means that real colonies are not put into the Proximity exclusion group if they are too close to debris. There are two discard groups by default; NC Irregular 1 and NC Irregular 2. These groups exclude debris based on compactness and axis ratio respectively. You can delete these groups if not required. Non-Cellular discard groups can be created by selecting a statistic from the drop-down list and drawing the discard line (icon denoted by a red cross) on the histogram.



You can create several of these Non-Cellular discard groups and these groups are listed on the Non-Cellular tab. By default all discard groups listed on the Non-Cellular tab are colored red but you can change this if needed. Discard groups can also be hidden by selecting the group and selecting the **Hidden** checkbox.

## Groups Tab

Use the Groups panel to visualize detected colonies and their corresponding groups and to create new groups for picking. You can optimize and setup groups at the Preview stage, but it is best to create groups after imaging. The grouping system is designed to generate exclusion groups for example, Too Small, Irregular, Proximity, and inclusion groups to pick specific populations, for example, High FITC, High Rhodamine. You can pick multiple groups in the same picking run and you choose which groups to pick later on the [Picking Review Tab on page 69](#).

Image thumbnails display to the left of the graphs where you can hover over the points on the graph to visualize the colonies.

There is the option to change between **Rank** and **Scatter** graph for plotting statistics and creating new groups.

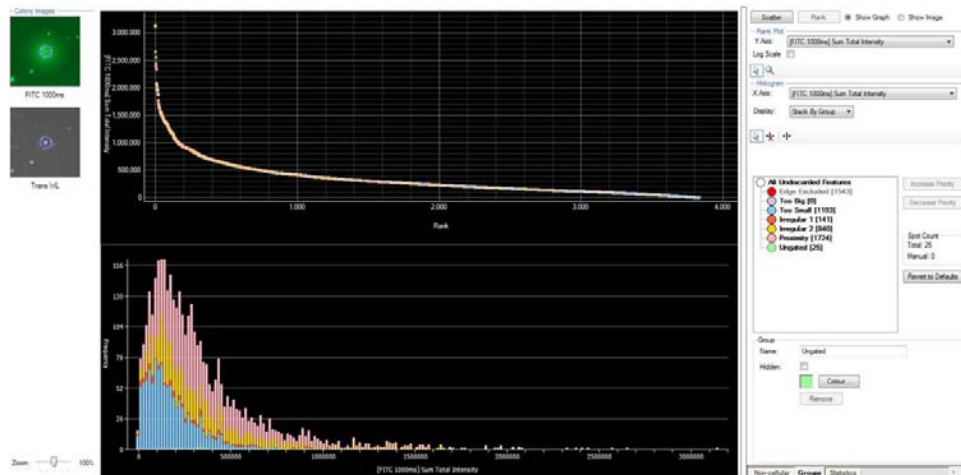
The grouping system works by priority with the group at the top having highest priority. For example, a colony that would fall into two groups is placed in the highest priority group.

Click **Increase Priority** and **Decrease Priority** to change the group priority. Colonies that do not fall into an exclusion group are retained in the bottom category. By default, this is named Ungated as these are 'good' colonies that have not been rejected by morphological characteristics or close neighbor proximity.

The bracketed number to the right of the group name indicates the number of colonies in the group.

## Spot Count

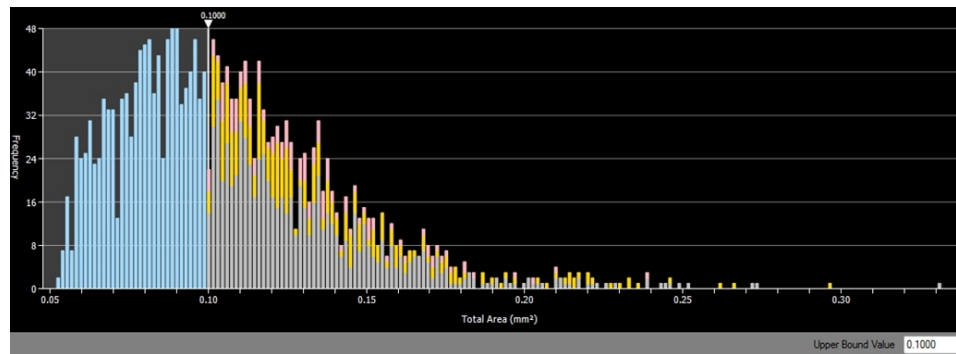
- **Total** - Displays the total number of colonies assigned to the groups you select.
- **Manual** - Displays the number of colonies you manually assign to the group.
- **Pickable** - When you select **All Undiscarded Features**, this displays the total number of pickable colonies.
- **Revert to Defaults** - Click to change all the group values back to the original default values.



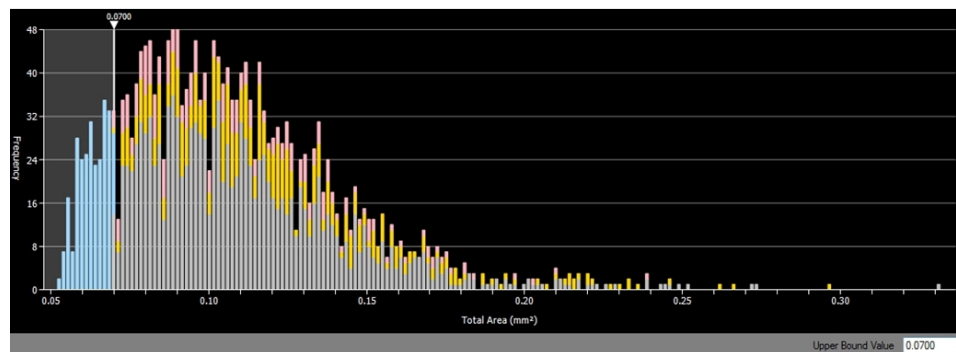
There are six default exclusion groups:

- **Edge Excluded** - Edge excluded colonies are outside the pickable region of the well and cannot be selected. This group has highest priority and cannot be moved to a lower priority.
- **Too Big** - Colonies greater in size than  $0.7\text{mm}^2$  fall into this group.
- **Too Small** - Colonies smaller in size than  $0.1\text{mm}^2$  fall into this group.
- **Irregular 1** - Colonies that have a Compactness value less than 0.6 fall into this group.
- **Irregular 2** - Colonies that have an Axis Ratio value less than 0.6 fall into this group.
- **Proximity** - Colonies that have a Proximity value less than 0.1 mm fall into this group.

Double-click the group name to display the group and its gate on the histogram where you can edit the group.



Move the slider on the histogram to a new value or enter the new value in the **Upper Bound Value** field in the lower right corner of the histogram (press **ENTER** to apply changes). The group gate displays the colonies that now fall into the altered group.



At the bottom of the Groups tab, you can enter a new name for the group and press **ENTER**. Select the **Hidden** checkbox to hide the group from view. This hides the group from view on all images and in the Data Table tab and Graphs tab. Click **Color** to change the group color.

Select a group and click **Remove** to delete the group. You cannot delete the Edge Excluded group or the Ungated group.


The bottom category (Ungated by default) is not a defined group and so does not have a gate on the histogram that can be changed.

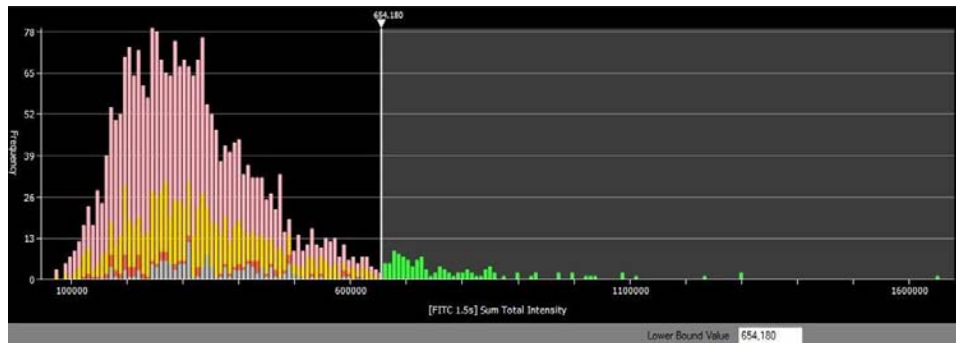
### Creating New Groups

Example: Creating a group for picking high FITC-expressing colonies.

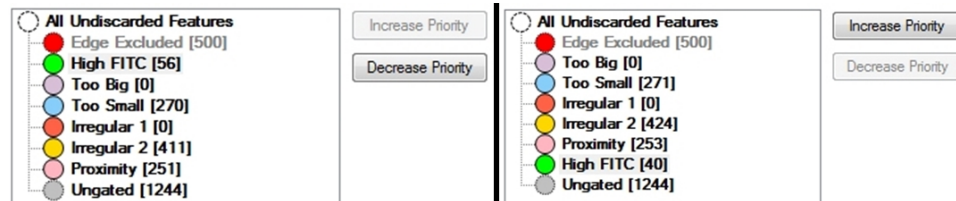
A white light Prime image and a FITC fluorescent image must be acquired to do this example.

To create a new group:

1. On the Graphs tab, click **Rank** to display the rank plot and histogram .
2. Select a fluorescent statistic from the drop-down list, for example, [FITC] Exterior Mean Intensity or [FITC] Sum Total Intensity. This displays on both graphs.
3. Click  **New Line Gate**, click the histogram where the group will start and drag the line off the end of the histogram to create the group.



4. Name the group, for example, High FITC, give it a color and click **OK**.  
The group displays below the Edge Excluded group.
5. Select the new group and decrease its priority so that it displays above the Ungated category.



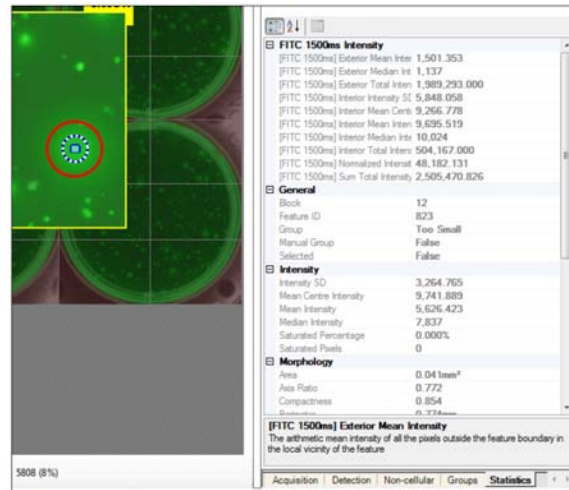
**Decrease Priority** - Only colonies with good morphological and proximity characteristics can be put into the new group.

All colonies in this group display in the Group color on the Images, Statistics, and Graphics tabs.

## Statistics Tab

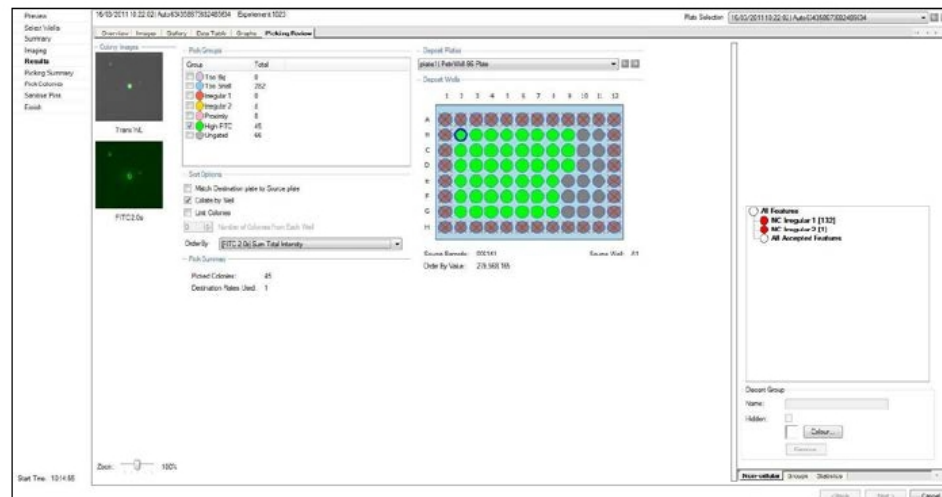
On the Statistics tab, you can view statistics for any colony on the Images, Data Table, Graphs, and Picking Review tabs.

Select the **Statistics** tab to display the statistics for a colony of interest. You can categorize the statistics in this list (Fluorescent Intensity, General, Intensity, Morphology, Position) or order the statistics alphabetically.



## Picking Review Tab

The Picking Review tab displays a summary of the colonies to be picked. Any number of groups can be selected to be picked here, and to be ordered using various criteria.



A summary of the picking review tab appears to the right of the main screen, including the following:

- **Colony Images** - When you select a well in Deposit Wells, the corresponding colony image displays on the left. Use the Zoom slider to zoom in and out of the colony images.
- **Picking Groups** - All the groups available to be picked are listed here, including exclusion groups such as Too Small and new groups that you create with specific characteristics.

- **Sort Options** - Provides several ways to sort colonies.
  - **Order by** - Allows colonies within a checked group to be picked in order of any of the parameters in the drop-down list, for example Fluorescent Exterior Mean Intensity. If you select several groups, you should order by Group.
- **Pick Summary** - Displays the number of colonies to be picked and the number of destination plates into which colonies are picked.
- **Deposit Plates** - Lists the number and type of destination plates (as previously defined in Modify Settings, Picking Settings) that are required for picking.
- **Deposit Wells** - Displays the wells that the colonies are going to be picked into. The wells are colored according to the corresponding colony that is picked into them. Any wells that have been excluded from picking display with red crosses over them. Underneath the well map, the following information displays when you hover the cursor over the colony assigned wells:
  - **Source Barcode** - Displays the barcode of the source plate where the picked colony comes from.
  - **Source Well** - Provides the source well number of the picked colony.
  - **Order By Value** - Provides the order value (assigned from the sort options drop-down list) of the colony to be picked.

## Picking Summary

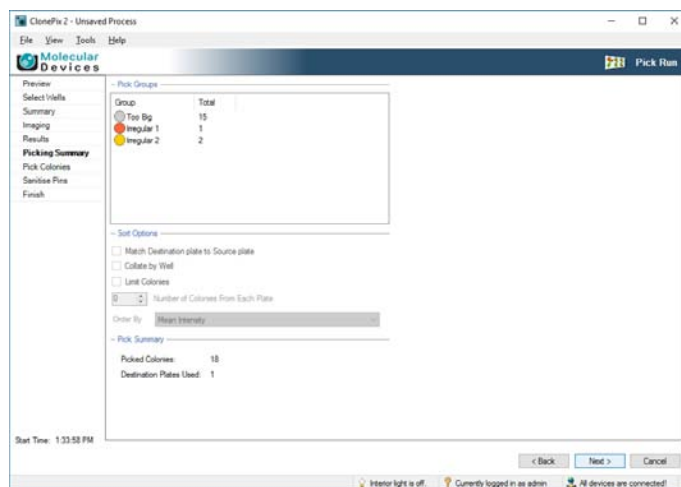
The Picking Summary screen displays the pick information before the picking starts.

Click **Next** to start picking.

To go back to the Results screen where you can make changes, click **Back**.



**CAUTION!** Selecting Back reprocesses all the images captured, which can take a while if several plates have been imaged.



When you start the picking, a message prompts you to load the destination plates into the destination stacker cassette. Raise or lower the destination stacker, if needed.

## Pick Colonies

The Pick Colonies screen displays the Picking Progress as the colonies are being picked. The automated picking step initiates with a pin wash.

**Picking Progress**

Microplates  
 Source Plate: Auto6342084869556353240  
 Destination Plate 1: Auto6342085253260786851

Picking colony 11 of 50

| Source                        | Destination                   | Thumbnail | Before | After |
|-------------------------------|-------------------------------|-----------|--------|-------|
| Auto6342084869556353240<br>A2 | Auto6342085253260786851<br>F3 |           |        |       |
| Auto6342084869556353240<br>B2 | Auto6342085253260786851<br>E3 |           |        |       |
| Auto6342084869556353240<br>A3 | Auto6342085253260786851<br>D3 |           |        |       |

Options  
 Change wash bottle  
 Pause  
 Pick: Safe  
 Terminate  
 Imaging  
 Image Before Pick  
 Image After Pick

Screen options include:

- **Microplates** - This section displays a schematic of the source plate and the destination plate. As colonies are being picked from a well of a source plate, the well is colored pink. As the colonies are being picked into the destination plate, the wells are colored pink. The barcodes of the plates being used appear and any wells not being used display a red cross over the well.
- **Colony Images** - As colonies are being picked, several pieces of information display; source plate barcode and well number, destination plate barcode and well number, a thumbnail image of the colony to be picked, a thumbnail image of the pin in place before the colony is picked and a thumbnail image of the pin in place after the colony has been picked. It is optional to have the before and after images display. Select or clear the checkbox before and after images.

- **Options** - Select from the options to stop the colony picking, including:
  - **Change Wash Bottle** - Enables sterile water to be replaced during the picking. To activate:
    - Click **Pause**
    - Select **Change Wash Bottle**, and then click **Continue**.  
Immediately before the next wash cycle, a prompt appears to swap over the bottle (if needed), and starts the long purge. The next cycle of pin washing reverts to the original purge settings.
  - **Pause** - Temporarily stops at any point during the picking.
  - **Park Safe** - When in Pause mode, you can move the picking head to the Park Safe position. Click **Continue** to restart the picking.
  - **Terminate** - When in Pause mode, the picking ends immediately. Use with caution.
- **Imaging** - When selected, the camera repositions and displays on-screen images of the pin tip before and after picking. Options include:
  - **Image Before Pick** - Select to display each colony immediately before being picked.
  - **Image After Pick** - Select to display if each colony has been successfully collected by the pin.

You can switch these options on and off at any time during the run.



**Note:** These images are not saved. If saving the images is required, see [Change Picking Settings on page 47](#) and select the Save Target and Aspirate Images.



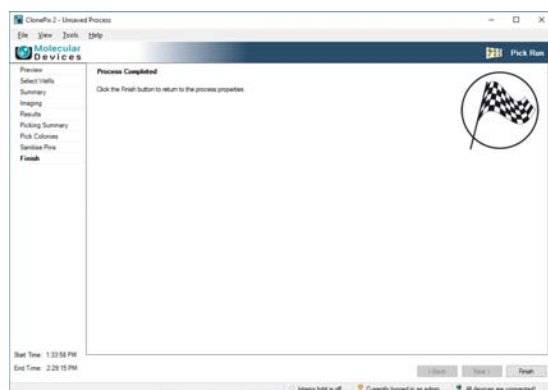
**Note:** Activating these images significantly slows the picking speed, and small colonies might not be visible.

## Sanitise Pins

In between each pick and at the end of each run, the pins are cleaned as specified in the Sanitise Pins Options set up. See [Change Sanitize Pin Options on page 49](#).

## Finish

When the picking ends, the Finish screen appears.



Click **Finish** to return to the Pick Run process Start screen. See [Start Pick Run on page 45](#)

To return to the Main Navigation screen click **Close Process**. If the process has not been saved, a prompt displays for the settings to be saved.

To view the results of the picking run, go to Review Results. See [Review Results Process on page 75](#).



## Imaging Run Process

Use the Imaging Run process for plate imaging without committing to a picking run. The default process includes all the parameters required for a standard imaging run, including image acquisition, colony detection, and group settings.



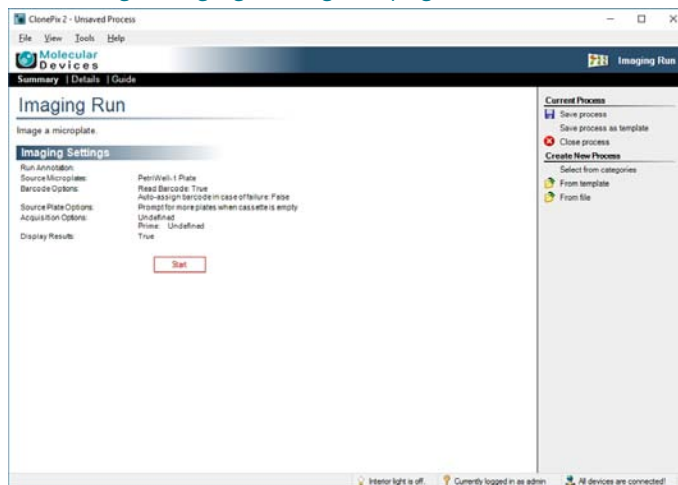
**Note:** Images acquired using this process cannot be used for picking later.

After imaging, source plates return to the stacker cassette, and then the process ends. Imaged microplate results are accessible in Review Results with data for all detected colonies. Imaging run data can be reprocessed in Review Results. See [Review Results Process on page 75](#).

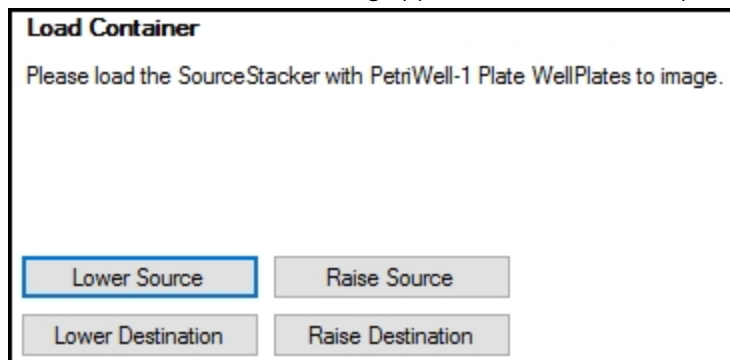
### Start Imaging Run

To start Imaging Run:

1. Before you continue, make setting changes. On the Imaging Run > Start screen, you can change the following settings:
  - [Change Imaging Settings on page 45](#)



2. Click **Start** to start the run process.
3. When the Load Container dialog appears, load the source plates into the source stacker.



4. Raise or lower the source stacker, if needed.
5. Click **Next**.

## Preview

The Preview screen starts the image acquisition and colony detection set up. This is the same process as Pick Run. See [Preview on page 49](#).

## Select Wells

Use the Select Wells screen to choose which wells to image. This is the same process as Pick Run. See [Select Wells on page 53](#).

## Summary

Use the Summary screen to confirm that all settings are correct before you proceed to imaging. This is the same process as Pick Run. See [Picking Summary on page 70](#).

## Imaging

The Imaging screen displays the images that are being captured as imaging progresses. The selected wells image in sequence for each of the acquisition options you select. The images are then processed together for colony detection according to the defined settings. This is the same process as Pick Run. See [Imaging on page 55](#).

## Results

If you select the Display Results option, see [Change Imaging Settings on page 45](#). The results appear after imaging. The Results screen has a similar format to the Preview screen. Use the Well Map to navigate the images and to view collated statistics for the entire plate. You can adjust Groups options. See [Groups Tab on page 66](#).

If you do not select the Display Results option, you can access the images and results from the Main Navigation screen > Review Results.

## Finish

The finish screen displays when the picking ends.

Click **Finish** to return to the Imaging Run process Start screen. See [Start Imaging Run on page 73](#).

Click **Close Process** to return to the Main Navigation screen. If the process has not been saved, a prompt displays for the settings to be saved.

To view the results of the picking run, go to Review Results. See [Review Results Process on page 75](#).

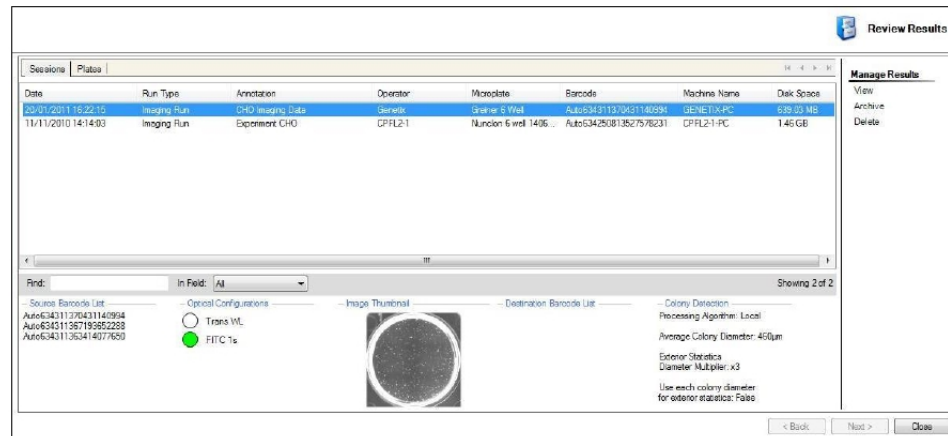
## Review Results Process

Use the Review Results process screen to review a completed Imaging Run and Picking Run.

By default, data is stored on the C: drive in a folder named Image Archive.

To open the Image Archive, click **Next** or to access data remotely, click **Browse**.

The Sessions tab displays a list of completed runs:



The runs appear in two tabs:

- **Sessions** tab - Results appear as one row even if the run consisted of several plates.
- **Plates** tab - Results appear as one row per plate processed so if there were three plates in the run, three individual rows appear.

Click the column headers to sort the results list or a double-click to invert the selection.

To view the data, do one of the following:

- On the **Sessions** tab, either double-click on a row or click once on a row and then click **View** to open the data set.
- On the **Plates** tab, do the same as in the Sessions tab, or to view multiple plates, use Shift-Click or Control-Click to highlight the rows, and then click **View**. You can also combine the data from multiple plates and view and process the plates as a batch.
- Search the data sets using **Find**. The following fields are available to search: **Date**, **Run Type**, **Annotation**, **Operator**, **Microplate**, **Barcode**, and **Machine Name**.

When you select a data set, the following information displays:

- **Source Barcode List** - Lists the barcodes on the source plates used.
- **Optical Configurations** - Lists the optical configurations used in the run and the colors associated with them.
- **Image Thumbnail** - Displays the prime configuration well image.
- **Destination Barcode List** - Lists the barcodes on the destination plates used.
- **Colony Detection** - Displays a summary of the parameters used to detect the colonies.

After you open data for viewing, depending on whether imaging or picking data is being opened, the tabs available are the same as seen throughout the Pick Run and Imaging Run stages. In a picking run data set, the Picking Review tab is visible. See [Picking Review Tab on page 69](#).



**Note:** For data integrity, Pick Run data is fully accessible in Review Results but the data are locked and cannot be changed. Imaging Run data can be changed but cannot be saved.

- **Archive** - Stores data to Image Archive or transfers data to CD or DVD. Shift-click or Control-click to highlight the required rows and then click Archive.
- **Delete** - In Image Archive, select a data set and click **Delete** to remove the data from the hard drive.

## Export Results Data

For Pick Runs, an Output Results tab displays destination plate, well and barcode information. Pick Run data are non-modifiable to provide an accurate record of the run.

For Imaging Run, data are modifiable on the Detection tab and Group tab, but your changes are not saved when you close the file.

You can export Graphs and Images or copy them to the Windows clipboard or saved as a file. You can paste them into presentations, or other applications for sharing.

You can export Statistics with a choice of data formats including comma-separated values (CSV), XML, and Excel-XML. On the Statistics tab, click **Export Statistics**.

Generate a Results Report to provide a summary of a run including its settings. Create it on the Review Results process screen or the Remote Data Viewer software in XML format so that it can be opened in a web browser and printed or stored electronically. To create a Results Report from the Statistics tab, click **Generate Report**.

## Remote Data Viewer

Remote Data Viewer software is an optional software to access ClonePix 2 software results remotely on another computer (order code: SL4950-A01). You can install it on one computer. The format is identical to the Review Results features in ClonePix 2 software, and allows you to access results from a ClonePix 2 system through a network or copies of runs transferred elsewhere.

## Data Tracking Logs

A Data Tracking Log file is created for each run. This is an XML file with the same schema used by other instruments. It includes some statistical information for each colony along with a 'replicate' element for each transfer from a source well to a destination well.

The Data Tracking log can be of interest for those wishing to automatically process the transfer of samples between plates. However the Automatic Logs option can be more useful.

Data Tracking Logs are generated in the 'Data Tracking' sub folder of the configuration folder. The configuration folder is stored under the local applications data folder.

On Windows 7 and Windows 10:

C:\ProgramData\Molecular Devices\Fusion\ClonePix 2\DataTracking

One file is generated for each imaging or picking run. The file name is derived from the data and time of the run, and the file extension is XML.

## Automatic Logs

---



**Note:** To enable the Auto Logs, ask Molecular Devices Technical Support.

---

Auto Logs can generate by the system for each run with some level of control of their content.

---



**Note:** Auto Logs do not generate by default. When enabled, one log is generated for each run.

---

The log is broadly the same format as that provided by the Export Statistics in Review Results or the Remote Data Viewer. See [Export Results Data on page 76](#) and [Remote Data Viewer on page 76](#).

When the Auto Logs function is active, the folder generates for each run with a name derived from the data and time of the run. A file saves in the folder with the name Auto.csv or Auto.xml, depending upon the file-type selection.

## Raw Data

The default location for raw data (TIF images and XML files) is: C:\image archive.

---



**CAUTION!** Viewing data from this location is not recommended.

---



Utility Processes provide several processes to control and maintain the hardware.

### Picking Head Management

Use Picking Head Management functionality each time the picking head is removed or replaced.

#### Removing the Head

To remove the picking head:

1. Click the **Remove Head** icon to bring the head to the removal position.
2. On the picking head, remove all of the tubing from the left side of the head.



**Tip:** Place a microplate lid under the picking head to collect any drops of water.

3. On the right side, loosen the thumbscrew that secures the head to the actuator assembly.



**CAUTION!** Do not lose the washer.

4. Slide the head out of the actuator while being careful not to damage the pins.
5. To return the actuator to a safe position, close the door and click **Next**.

For picking head cleaning procedures, see [Cleaning the Picking Head on page 88](#).

#### Replacing the Head



**CAUTION!** When handling the picking head take care that the pins are not knocked. Damage to pins can produce unreliable results and can cause serious damage to the instrument. If you suspect any pin damage, replace them immediately.

To replace the picking head:

1. Click **Replace Head** to bring the head to the replacement position.
2. Slide the head on to the actuator while being careful not to damage the pins.
3. On the picking head, attach the picking tubing, and on the right side, tighten the thumbscrew that secures the head to the actuator assembly.
4. Replace the tubes starting from the back, pushing on the picking tubing.
5. Click **Next** to continue to Pin Fire Test. See [Pin Fire Test on page 35](#).
6. Click **Next** to continue to Alignment for plate and primary pin. See [Alignment on page 36](#).
7. Click **Next** to continue to Pin Offsets. See [Adjusting Pin Offsets](#).

## Adjusting Pin Offsets

The Pin Offsets process aligns the remaining seven pins relative to the primary pin. Run this whenever you remove or replace the pins compensate for  $\mu\text{m}$  differences between pins.



**CAUTION!** If you touch the pins or move the pins in any way, you must adjust the pin offsets for pin one and pins two through eight.

To adjust pin offsets:

1. Following on from Alignment, see [Alignment on page 36](#), you are prompted about using the same plate. To continue, click **Next**, or to cancel, **Close**.
2. To continue and extend the Primary Pin into well A1, click **Next**.

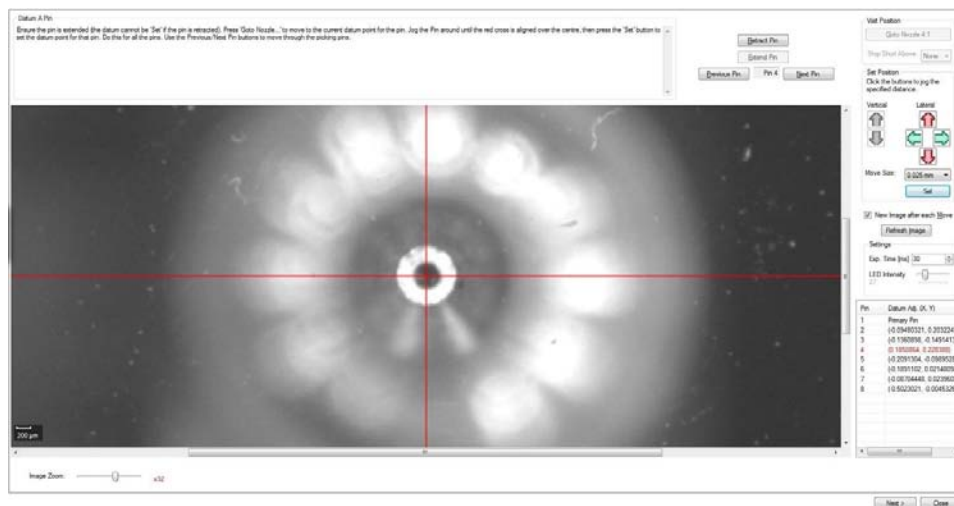


**Note:** The datum point for the primary pin cannot be set here. It should already have been set in the Alignment procedure. If the primary pin is misaligned, exit the Replace Head process and start again.

3. Click **Next Pin** to extend the next pin into well A1.
4. Align the pin using the X and Y axis Jog buttons until the pin hole is centered on the red cross.
5. When the pin is in the required position, click **Set**. The new pin coordinates appear in red.
6. To move through and set the remaining pins, click **Next Pin**.



**Note:** If you click Previous Pin, the pin might appear not to be aligned because the actual position is shown rather than the offset on the assumption that you want to create the offset.



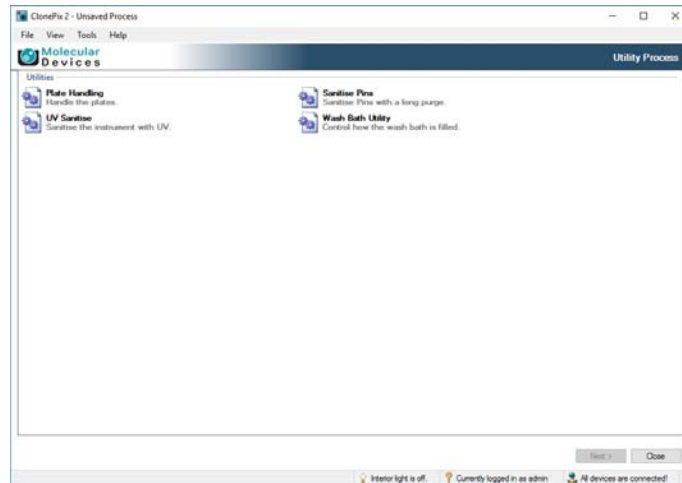
**Pin Correctly Aligned to Camera and Pins 2 and 3 Already Aligned Shown in red**

7. Click **Next** for Wash Bath Utility. See [Wash Bath Utility on page 40](#).
8. Click **Next** for Sanitize Pins in 70% Ethanol. See [Alternative: Sanitize Pins in 70% Ethanol on page 41](#).
9. Click **Next** for Ultra Violet Sanitize. See [Ultra Violet Sanitize on page 41](#).
10. Click **Next** for Sanitize Pins to purge the system with sterile water. See [Sanitize Pins on page 42](#).
11. Click **Next** twice to automatically park the head safely, return the source plate, and close the illumination cover.



## Utility Process

The Utility Process screen includes much of the same functionality as the Prepare for Pick Run process, with the exception of Plate Handling settings. See [Prepare for Pick Run on page 35](#).



- **Plate Handling** - Set Source Plate and Destination Plate handling.

To set source plate handling:

- For **Plate Station**, select **Source Stacker**.
- For **Microplate**, select the plate type to use.
- Click **Get Plate** and confirm that the plate collects from the Source Plate Station.
- Click **Return Plate** and confirm that the plate returns to the Source Plate Station.

To set destination plate handling:

- For **Plate Station**, select **Destination Stacker**.
- For **Microplate**, select the plate type to use.
- Click **Get Plate** and confirm that the plate collects from the Destination Plate Station.
- Click **Return Plate** and confirm that the plate returns to the Destination Plate Station.

- **Sanitize Pins** - See [Sanitize Pins on page 42](#).
- **UV Sanitize** - See [Ultra Violet Sanitize on page 41](#).
- **Wash Bath Utility** - See [Wash Bath Utility on page 40](#).



Before operating the instrument or doing maintenance operations, make sure that you are familiar with the safety information in this guide. See [Safety Information on page 6](#).



**WARNING!** Service or maintenance procedures other than those specified in this guide can be done only by Molecular Devices qualified personnel. When service is required, contact Molecular Devices Technical Support. See [Obtaining Support on page 96](#).

### Doing Preventive Maintenance

You should do daily and weekly maintenance. We strongly recommend that a complete instrument preventative maintenance service be done annually by Molecular Devices.

#### Daily Maintenance

- Check that the instrument interior is clean and free from any obstruction.
- Ensure sterile water feed bottle is connected to the system fluid supply line. Empty the waste bottle every time that a new sterile water bottle is connected.
- If the system has been sanitized, purge the system fluid supply line on first use to remove air bubbles that develop overnight. If there is doubt about sterility, run the Prepare for Pick Run routine that includes a purge of the system fluid supply line.
- During the purge, check that there are no air bubbles in the supply line coming from the sterile water bottle, which will indicate a fault in the connection, and that there are no leaks around the picking head, which will indicate a blocked pin. If there is a leak, or if a tube detaches from the head, replace blocked pin with a new one or follow the pin cleaning procedure. See [Cleaning the Head and Pins on page 87](#).
- Fill ethanol wash bath with 70% ethanol. For systems fitted with a replenishing ethanol wash bath, check that there are no kinks in the tubing and then run through with 70% ethanol. Check that the bath fills with ethanol and that the waste ethanol flows freely into the waste bottle. On first use, it might be necessary to run through several times.

#### Weekly Maintenance

- Any week that the ClonePix 2 system is not used, it is important to purge the system fluid supply line with sterile water and sporicide. Failure to keep the pumps regularly primed may lead to system malfunction.
- Sonicate the picking pins.
- Check operation of interlock switch and emergency stop.
- Check the compressed air filter for signs of moisture. If necessary, push the drain upwards to force moisture out.
- If you use the external air compressor, drain the accumulated moisture from the internal air tank weekly. If you use the external air compressor, drain the accumulated moisture from the internal air tank.

#### Monthly Maintenance

- Sanitize the picking head and silicon fluid tubing.
- Sanitize the source plate illumination cover.

### Annual Maintenance

A complete instrument maintenance should be done every six months by an approved service engineer. To obtain a maintenance contract or schedule a service visit, contact your representative or technical support.

### As needed


- If the instrument is used in dusty conditions, check the HEPA filter or pre-filter and have Molecular Devices replace it more frequently than recommended.

## Changing Pins

The pins need to be removed from the picking head for cleaning and to change to a different pin type.



---


 **CAUTION!** If you touch the pins or move the pins in any way, you must adjust the pin offsets for pin one and pins two through eight. See [Adjusting Pin Offsets on page 80](#).

---

To change pins:

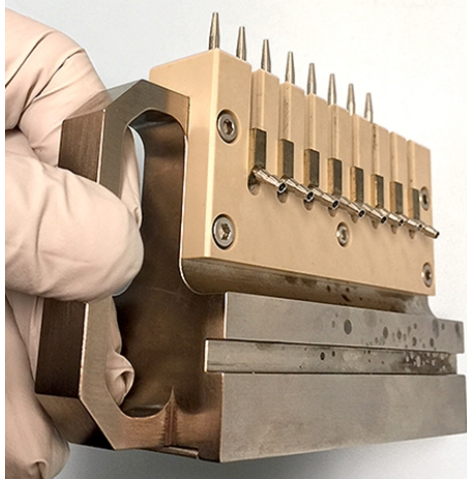
1. Click **Replace Head** to bring the head to the replacement position.
2. On the picking head, remove all of the tubing from the left side of the head.
3. On the right side, loosen the thumbscrew that secures the head to the actuator assembly.

---

 **CAUTION!** Do not lose the washer.

---

4. Slide the head off of the actuator and continue working over a workbench



5. On the picking head, use the Pin Removal Key (X4948) to turn counter-clockwise and loosen each of the pins.



6. Before installing the new pins, verify that each pin has an O-ring (X1036) on the thread.



7. When installing each pin into the picking head, use the Pin Removal Key to turn the pin clockwise until you feel some resistance.



**CAUTION!** Do not over-tighten.

---

8. Slide the head onto the actuator while being careful not to damage the pins.
9. On the picking head, attach the picking tubing, and on the right side, tighten the thumbscrew that secures the head to the actuator assembly.
10. Replace the tubes starting from the back, pushing on the picking tubing.
11. Click **Next** to continue to Pin Fire Test. See [Pin Fire Test on page 35](#).
12. Click **Next** to continue to Alignment for plate and primary pin. See [Alignment on page 36](#).
13. Click **Next** to continue to Pin Offsets. See [Adjusting Pin Offsets on page 80](#).

## Cleaning the Instrument

Follow all of the cleaning procedures outlined in this guide for the instrument.

Do the following before you clean equipment that has been exposed to hazardous material:

- Contact the applicable Chemical and Biological Safety personnel.
- Review the Chemical and Biological Safety information contained in this guide. See [Chemical and Biological Safety on page 10](#).

Avoid spilling liquids on the system. Fluid spilled into internal components creates a potential shock hazard. Wipe up all spills immediately. Do not operate the system if internal components have been exposed to spilled fluid. Unplug the instrument if there is a fluid spill in the instrument and contact Technical Support.

Perform only the maintenance tasks described in this guide. Contact a Molecular Devices service engineer to inspect and perform a preventive maintenance service on the instrument each year. See [Obtaining Support on page 96](#).



**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.

---

Clean all components that come into close contact with biological materials. For efficient decontamination of pathogenic micro-organisms, wipe all non-removable parts within the instrument with a cloth using 70% ethanol.



**CAUTION!** Molecular Devices recommends that you always use ethanol for cleaning, because autoclaving is not compatible with anodized parts. Do not use abrasive cleaners, because it damages the surface of the bed.

---

The instrument can be left in a laboratory during formaldehyde vapor fumigation at an appropriate concentration.



**CAUTION!** Excessive formaldehyde treatment damages sensitive electrical and optical components.

---

## Sterility Validation Tests

To test the sterility of your environment:

### Chamber Sterility Test

To test chamber sterility:

1. Prepare the instrument using the cleaning procedures provided.
2. Place a Trypticase Soy Agar\* plate open on the process bed during a prolonged mammalian cell picking run.  
\*A general purpose medium which supports the growth of fastidious and non-fastidious microorganisms.
3. After 20 minutes, cover the plate, remove the plate, and replace the plate with another open plate.
4. Repeat steps 1 and 2 six times spanning two hours.
5. Incubate the plates at 37°C for two weeks to encourage any potential contaminant growth. The chamber is sterile when after this period there is no contamination on any of the agar plates.

### Mammalian Cell Picking Sterility Test

To test mammalian cell picking sterility:

1. Prepare the instrument using the cleaning procedures provided.
2. Plate CHO-S cells in chemically defined medium without antibiotics and render them semi-solid using CloneMatrix.
3. Incubate the plates at 37°C for 10 days to grow the cells into clonal colonies.
4. Select colonies (n=175) using the instrument into 96-well plates prefilled with 200RI liquid medium without antibiotics.
5. Grow the picked cells for 12 days, replenish with fresh liquid medium without antibiotics, and then expanded to exhaustion for four weeks.  
Sterility is confirmed when after four weeks there is no visible contamination in any of the wells.

## Cleaning the System Fluid Supply

The cell collection process relies on the presence of non-expandable system fluid between the pumps and the picking pins. The ClonePix 2 system uses sterile water for the system fluid supply. The system fluid is fed in from a detachable wash bottle and the waste fluid is voided directly to the waste bottle. The system fluid is also used to purge out the picking pins between each cycle of picking.

Change the Wash and Waste bottles using the quick-release connectors with the check valves that close when the connection is broken. Molecular Devices recommends liberally spraying the connectors with Spor-Klenz Ready-To-Use Cold Sterilant before re-connecting.

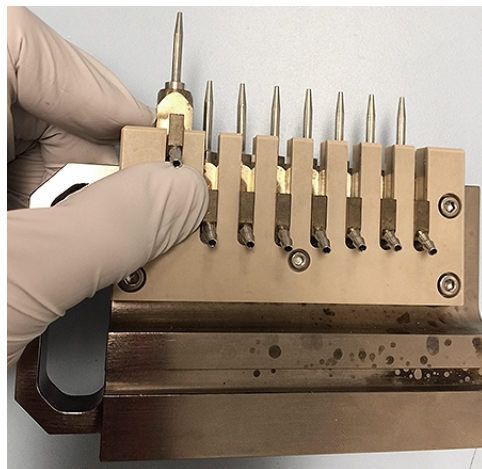
To ensure a sterile supply of system fluid, sanitize the system fluid lines once a week, and after any length of non-use. Follow the software procedures in [Prepare for Pick Run on page 35](#).

## Cleaning the Head and Pins

The reusable pins in the head are sanitized when you run a process that includes a Sanitise profile. They are cleaned before the first pick, between each cycle of picking, and at the end of the run.

You can leave the picking head installed and spray it with 70% ethanol and let it dry in the HEPA laminar air flow.

For a more thorough cleaning of the head and pins, remove the head from the actuator. See [Removing the Head on page 79](#).



The head is housed in an actuator system that permits easy exchange and set-up of the head. Although it is possible to manually move the actuator assembly, Molecular Devices recommends using the software to safely move the actuator into position to remove or install the head. See [Picking Head Management on page 79](#).



**WARNING!** The actuator assembly has moving parts that can cause pinch injuries if it is moved manually. To prevent pinch injuries, use the software Change Head process to safely move the actuator.

Before storing the head, or whenever it needs cleaning, wipe the exterior of the head with a cloth using 70% ethanol.

To store the head, slide it into its metal cover with the pins facing inward and then secure it in place with the thumbscrew and washer.



**Note:** Automated cleaning with a Sanitise profile does not replace sonicating the pins. See [Sonicating the Pins on page 90](#).

## Cleaning the Picking Head

You can thoroughly clean the picking head with either ethanol or a sterilizing agent.



**CAUTION!** Autoclaving the head is not recommended because it can cause functional deterioration. Molecular Devices recommends at least one picking head for each project.



### Picking Head Without Pins

#### Cleaning the Picking Head With Ethanol

To clean the picking head with ethanol:

1. Remove picking head from the instrument and remove the pins and the picking tubing. See [Removing the Head on page 79](#).



2. Submerge the head and O-rings into 100% ethanol.
3. Press the piston actuator to allow the interior to fill.



4. Soak for 10 minutes.
5. In a sterile environment, lift the head out of the ethanol.
6. Press the piston actuator to allow the interior to empty.
7. Leave to air dry.
8. Replace the picking head. See [Replacing the Head on page 79](#).



## Cleaning the Picking Head With Sterilizing Agent

To clean the picking head with sterilizing agent:

1. Remove picking head from the instrument and remove the pins and the picking tubing. See [Removing the Head on page 79](#).
2. Prepare the sterilizing agent in 1 liter of warm water.
3. Submerge the head into the sterilizing agent.
4. Press the highest piston to allow the interior to fill.
5. Soak for 10 minutes.



**CAUTION!** Do not leave the head in the sterilizing agent for longer than 10 minutes because acetic acid generated as a by-product can cause discoloration or functional deterioration.

6. In a sterile environment, lift the head out of the sterilizing agent.
7. Press the piston actuator to allow the interior to empty.
8. Submerge the head into a beaker of sterile water.
9. Press the highest piston to allow the interior to fill.
10. Remove the head from the sterile water.
11. Press the piston actuator to allow the interior to empty.
12. Repeat steps 8-11 multiple times to remove the sterilizing agent.
13. Hold the head over an empty beaker and spray it liberally with 100% ethanol including the interior.
14. Press the piston actuator and ensure that the interior is well drained.
15. Leave to air dry
16. Replace the picking head. See [Replacing the Head on page 79](#).

## Cleaning the Picking Pins



**Note:** Molecular Devices recommends at least two sets of Picking Pins for each project.

The reusable Picking Pins are frequently cleaned automatically during a picking run. Specifically, they are cleaned before the first pick, between each cycle of picking, and at the end of the run.

There are four components to the automated cleaning process:

- Sterile water unidirectional purge to clean the internal bores of the picking pins.
- Ethanol Wash Bath brushing of the picking pins for effective mechanical cleaning of the pin exterior.
- Halogen dryer to heat the picking pins and remove residual ethanol.
- Sterile air blower blows over the picking pins to cool them and ready them for the next pick.



**Note:** The length of each of the above steps can be user defined according to the application, but the halogen dryer should not be run for less than 10 seconds.



**CAUTION!** The automated cleaning is not designed to replace the sterilization of picking pins by sonication.

Sterilize the Picking Pins, O-rings, and Pin Removal Key regularly through sonication. See [Sonicating the Pins](#).

### Sonicate the Pins

You should sonicate the pins weekly. Before sonicating the pins, remove the pins from the head and sonicate the pins only.

To clean the head, wipe it with a cloth using 70% ethanol. If necessary for further decontamination, soak the head in 100% ethanol. Before reassembly, make sure that all parts are thoroughly dry.

If necessary, you can soak the springs and screws in 100% ethanol and then let them thoroughly dry.

To sonicate the picking pins:

1. Prepare a 2% solution of aQu Clean Microarray Pin Cleaning Solution.
2. Sonicate the pins in the 2% aQu Clean solution for 10 minutes.
3. Rinse the pins with deionized water.
4. Sonicate the pins in deionized water for 10 minutes.
5. Rinse the pins with 100% ethanol, or autoclave.
6. Allow the pins to thoroughly dry in a biosafety hood for a few hours or overnight.

### Cleaning the Wash and Waste Bottles

We recommend at least three Wash Bottles and one Waste Bottle for each project. One Wash Bottle should be kept for cleaning the system fluid supply with 70% ethanol or sterilizing agent. The Wash Bottles that are connected to the Feed tubes must be sterile before use.

To clean the wash bottles:

1. Fill the Wash Bottle with deionized water.
2. Replace the bottle cap loosely.



**CAUTION!** Failure to leave the bottle cap on loosely can damage the bottle.

---

3. Seal the air inlet pipe with an autoclavable 0.2  $\mu\text{m}$  (microns) disposable filter.
  4. Autoclave using standard conditions: 121°C at 15 psi (103 kPa) for the required time. The time required to ensure sterility depends on your autoclave system.
- 



**CAUTION!** Do not autoclave the bottles filled with alcohol or sterilizing agent.

---

### Cleaning the Wash and Waste Tubing

The Wash and Waste Tubing are silicon tubes that connect the Wash Bottle and Waste Bottle.

Remove the tubes and autoclave as required using standard conditions: 121°C at 15 psi (103 kPa) for 20 minutes.

## Cleaning the Picking Tubing

The Picking Tubing consists of eight silicon tubes that connect between the picking head and the actuator.

The cell collection process relies on sterile water as non-expandable system fluid between the pumps and the picking pins. The system fluid is fed in from a detachable wash bottle and the waste fluid is voided directly to a detachable waste bottle. The system fluid is also used to purge out the picking pins between each cycle of picking. See [Cleaning the Picking Pins on page 89](#).



**Note:** To separate the sample from the system fluid, at the tip of the picking pins, is a small amount of air, default 2RI, drawn in from the sterile interior immediately before sample collection.

These tubes can be cleaned along with the picking head.



**Tip:** Spor-Klenz Ready-To-Use Cold Sterilant (VZR #21899-864) sterilizing agent can be used to stop fungal growth, while ethanol will not. Thoroughly flush out all Spor-Klenz sterilizing agent before using the instrument again.



**CAUTION!** Never clean the system fluid tubes with bleach.



**Note:** If you autoclave, use standard conditions: 121°C at 15 psi (103 kPa) for 20 minutes.

## Cleaning the Ethanol Wash Bath

The pins are cleaned during the process in the stainless steel Ethanol Wash Bath. The wash bath includes a nylon bristle insert. During normal use, the wash bath is filled with 70% ethanol.



**Tip:** When the instrument is not in use, Molecular Devices recommends that you remove and rinse the nylon bristle insert with distilled water.

If contamination is suspected, the wash bath and nylon bristle insert can occasionally be autoclaved using standard conditions: 121°C at 15 psi (103 kPa) for 20 minutes. Slight warping of the nylon bristle insert after autoclaving is normal, which you can force back to straightness.



**CAUTION!** A severely warped block must not be used because it can damage pins and the instrument.

## Cleaning the Source Plate Imaging Cover

You can remove the Source Plate Imaging Cover and clean it with ethanol.



### Source Plate Imaging Cover Upside Down



**Note:** Autoclaving will discolor the cover but it will not change the functionality.

## Draining the Compressor

The following procedure is for the optional DynAir DA7001CS air compressor. The compressor collects water in the internal air tank. This water must be emptied into the externally mounted fluid collection bottle. Draining frequency depends on instrument use:

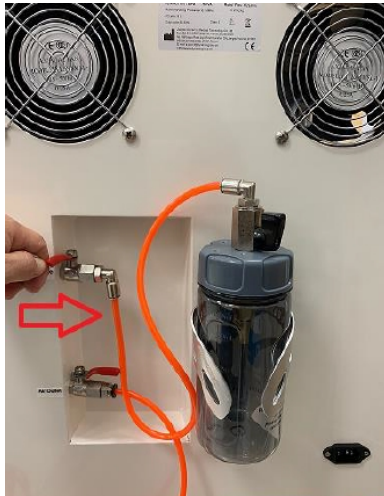
- High use - drain once a day
- Low use - drain once a week



To drain the air tank:

1. Verify that the compressor pressure is 105 psi.
2. Turn off the air compressor.

3. Slowly open the valve by turning the lever on the Drain Valve to the right about a quarter of the way open until you hear the sound of high pressure blowing air.



4. When the sound of air stops and fluid appears in the fluid collection bottle, close the valve by turning the lever all the way back to the left.



To empty the external fluid collection bottle:

1. Remove the fluid collection bottle from the metal holder, but leave the bottle cap lever vertical (open) and tube attached.
2. Unscrew the fluid collection bottle cap, and dispose of the fluid as required by your lab protocol.
3. Screw the fluid collection bottle cap back on, and return the bottle to the metal holder.

## Moving and Storing the Instrument



**Note:** We recommend not moving the instrument after installation. If relocation is required, standard lifting gear is adequate but only do the move with Molecular Devices supervision.

---

The ClonePix 2 system has caster wheels for ease of movement.

If you use a forklift or dolly trucks to move the instrument, it must be properly balanced before lifting.

---



**CAUTION!** Do not use any part of the exterior bodywork to lift the instrument as this can cause irreparable damage.

---

The instrument must be stored and transported in temperatures within the range -13°F (-25°C) to 131°F (55°C).

## Commissioning Process

The Commissioning Process must be done by Molecular Devices field service engineers. Contact Technical Support. See [Obtaining Support on page 96](#).

## Troubleshooting

This section describes some common problems that can occur and possible solutions.

Generally, crash recovery involves some form of the following procedure:

1. An error is identified in the software or the instrument stops.
2. If the error condition is caused by a receptacle or other item, then remove the item from the instrument.
3. Reactivate the software, generally by clicking **Next**.

### Troubleshooting Barcodes

If you use barcodes and the instrument does not detect them correctly, try the following:

- If one out of 10 plates fails to be read, check the placement of the barcode label on the plate. This might be fixed with a small change in the plate configuration.
- Add more white space at beginning or end.
- Restrict numbers to a maximum of six (6) digits.
- Verify that the codes meet the requirements. See [Using Barcodes on page 29](#).

If detection is still a problem, contact Molecular Devices Technical Support. See [Obtaining Support on page 96](#).

## Common Problems and Possible Solutions

### Machine will not turn on.

- Check that main power switch is turned on and the emergency stop button is released (pulled out).

### Computer will not start.

- Check that computer is powered on.

### One or more of the axes will not move.

- Check that the main power is turned on and the START button has been pressed.
- Manually move the actuator to the middle of the bed.
- Disconnect and reconnect the control cable between the computer and right side of the instrument. A small screwdriver might be needed.
- Re-start the instrument.

### Interior light or UV germicidal lamp will not turn on.

- Check that the bulbs are not burned out.

### Picking alignment is incorrect.

- Ensure that the pins are correctly aligned.
- Check for any bent pins and replace as needed.

### Silicon tubing becomes detached from a pin during the purge.

- The pin might be blocked. Replace with a new pin.

### Water is leaking from picking head or actuator.

- The pin might be blocked. Replace with a new pin.

### Poor picking results.

- Air in the system. Purge the supply line.
- Check Z height of the pin.

### Not depositing.

- The pin is blocked. Run the Purge option or replace the pin.
- Dispense volume is too low.

### No liquid flow to Wash Station.

- System fluid supply bottle is empty
- Check the air supply and the liquid tubes for kinks.
- Check that the outputs in the configuration are correct. Contact Molecular Devices Technical Support [support.moleculardevices.com/](http://support.moleculardevices.com/).

### Wash Station is overflowing.

- Check that no kinks are in the ethanol waste tubing.
- Waste bottle full and/or vacuum pump is broken.

### Pins are catching on the holes of the wash station.

- Datum point is not set accurately.
- The pins are damaged. Replace the pins.
- The pins have two O-rings. Remove one.

### Water is leaking from the top of pins.

- Make sure each pin is fitted with one o-ring.

**Pins are not firing correctly.**

- Check that air compressor is working.
- Pistons might have accumulated residue. Run Pin Fire Test for up to 1 hour.



**CAUTION!** Do not run Pick Run Fire Test if the pins are not firing correctly.

---

**Illumination cover is not opening or closing correctly.**

- Check that the air compressor is working.

**No ethanol flow to the Replenishing Ethanol Wash Bath.**

- Check that no kinks are in the ethanol waste tubing.
- Peristaltic pump tubing has become blocked, usually after period of inactivity, and needs clearing.

**Replenishing Ethanol Wash Bath is overflowing.**

- Check that no kinks are in the ethanol waste tubing.
- Peristaltic pump tubing has become blocked, usually after period of inactivity, and needs clearing.

**Error: One or more drivers on the machine failed to home correctly.**

- See Knowledge Base Answer number [20364](#).

## 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—[support.moleculardevices.com/](http://support.moleculardevices.com/)—describes the support options offered by Molecular Devices, including service plans and professional services. It also has a link to the Molecular Devices Knowledge Base, which contains documentation, technical notes, software upgrades, safety data sheets, and other resources. If you still need assistance, you can submit a request to Molecular Devices Technical Support.

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.

---



## Generating Error Reports

When the instrument is in an error state, you can generate an error report. This report is required when you contact Molecular Devices Technical Support.

To generate an error report:

1. Select **Tools > Configuration**.
2. Select **System Log**.
3. In the **System Log** list, select the most recent Log.
4. In the **Event Severity** section, select the **Error** checkbox and the **Fatal** checkbox.
5. To review the details, highlight the results lines.
6. To generate a report of the results listed, select **Tools > Prepare Error** report to display the Error Reporting wizard.
7. Click **Next**.
8. Specify the location to save the error logs.
9. Send the report file to Molecular Devices Technical Support.



## Appendix A: Technical Specifications



If you use the air compressor, see the documentation that comes with the compressor for details.

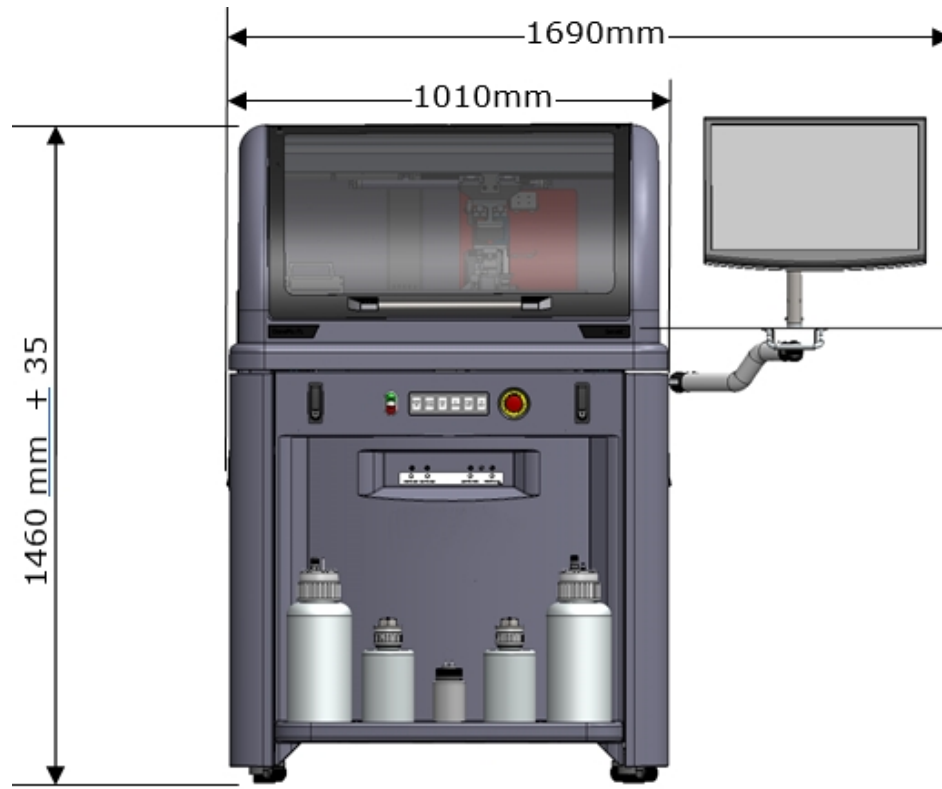
### ClonePix 2 Mammalian Colony Picker Technical Specifications

| Item                       | Description   |
|----------------------------|---|
| <b>Imaging</b>             |   |
| Software                   | Dedicated imaging software pre-installed on a high-specification PC, Microsoft Windows 7, 32-bit, and Windows 10 64-bit   |
| White light imaging        | Trans-illumination for imaging low contrast colonies such as adherent monolayers or small colonies in suspension<br>Epi-illumination for imaging colonies as they are collected |
| Fluorescence imaging       | Software-controlled switching between up to 5 excitation /emission filter pairs (we recommend no more than 3 filters be multiplexed for optimal performance)                    |
| Data tracking              | Internal barcode reader for source and destination plates enables data tracking for each run  |
| Camera                     | Integrated 16-bit cooled CCD camera   |
| Imaging speed              | 6-well microplate: 5 min for 2 wavelengths (standard conditions)  |
| Resolution                 | Standard: 28 micron   |
| <b>Instrumentation</b>     |   |
| Containment                | Fully enclosed working environment with Class 100-type, HEPA filtration   |
| Source plate type          | See <a href="#">Plate Compatibility on page 103</a>   |
| Destination plate type     | See <a href="#">Plate Compatibility on page 103</a>   |
| Source plate capacity      | 10 plates   |
| Destination plate capacity | 10 plates   |
| Picking head               | 8 x picking pins – each pin independently controlled  |
| Picking pin size           | Diameter of picking pins is application specific –F1: suspension cells (0.4 mm), F2: adherent cells (0.7 mm)  |
| Picking speed              | > 200 clones per hour   |
| Wash bath                  | Ethanol wash bath, automatically refilled   |
| Picking system fluids      | 5 L sterile water supply, 5L waste bottle   |
| Pin drying                 | Proprietary halogen pin drying station  |
| Instrument dimensions      | 1010 mm (width) x 900 mm (depth) x 1490 mm (height)   |
| Instrument weight          | 350 kg  |

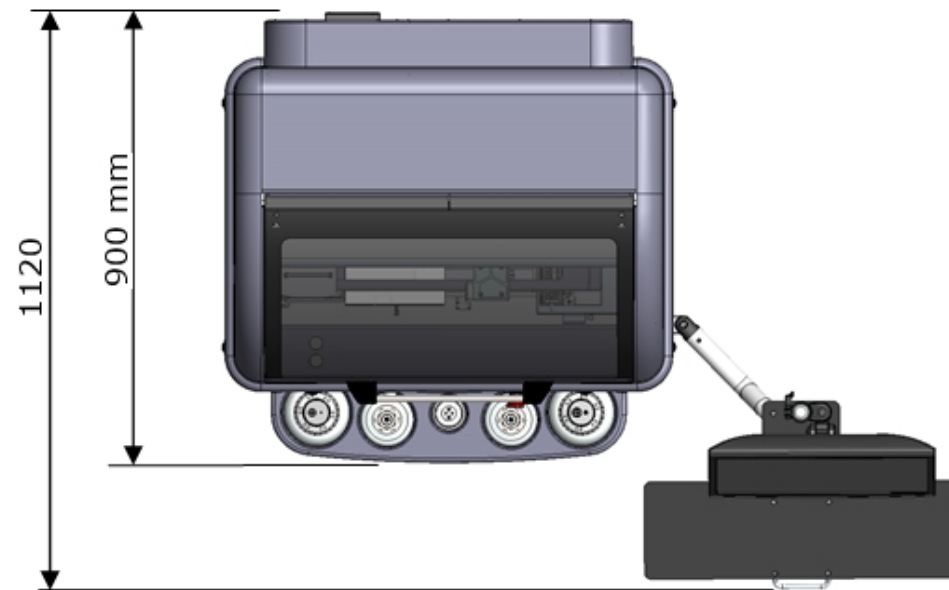
**ClonePix 2 Mammalian Colony Picker Technical Specifications (continued)**

| Item   | Description  |
|--|--|
| <b>Compressor air requirements</b>                               |  |
| Air  | Clean, oil-free with sub-micron filtration                             |
| Minimum operating pressure                                       | 6 bar (~90psi)   |
| Minimum operating volume   | 80 L/min   |
| <b>Optional air compressor</b>                                   |  |
| Compressor unit  | DynAir DA7001CS, Clean, oil-free compressor with sub-micron filtration |
| Dimensions   | 440 mm (width) x 440 mm (depth) x 613 mm (height)                      |
| Weight   | 84 lbs (38 kg)   |
| Power  | 0.75 KW  |
| Tank size  | 9 L  |
| Pump head number   | 2  |
| Rated output free air  | Max 5.37CFM/152L.Min   |
| Noise level  | 56 ±5 dB/A   |
| Drain  | Manual / buffer bottle   |
| <b>Regulatory approval</b>                                       |  |
| Compliance   | CE   |
| Quality  | ISO9001:2008 certified   |
| <b>Electrical Ratings</b>  |  |
| 100V   | 50/60 Hz, 1250W  |
| 115V   | 60 Hz, 1250W   |
| 230V   | 50 Hz, 1250W   |
| Factory set and changed by service personnel during installation |  |

### Instrument Dimensions



Front View of the ClonePix 2 Mammalian Colony Picker



Top View



# Appendix B: Plate Compatibility



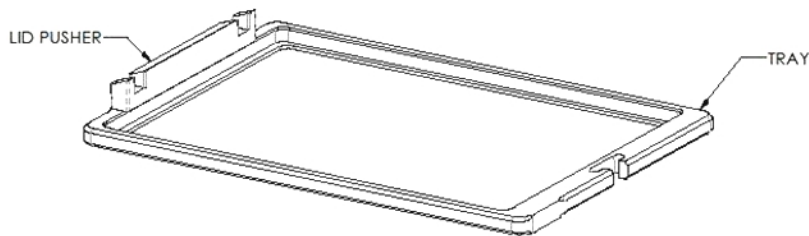
The ClonePix 2 stacker system uses plate holders and a lid lift system to transport the plate into the correct position ready for picking and depositing.

## Plate Holders

There are two types of plate holders that are designed for various plate types. The Lid Pusher at the back of the plate holder differs between the two types. This is an important part of the plate holder because it pushes the lid onto the lid lift.



**CAUTION!** If you use the wrong plate holder, the lid is not pushed onto the lid lift correctly and result in stacker failure.



### Plate Holders

| Part Number                  | Lid Pusher Color | Part Number |
|------------------------------|------------------|-------------|
| PLATE CARRIER ASSY (13.4 MM) | Green            | ME004233    |
| PLATE CARRIER ASSY (8.7MM)   | Blue             | ME004232    |

### Source Plates

| Part Number              | Lid Pusher Color | Lid Lifter Assembly |
|--------------------------|------------------|---------------------|
| Greiner 6 well - #657185 | Green            | Red                 |
| Nunc Omni - #242811      | Blue             | Silver              |
| Nunc – 6 Well -#140675   | Green            | White               |

**Destination Plates**

| Part Number                             | Lid Pusher Color | Lid Lifter Assembly |
|---|------------------|---------------------|
| Costar 96 FB - #3300                    | Blue             | Silver              |
| Costar 96 FB - #3595                    | Blue             | Silver              |
| Costar 96 FB - #3596                    | Blue             | Silver              |
| Costar 96 FB - #3598                    | Blue             | Silver              |
| Falcon 96 FB - #353072                  | Blue             | Silver              |
| Greiner 96 FB - #655185                 | Blue             | Silver              |
| Matrix 96 FB - #4942                    | Blue             | Silver              |
| Nunc 96 FB - #167008                    | Blue             | Silver              |
| Nunc 96 FB - #143761 (Lids from 163320) | Blue             | Silver              |

**Lid Lifter Assembly**

There are three different types of Lid Lifter Assembly used in the ClonePix 2 system. These assemblies are designed to remove the lids of specific plate types.



**CAUTION!** To avoid plate jams, always use the correct lid lifter assembly with the correct plate type.




# Appendix C: Replacement Parts, Accessories, and Consumables



## Replacement Parts

### ClonePix 2 Instrument Replacement Parts

| Part Number | Description  |
|-------------|--|
| X4941       | Additional picking head populated with eight X4961 (F1 Pins)<br> <b>Note:</b> The standard instrument comes with one picking head with F1 pins. |
| X4942       | Additional picking head populated with eight X4962 (F2 Pins)   |
| X4949       | Additional picking head without pins   |
| X4961       | F1 – ClonePix colony picking pin, 0.40 mm diameter (Single pin)  |
| X4962       | F2 – ClonePix colony picking pin, 0.70 mm diameter (Single pin)  |
| X1036       | O Rings for picking pins (Pack of 8)   |
| X4970       | Replacement picking tubing (Pack of 8)   |
| X4948       | Picking Pin Removal Key  |
| X4972       | 2L Bottle and Fittings (for either wash or waste)  |
| X4975F      | 5L Feed Bottle and Fittings (for wash)   |
| X4975W      | 5L Waste Bottle and Fittings   |
| X4976       | Nylon bristle ethanol bath insert  |

## Accessories and Consumables

### ClonePix 2 System Optional Extras

| Part Number | Description  |
|-------------|--|
| SL4950-A01  | ClonePix Remote Data Viewer Software   |
| X4990       | ClonePix filter set ex 440 em 505 (CFP)  |
| X4992       | ClonePix filter set ex 500 em 550 (YFP)  |
| X4993       | ClonePix filter set ex 530 em 590 (CloneDetect 549, Rhodamine)                         |
| X4992       | ClonePix filter set ex 622 em 700 (CloneDetect 622, Cy5)                               |
| SLLG025SS   | Millex-LG filter unit 0.20 µm hydrophilic 25 mm ethylene oxide sterilized              |
| K2505       | aQu Clean pin cleaning solution (1L)   |
| K8150       | CaliBeads: Fluorescent beads for ClonePix 2. 200 µm diameter. Pan-wavelength. (100 ml) |

## Reagents and Media

For an up-to-date list of reagents and media, see the website at: [www.moleculardevices.com](http://www.moleculardevices.com).



## Appendix D: Imaging Definitions



**Statistics for Prime Configurations**

| Statistic             | Description   | Category    | Unit            |
|-----------------------|---|-------------|-----------------|
| Actual X              | The X co-ordinate of the center of the feature in mm  | Position    | mm              |
| Actual Y              | The Y co-ordinate of the center of the feature in mm  | Position    | mm              |
| Area                  | The area covered by the feature (excluding 'child' features) in square mm   | Morphology  | mm <sup>2</sup> |
| Axis Ratio            | The ratio of the minimum and maximum radiuses of the feature measured from 0 (very elongated) to 1 (a perfect circle) | Morphology  |                 |
| Block                 | The identifier of the block of features this feature is associated with   | General     |                 |
| Compactness           | A measure of how compact the feature is, measured from 0 (not compact) to 1 (a perfect circle)                        | Morphology  |                 |
| Deposit Barcode       | Barcode of the plate the colony has been deposited in   | Information |                 |
| Deposit Well          | The well the colony has been deposited in   | Information |                 |
| Edge Excluded         | Whether the colony center lies within the exclusion zone  | Position    |                 |
| Feature ID            | A unique identifier   | General     |                 |
| Group                 | Group the feature is assigned to  | General     |                 |
| Image Column          | The column the image containing the feature is in   | Position    |                 |
| Image Row             | The row the image containing the feature is in  | Position    |                 |
| Intensity SD          | The standard deviation of the intensity of the pixels in the feature  | Intensity   |                 |
| Manual Group          | Whether the feature was manually added to its current group   | General     |                 |
| Mean Center Intensity | The mean intensity of the nine pixels at the center of the feature  | Intensity   |                 |
| Mean Intensity        | The mean intensity of all the pixels in the feature   | Intensity   |                 |
| Median Intensity      | The median intensity of all the pixels in the feature   | Intensity   |                 |
| Perimeter             | The length of the perimeter of the feature in mm  | Morphology  | mm              |
| Picked                | Flag to signify if the colony has been picked.  | Information |                 |
| Pixel Area            | The area covered by the feature (excluding 'child' features) in pixels  | Morphology  | pixels          |

**Statistics for Prime Configurations (continued)**

| Statistic            | Description  | Category   | Unit            |
|----------------------|--|------------|-----------------|
| Pixel Perimeter      | The length of the perimeter of the feature in pixels   | Morphology | pixels          |
| Pixel Radius Max     | The maximum radius of the feature in pixels  | Morphology | pixels          |
| Pixel Radius Min     | The minimum radius of the feature in pixels  | Morphology | pixels          |
| Pixel Radius SD      | The radius standard deviation of the feature in pixels   | Morphology | pixels          |
| Pixel Total Area     | The total area of the feature (including 'child' features) in pixels                           | Morphology | pixels          |
| Pixel X              | The X co-ordinate of the center of the feature in pixels relative to the top left of the image | Position   | pixels          |
| Pixel Y              | The Y co-ordinate of the center of the feature in pixels relative to the top left of the image | Position   | pixels          |
| Proximity            | The distance to the closest neighboring colony in the same image                               | Morphology | mm              |
| Radius Max           | The maximum radius of the feature in mm  | Morphology | mm              |
| Radius Min           | The minimum radius of the feature in mm  | Morphology | mm              |
| Radius SD            | The radius standard deviation of the feature in mm   | Morphology | mm              |
| Saturated Percentage | The percentage of saturated pixels in the feature  | Intensity  | %               |
| Saturated Pixels     | The number of saturated pixels in the feature  | Intensity  |                 |
| Selected             | Whether the feature is currently selected  | General    |                 |
| Source Barcode       | The barcode of the plate that the feature is in  | Position   |                 |
| Source Well          | The well the feature is in   | Position   |                 |
| Total Area           | The total area of the feature (including 'child' features) in square mm                        | Morphology | mm <sup>2</sup> |
| Volume Equivalent    | The volume of a sphere having the same cross-sectional area as the feature in cubic mm         | Morphology | mm <sup>3</sup> |
| Well Index           | A numerical annotation of the Source Well for graphical presentation                           | Position   |                 |

**Statistics for Fluorophore Configurations**

| Statistic                      | Description   | Category   | Unit            |
|--------------------------------|---|------------|-----------------|
| Exterior Area                  | The area of the pixels outside the feature boundary in the local vicinity of the feature                          | Morphology | mm <sup>2</sup> |
| Exterior Mean Intensity        | The arithmetic mean intensity of all the pixels outside the feature boundary in the local vicinity of the feature | Intensity  |                 |
| Exterior Geo Mean Intensity    | The geometric mean intensity of all the pixels in the feature   | Intensity  |                 |
| Exterior Median Intensity      | The median intensity of all the pixels outside the feature boundary in the local vicinity of the feature          | Intensity  |                 |
| Exterior Total Intensity       | The total intensity of all the pixels outside the feature boundary in the local vicinity of the feature           | Intensity  |                 |
| Interior Intensity SD          | The standard deviation of the intensity of the pixels in the feature  | Intensity  |                 |
| Interior Mean Center Intensity | The mean intensity of the nine pixels at the center of the feature  | Intensity  |                 |
| Interior Mean Intensity        | The arithmetic mean intensity of all the pixels in the feature  | Intensity  |                 |
| Interior Geo Mean Intensity    | The geometric mean intensity of all the pixels in the feature   | Intensity  |                 |
| Interior Median Intensity      | The median intensity of all the pixels in the feature   | Intensity  |                 |
| Interior Total Intensity       | The total intensity of all the pixels in the feature  | Intensity  |                 |
| Normalized Intensity           | The total intensity of the feature divided by the primary area  | Intensity  |                 |
| Sum Total Intensity            | The sum total intensity of the feature (interior and exterior)  | Intensity  |                 |



## Appendix E: Guide to Fluorophore Configurations



The ClonePix 2 system acquires images of fluorescing colonies and generates a range of statistics from the images in order that the colonies can be screened to select the most 'fit'. The user can choose which parameters to use to meet their definition of 'fit' – the software does not enforce any particular definition of what the 'best' colonies are.

Most users want to select either the highest expressing colonies or those that are positive for a specific parameter. The software generates various statistics to facilitate this. This appendix explains how these statistics are generated and provides advice on how they might be used.

### Measuring Fluorescence

The ClonePix 2 software works on the principal that all colonies can be detected under white light. All the detection algorithms in the software process the white light (Prime Probe) images to detect colony boundaries.

There are two principal advantages to this approach:

- Colonies that are not fluorescing can still be detected and considered when deciding which colonies to select.
- Knowing the boundaries of all colonies allows for a more rigorous approach to processing the fluorescent images.

Fluorescent colonies can be divided into two types:

- Those whose fluorescence is purely 'internal', that is, within the colony boundary as determined from the white light image.
- Those whose fluorescence is also present outside the colony boundary.

Generating statistics for the first type of fluorescent colony is straightforward because for any given pixel in the image the pixel is either within the boundary of one colony or it is a background pixel and can be ignored. The only complication is determining what proportion of the intensity of the pixel is 'real' fluorescence and what proportion is 'background' fluorescence. Background fluorescence would exist even in the absence of any colonies since fluorescence images rarely have a background of zero – pure black.

Generating statistics for the second type of fluorescent colony (those whose fluorescence is also present outside the colony boundary) is more challenging. This is because when two or more of these colonies are close together, their fluorescence mingles together and there is no deterministic mechanism to accurately decide what proportion of intensity of a given pixel can be attributed to each colony. The problem is compounded as the density of colonies increases (that is, when there are more colonies in each plate). Also, fluorescence from the second type of colony can affect the first type of colony, which can therefore no longer be assumed to be independent of other colonies.

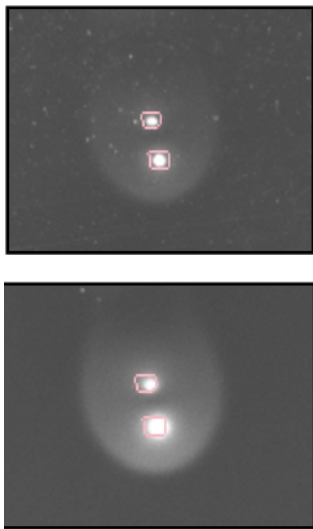
Because of this difficulty, the ClonePix 2 software does not attempt to determine what proportion of intensity of pixels outside of colony boundaries comes from any particular colony. Instead, all the pixels surrounding each colony are included. This results in some pixels being included in the total for several colonies, but this is not a shortcoming.

It is important to know that the ClonePix 2 system measures relative fluorescence, not absolute fluorescence. This means that fluorescent statistics can only meaningfully be compared when generated from the same plate or batch of identical plates at the same time. It is not necessarily meaningful to compare fluorescent statistics between different plates imaged on different days. This is because several things can vary that affect fluorescence, for example exposure time. Varying the exposure time used when acquiring fluorescent images increases or decreases the absolute values of the fluorescent statistics. However, all colonies in general get brighter or darker so their brightness relative to each other is unaffected.

## Example of Fluorescent Statistic Generation

The white light image (top) shows two colonies close together. Their approximate boundaries, as determined by the colony detection algorithm are shown in pink.

The corresponding fluorescent image is shown (bottom) with the colony detection algorithm in pink overlaid. Several fluorescent statistics can be generated from this directly.



**Note:** A background subtraction algorithm is applied to all fluorescent images before intensity statistics are generated. This algorithm seeks to determine what level of fluorescence there would be in the absence of any colonies, and removes this from the image so that the intensity statistics include only 'real' fluorescence.

## Interior Intensity Statistics

The simplest statistic is **Interior Total Intensity**. This is the sum of the intensities of all the pixels of the fluorescent image within the colony boundary from the matching white light image. Larger colonies have larger totals because they are larger. This statistic is not therefore especially useful on its own but several other statistics are derived from it.

**Interior Mean Intensity** is Interior Total Intensity divided by colony area (in pixels). It is the 'average' brightness of all pixels within the colony, or the arithmetic mean. Since it does take account of area, this statistic can be used to compare colonies with different areas. In fact, this is probably the most powerful statistic (along with Interior Median Intensity) since colonies with a high Interior Mean Intensity are overwhelmingly likely to be bright colonies – rather than faint colonies that happen to be near a bright colony that is dispersing fluorescence into the surrounding medium.



**Interior Geo Mean Intensity** is similar to the Interior Mean Intensity but is calculated as the geometric mean. This statistic can never be greater than the Interior Mean Intensity, since the way it is calculated tends to reduce the influence of the largest values. Whether the geometric mean is a 'better' statistic to use than the arithmetic mean depends entirely on the colonies in question.

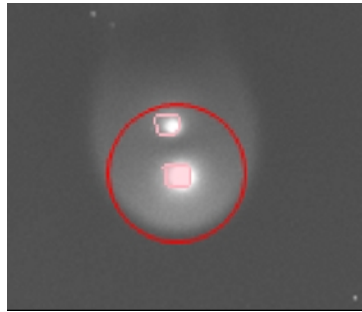
**Interior Intensity SD (Standard Deviation)** is a measure of how the intensity of pixels within a colony boundary varies. Colonies with a low value for this statistic must have pixels that all have a similar brightness. Conversely, those with high values have both bright pixels and faint pixels with their colony boundaries. This statistic can be used to differentiate between colonies that have similar Interior Mean Intensity Values.

**Interior Median Intensity** is the intensity of the pixel midway between the faintest pixel and the brightest pixel within the colony boundary. This is similar to Interior Mean Intensity, but tends to be more robust since its value is less affected by a few bright or a few faint pixels. As with Interior Mean Intensity, colonies with a high Interior Median Intensity are likely to be bright colonies.

**Interior Mean Center Intensity** is the mean intensity of the nine pixels closest to the center of the colony (that is, the center pixel and its immediate eight neighboring pixels). Often, colonies are brightest at their center so this statistic provides an unbiased measure of the brightness of a colony.

## Exterior Intensity Statistics Calculation

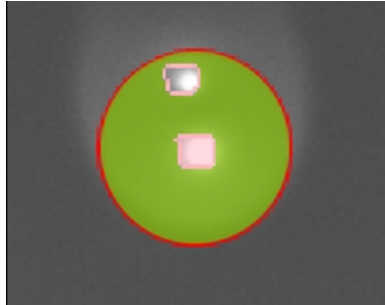
The fluorescent image is shown again with the lower colony highlighted (its interior is shaded pink). Around the colony is a red circle, which is visualized by moving the mouse pointer over the colony. The diameter of the red circle is derived from a combination of the Average Colony Diameter and the Exterior Statistics Diameter Multiplier provided by the user in the Detection tab when the image was originally processed.



In this example, the user had selected an average colony diameter of 0.4 mm and an Exterior Statistics Diameter Multiplier of x3 (left). Therefore the red circle has a diameter of 1.2 mm. It is important to appreciate that the size of this circle is the same for every colony in the image even though the colonies themselves are different sizes.

The significance of the red circle is that it defines the boundary within which pixels for a given colony are processed to generate the Exterior Intensity Statistics.

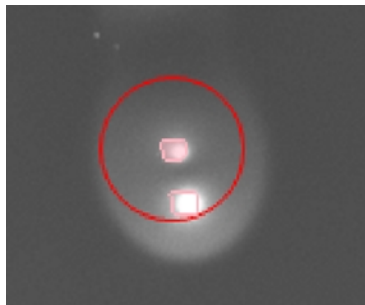
Exterior Intensity Statistics are generated by first considering all the pixels within the red circle. Following this, any pixels within boundary if the colony of interest are excluded (the pixels shaded in pink). Next, all the pixels that lie within any other colonies are excluded. So in the image on the right, all the pixels within the boundary of the second unshaded colony would be excluded. This is because any fluorescence within the boundary of a neighboring colony is likely to have come from the neighboring colony.



What remains are the pixels that are local to the colony of interest but not within any other colony boundary (highlighted in green in the image). These are the pixels that are used to generate the 'exterior intensity' statistics.

If the colony is larger than the Average Colony Diameter specified by the user or if the colony is an irregular shape, it is possible that some of the pixels within the colony boundary would lie outside the red circle. In those circumstances, the red circle is not a good approximation of the area 'local' to the colony. Also it is not clear how to meaningfully compare irregular shaped colonies with the more circular shaped colonies because it is unlikely that the 'local' area values produced can be used for comparative purposes.

Neighboring colonies are treated in exactly the same way. The top colony (shaded in pink) has been selected. Again, the red circle overlaps the lower colony. This means that many pixels are used to generate exterior statistics for both colonies.



In this example, both colonies are fluorescing and dispersing their fluorescence into the local medium so, for the pixels close to both colonies, it is not clear what proportion of the intensity of each pixel is derived from which colony. The area of fluorescence around the colonies looks uniform implying that the fluorescence from each colony is not cumulative. This suggests that counting the intensities of the pixels twice (once from each colony) would be a reasonable approach.

## Exterior Intensity Statistics

**Exterior Area** (mm<sup>2</sup>) is the area of the 'exterior' pixels (those shaded green). It could be used to exclude colonies that are in crowded regions of the plate. A low value for a given colony compared to other colonies in the image indicate that the exterior intensity statistics for that colony might not be accurate. It is not in any way a measure of the actual size of the halo of fluorescence for any colony.

**Exterior Total Intensity** is the sum of the intensities of all the 'exterior' pixels. If there are neighboring colonies within the red circle, this value is lower. On its own, this statistic is not very useful since it cannot be used comparatively with other colonies.

**Exterior Mean Intensity** is the Exterior Total Intensity divided by the number of pixels in the exterior area (arithmetic mean). This statistic provides a good approximation of how bright the area immediately surrounding a given colony is, though there is no implication that the brightness is coming from the colony as opposed to one or more of its neighbors.

**Exterior Geo Mean Intensity** is similar to the Exterior Mean Intensity but is the geometric mean. See the explanation of Interior Geo Mean Intensity for further details.

**Exterior Median Intensity** is the intensity of the pixel midway between the brightest and the faintest pixel within the exterior pixels. Just as the Interior Median Intensity is more robust than the Interior Mean Intensity, so the Exterior Median Intensity is more robust than Exterior Mean Intensity.

**Sum Total Intensity** seeks to approximate the sum of the Exterior Total Intensity and Interior Total Intensity if there were no neighboring colonies within the red circle. Instead of using the actual Exterior Total Intensity it uses the Exterior Mean Intensity multiplied by the number of pixels within the red circle. The calculation also excludes the pixels within the boundary of the colony of interest but includes any pixels in neighboring colonies in the red circle. If this calculation were not taken into account, it would not be meaningful to compare Sum Total Intensity between colonies because the value would vary depending on the presence of neighboring colonies.

**Normalized Intensity** is Sum Total Intensity divided by colony area (in pixels). The highest secreting colonies have the highest values here but unfortunately in practice normalizing by area biases the statistic in favor of small colonies rather than bright colonies. This is because the interior intensity is always much brighter than the exterior intensity which decreases exponentially from the colony boundary.



## Appendix F: ClonePix 2 Data Viewer Software



**Note:** The ClonePix 2 Data Viewer software is optional. Contact Molecular Devices Technical Support to purchase.

The ClonePix 2 Data Viewer software enables you to view results generated by the ClonePix 2 instrument on a computer that is remote from the instrument.

You can either access a copy of the results files or you can access the folder on the instrument computer using Windows folder-sharing if the two computers are networked together. The default folder location is C:\Image archive.



**Tip:** Any time you install and use the ClonePix 2 Data Viewer software on a computer that is detached from the instrument, it functions as a data viewer.

### Software Prerequisites

The ClonePix 2 Data Viewer software requires the installation of Microsoft .NET™ Framework 4.5.2 or newer. Your computer might already have it installed.

To check for .NET Framework 4.5.2:

1. Go to the Control Panel and click **Add/Remove Programs**.
2. Look in the list for **Microsoft.Net 4.5.2** or newer.



**Tip:** Windows 10 computers should already have Microsoft .Net 4.6 or newer installed.

3. If Microsoft .Net 4.6 or newer is installed, continue to the next step, otherwise install it now.

### Installing ClonePix 2 Data Viewer Software

After verifying that the .NET Framework 3.5 SP1 is installed, you can install the ClonePix 2 Data Viewer software. See [Software Prerequisites](#).

To install the software:

1. Log in to the Molecular Devices Support website, go to the [ClonePix™ System Software Download Page](#), and in the table, click the **DataViewer > Installer** link.
2. Double-click ClonePix 2 **Data Viewer v.x.x.x**, where x.x.x is the software version number, to start the installation.  
To start the installation, if you are using the *Recovery DVD*, double-click **Setup\_ClonePix2\_Viewer.exe**.
3. Follow the on-screen instructions.

### Licensing ClonePix 2 Data Viewer Software

When the ClonePix 2 Data Viewer software is installed, you must license it.

The first time that you start the ClonePix 2 Data Viewer software after completing the installation, the software prompts you for a license.

## Using ClonePix 2 Data Viewer Software

When the ClonePix 2 Data Viewer software starts, the main process screen appears.

To view the results, locate where they are stored. On the right side of the screen, click **Results Location** and browse for the location of the results. The location can be a local folder on the computer, a CD-ROM or DVD, or, if the two computers are networked together, it can be a shared folder on the instrument computer.

After you select the location, the software catalogs all of the results and displays a list. You can then view the results the same way as you do on the instrument computer.

## Contact Us

Phone: [+1-800-635-5577](tel:+18006355577)  
Web: [moleculardevices.com](http://moleculardevices.com)  
Email: [info@moldev.com](mailto:info@moldev.com)

Visit our website for a current listing of worldwide distributors.