



# IonWorks Barracuda<sup>®</sup>

Automated Patch Clamp System

## Software Release Notes

Version 2.6  
August 2016



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 perform 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.

Product manufactured by Molecular Devices, LLC.

1311 Orleans Drive, Sunnyvale, California, United States of America 94089.

Molecular Devices, LLC is ISO 9001 registered.

© 2016 Molecular Devices, LLC.

All rights reserved.

# Contents

<b>Chapter 1: IonWorks Barracuda® Software General information</b>	<b>5</b>
System Requirements . . . . .	5
Installing IonWorks Barracuda Software . . . . .	6
Obtaining Support . . . . .	7
<b>Chapter 2: IonWorks Barracuda Software Version 2.6 Software Release Notes</b>	<b>9</b>
New in IonWorks Barracuda Software Version 2.6 . . . . .	9
<b>Chapter 3: IonWorks Barracuda Software Version 2.5.4 Software Release Notes</b>	<b>11</b>
Resolved Issues . . . . .	11
<b>Chapter 4: IonWorks Barracuda Software Version 2.5.3 Software Release Notes</b>	<b>13</b>
Modified . . . . .	13
Known Issues . . . . .	13
<b>Chapter 5: IonWorks Barracuda Software Version 2.5.2 Software Release Notes</b>	<b>15</b>
New in IonWorks Barracuda Software Version 2.5.2 . . . . .	15
Modified . . . . .	15
Resolved Issues . . . . .	16
<b>Chapter 6: IonWorks Barracuda Software Version 2.5.1 Software Release Notes</b>	<b>17</b>
New in IonWorks Barracuda Software Version 2.5.1 . . . . .	17
<b>Chapter 7: IonWorks Barracuda Software Version 2.5 Software Release Notes</b>	<b>19</b>
New in IonWorks Barracuda Software Version 2.5 . . . . .	19
Modified . . . . .	20
Resolved Issues . . . . .	21

Known Issues . . . . .	21
<b>Chapter 8: IonWorks Barracuda Software Version 2.3 Software Release Notes . . . . .</b>	<b>23</b>
New in IonWorks Barracuda Software Version 2.3 . . . . .	23
Modified . . . . .	23
Resolved Issues . . . . .	24
Known Issues . . . . .	25
<b>Chapter 9: IonWorks Barracuda Software Version 2.1.1 Software Release Notes . . . . .</b>	<b>27</b>
New in IonWorks Barracuda Software Version 2.1.1 . . . . .	27
Modified . . . . .	27
<b>Chapter 10: IonWorks Barracuda Software Version 2.1 Software Release Notes . . . . .</b>	<b>29</b>
New in IonWorks Barracuda Software Version 2.1 . . . . .	29
Modified . . . . .	29
Resolved Issues . . . . .	30
Known Issues . . . . .	31
<b>Chapter 11: IonWorks Barracuda Software Version 2.0 Software Release Notes . . . . .</b>	<b>33</b>
New in IonWorks Barracuda Software Version 2.0 . . . . .	33
Resolved Issues . . . . .	35
Known Issues . . . . .	37

## information

IonWorks Barracuda® Software controls the functions of the IonWorks Barracuda instrument, displays acquired data, and provides data analysis and export functions after an assay completes. The software comes pre-installed on new system computers.

The following topics are included in this chapter:

- [System Requirements on page 5](#)
- [Installing IonWorks Barracuda Software on page 6](#)
- [Obtaining Support on page 7](#)

## System Requirements

IonWorks Barracuda Software can be installed on additional computers to work in Analysis only mode. Before installing the software, make sure the computer meets the following minimum system requirements:

- CPU: Quad core 2.4 GHz
- RAM: 8 GB
- OS: Windows 7 (64-bit) and Windows 10 (64-bit)
- Graphics card supporting a resolution of 1920 x 1200



---

**Note:** The software is designed for a 24-inch monitor with a display resolution of 1920 x 1200, and default font setting. If you are installing on a computer with a lower resolution or screen size, or change the computer font size to larger than the 100% (default) setting, you might experience minor display issues.

---

## Installing IonWorks Barracuda Software

1. Double-click the **BarracudaInstaller\_2.6.x.exe** IonWorks Barracuda Software installation file. A Welcome to the IonWorks Barracuda Software Setup Wizard dialog is displayed.
2. When updating your software, you must first uninstall the current IonWorks Barracuda Software. When prompted, follow the uninstall wizard instructions.
3. Again, double-click the **BarracudaInstaller\_2.6.x.exe** IonWorks Barracuda Software installation file. A Welcome to the IonWorks Barracuda Software Setup Wizard dialog is displayed.
4. Click **Next**.
5. In the **License Agreement** dialog box, select **I accept the terms of the license agreement**, and click **Next**.
6. In the **Setup Type** dialog, designate the default mode in which you want the software to start.
  - ◆ In **Instrument**, IonWorks Barracuda Software connects to the IonWorks Barracuda instrument when the software is started. Use this mode for software installed on the system computer.
  - ◆ In **Simulator**, IonWorks Barracuda Software is not connected to the IonWorks Barracuda instrument. Use this mode for software installed on an analysis computer.
7. Click **Next**.
8. In the **Destination Folder** dialog, the **Install IonWorks Barracuda Software to** field displays the default installation directory. To change the installation directory, click **Change**, navigate to the desired directory, then click **OK**.
9. Click **Next**.
10. In the **Select Program Folder** dialog, leave the displayed default **Program Folder** settings. Select **Anyone who uses this computer** to make IonWorks Barracuda Software available to all users on the IonWorks Barracuda Automated Patch Clamp System computer, then click **Next**.
11. The **Completing the setup for IonWorks Barracuda** dialog appears. Click **Next**. The software installation begins.



---

**Note:** If an **Error** dialog appears, click **OK**, then click **Finish**. Repeat steps 1 through 9. At step 10, select **Only for me (current user)**, then click **Next**. Repeat step 11.

---

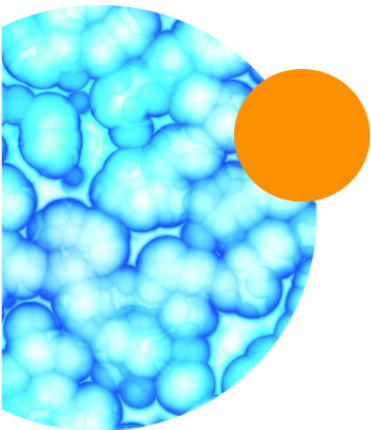
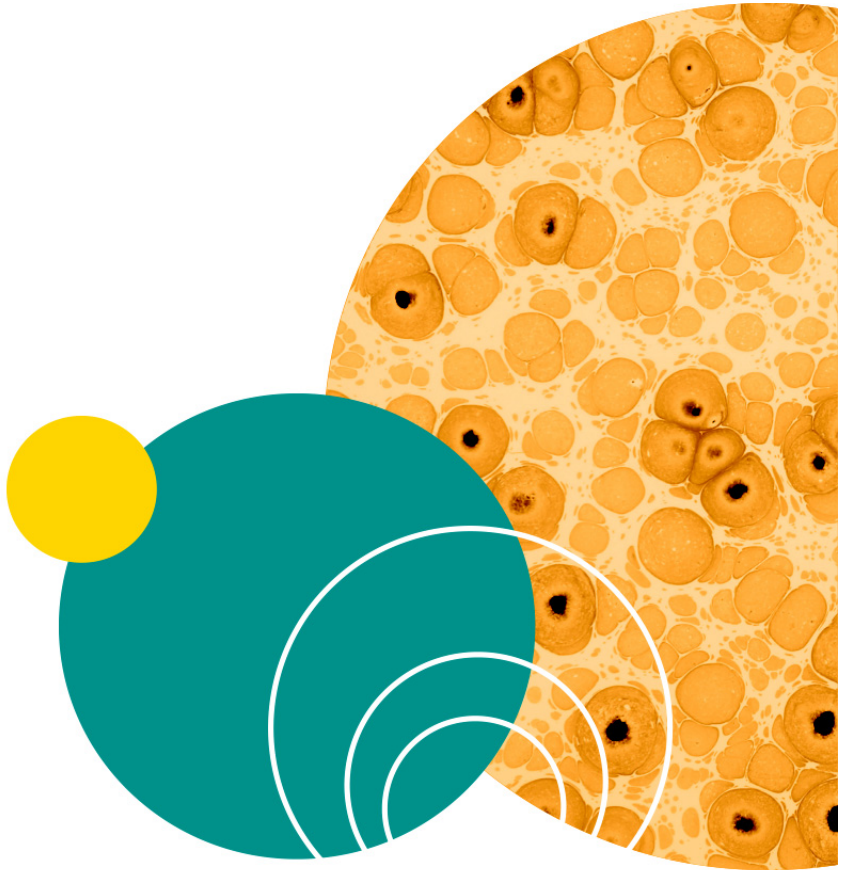
12. Click **Finish** to exit the wizard.

## Obtaining Support

Molecular Devices is a leading worldwide manufacturer and distributor of analytical instrumentation. We are committed to the quality of our products and to fully supporting our customers with the highest possible level of technical service.

Our support web site, [www.moleculardevices.com/support.html](http://www.moleculardevices.com/support.html), has a link to the Knowledge base with technical notes, software upgrades, and other resources. If you do not find the answers you are seeking, follow the links to the Technical Support Service Request Form to send an email message to a pool of technical support representatives.

You can contact your local representative or contact Molecular Devices Technical Support by telephone at 800-635-5577 (U.S. only) or +1 408-747-1700. In Europe call +44 (0) 118 944 8000. Please have the system ID number, system serial number, software version number, and the system owner's name available when you call.





# IonWorks Barracuda Software Version 2.6 Software Release Notes

## 2

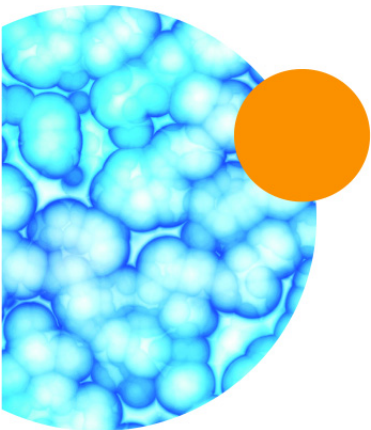
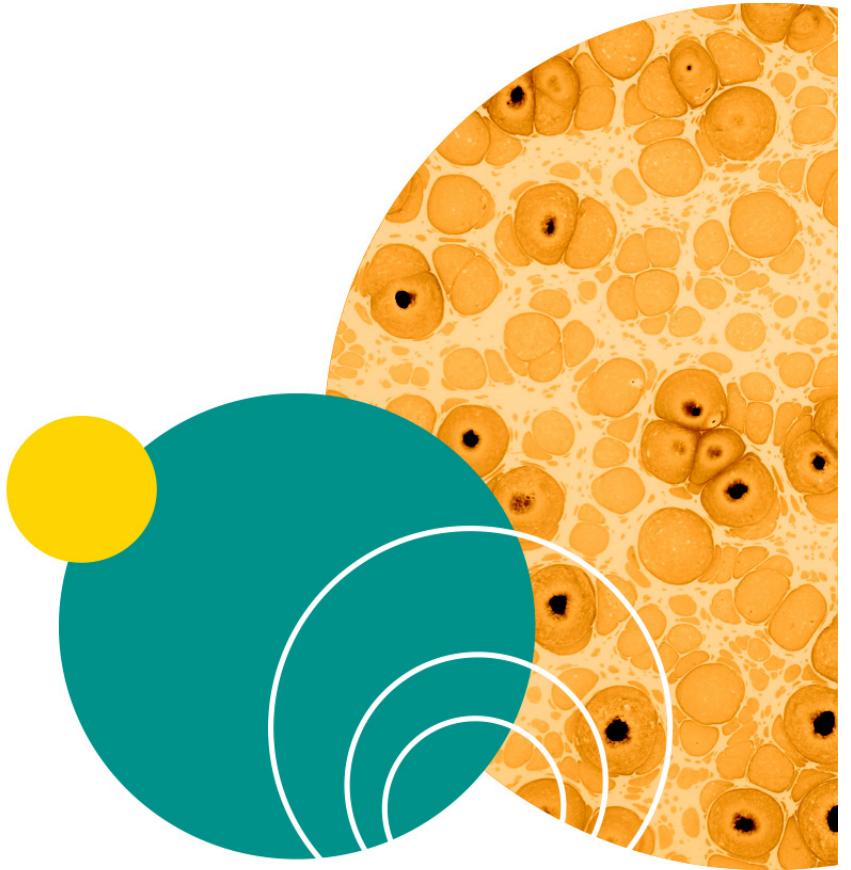
This chapter provides release notes for version 2.6 of the IonWorks Barracuda® Software. Read this entire document before you use the software. The following is a summary of the changes incorporated in this update as compared to version 2.5.4, the last general release of the IonWorks Barracuda® Software.

- [New in IonWorks Barracuda Software Version 2.6](#)

## New in IonWorks Barracuda Software Version 2.6

### Windows 10 Operating System Support

Adds software support for the Windows 10 (64-bit) operating system.



# IonWorks Barracuda Software Version 2.5.4 Software Release Notes

## 3

This chapter provides release notes for version 2.5.4 of the IonWorks Barracuda® Software. Read this entire document before you use the software. The following is a summary of the changes incorporated in this update as compared to version 2.5.3, the last general release of the IonWorks Barracuda® Software.

- [Resolved Issues](#)

## Resolved Issues

### Data File Naming Error In Multiple Protocol Mode

Tracking ID: FB3396

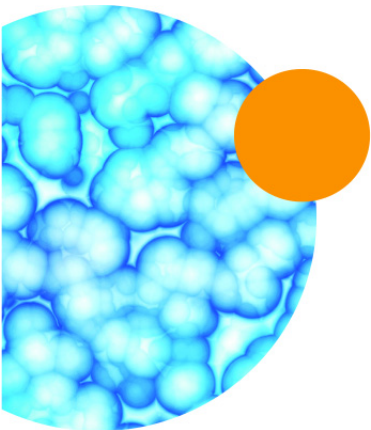
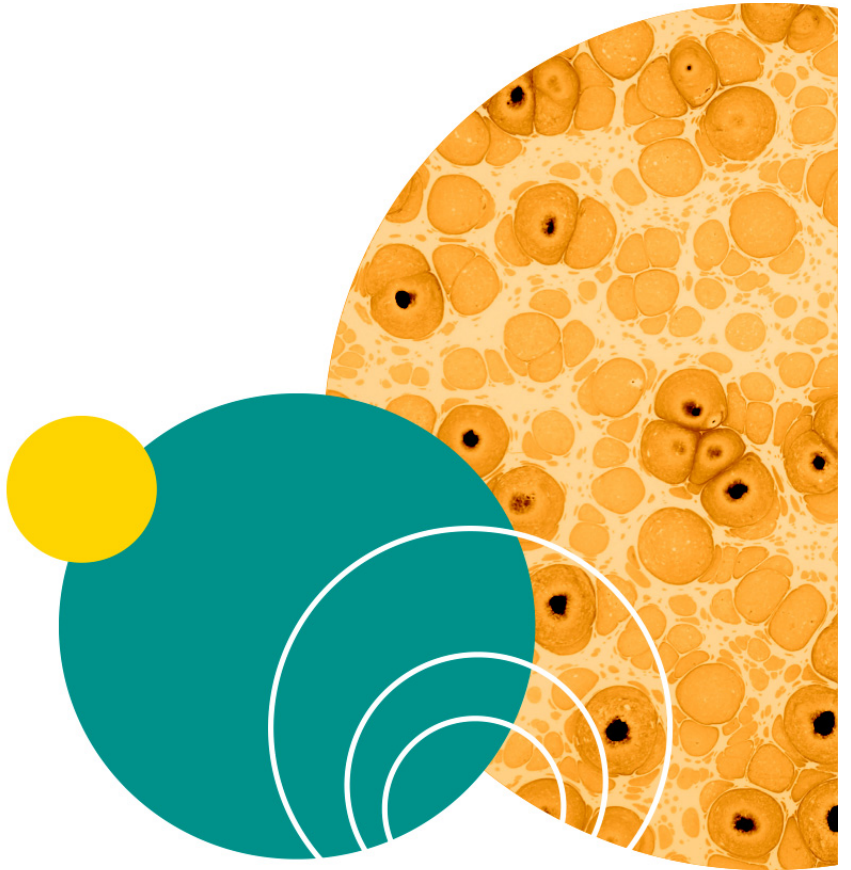
In Multiple Protocol Mode, the data files sometimes named as files from the previous experiment run.

**Resolution:**

The issue is now resolved in the software.

**Impact of the fix:**

This fix has no impact on data.



# IonWorks Barracuda Software Version 2.5.3 Software Release Notes

## 4

This chapter provides release notes for version 2.5.3 of the IonWorks Barracuda® Software. Read this entire document before you use the software. The following is a summary of the changes incorporated in this update as compared to version 2.5.2, the last general release of the IonWorks Barracuda® Software.

- [Modified](#)
- [Known Issues](#)

### Modified

#### Revised Help File and User Guide

The Help file and user guide have been revised to include supporting information and usage procedures for the new Replaceable Ground Electrodes (REGEs) and supporting plenum.

### Known Issues

#### Export Metric Dialog Should Close when OK Clicked

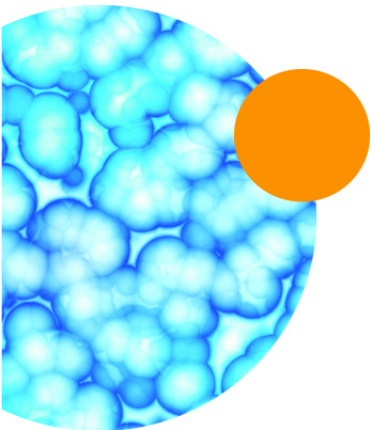
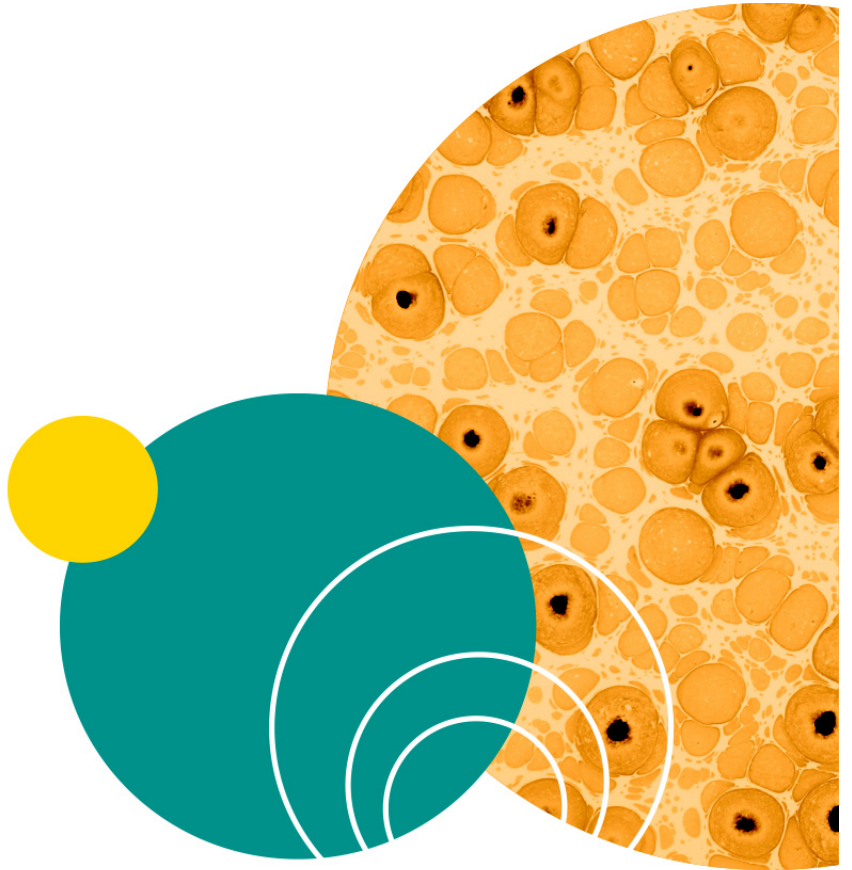
Defect ID:15136

##### Impact

When performing the **Export Metric** procedure **Edit > Channel Protocol Editor > Specify Metric Parameters > Configure Metric Export > OK**, the **Export Metric** dialog stay open when it should close.

##### Workaround

Click the **X** in the top right corner to close the dialog manually. Your changes save during the closing.



# IonWorks Barracuda Software Version 2.5.2 Software Release Notes

## 5

This chapter provides release notes for version 2.5.2 of the IonWorks Barracuda® Software. Read this entire document before you use the software. The following is a summary of the changes incorporated in this update as compared to version 2.5.1, the last general release of the IonWorks Barracuda® Software.

- [New in IonWorks Barracuda Software Version 2.5.2](#)
- [Modified](#)
- [Resolved Issues](#)

## New in IonWorks Barracuda Software Version 2.5.2

### RA1 Electrode-Plate Support and Usage Tracking

The PatchPlate Station supports the new RA1 Electrode-Plate. The **Electrode-Plate Type** field is now available in the **Add Experiment Notes** dialog for manually tracking which electrode-plate gets used in your experiments.

### Low Profile Reservoir Plate Type Support

The Seahorse Low Profile Reservoir replaces the deep buffer boat reservoir. The Seahorse Low Profile Reservoir option is available in the **Edit Channel Protocol > Plate Type** field.

## Modified

### Expanded Maintenance and Troubleshooting Information

The Maintenance and Troubleshooting information in the software-based Help file and the PDF file User Guide is expanded to include more hardware cleaning and user-based testing procedures.

## Resolved Issues

### Interrupted Amphotericin Filling

Defect ID: 12777

The software checks the status of every sub-system before proceeding to the next action to prevent the occasional occurrence of the data acquisition starting before the amphotericin filling completes.



# IonWorks Barracuda Software Version 2.5.1 Software Release Notes

## 6

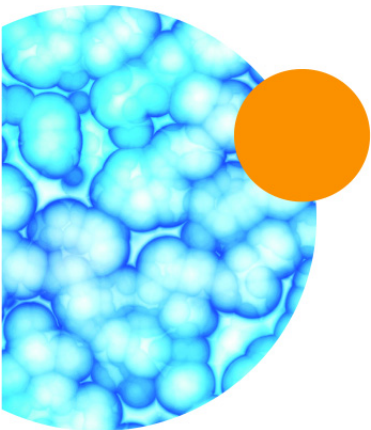
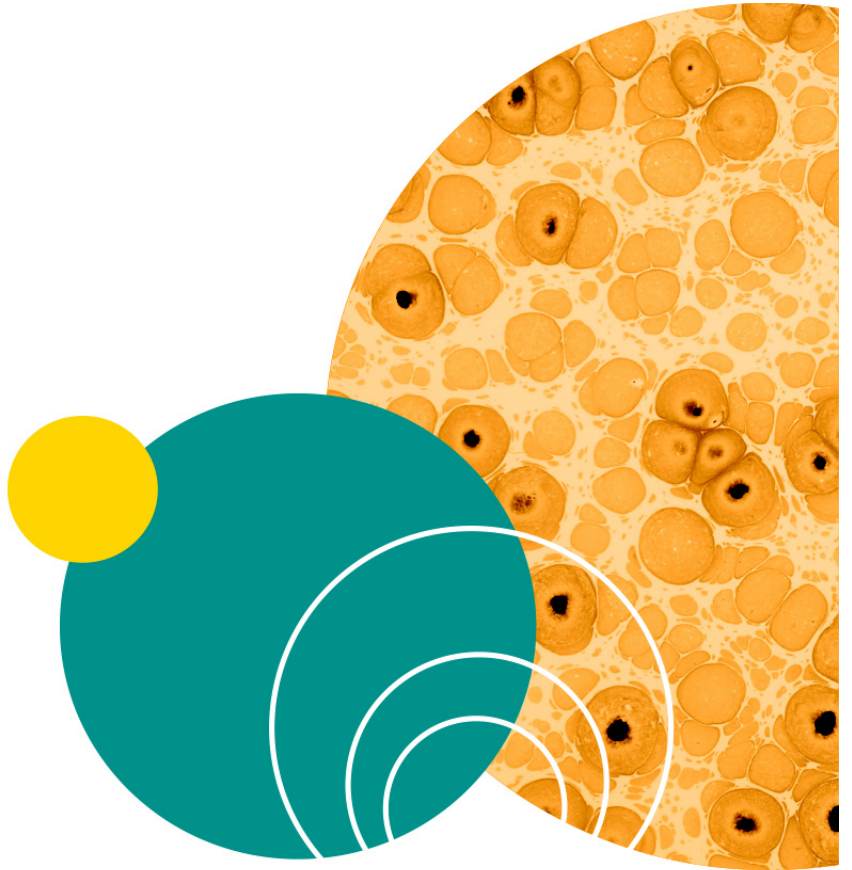
This chapter provides release notes for version 2.5.1 of the IonWorks Barracuda® Software. Read this entire document before you use the software. The following is a summary of the changes incorporated in this update as compared to version 2.5, the last general release of the IonWorks Barracuda® Software.

- [New in IonWorks Barracuda Software Version 2.5.1](#)

## New in IonWorks Barracuda Software Version 2.5.1

### Flush Tip Washer Lines Option

The **Utilities > Flush Tip Washer Lines** functionality is added as an instrument maintenance process. A 2% bleach solution is run through the lines to clean out growths and debris that can develop in the tubes, which cause blockages. Molecular Devices recommends running this process once every three months.



# IonWorks Barracuda Software Version 2.5 Software Release Notes

## 7

This chapter provides release notes for version 2.5 of the IonWorks Barracuda® Software. Read this entire document before you use the software. The following is a summary of the changes incorporated in this update as compared to version 2.3, the last general release of the IonWorks Barracuda® Software.

- [New in IonWorks Barracuda Software Version 2.5](#)
- [Modified](#)
- [Known Issues](#)

## New in IonWorks Barracuda Software Version 2.5

### Prime Pipettor Tips Option

In order to reduce the risk of dispensing air bubbles when you reuse pipettor tips, you can now select the **Prime Pipettor Tips** option to run at the beginning of your Setup Protocol.

### Create, Edit and Run Multiple Channel Protocols

Selecting **Experiment Mode > Multiple Protocol Mode** allows you to save a sequence of channel protocols as a Multiple Channel Protocol. This ability is available in the Acquisition as well as the Analysis Only instances of the software.

### Barcode Position Settings

When barcodes are used in experiments, the position of the barcode on a compound plate must now be specified as part of each compound definition in the **Edit Channel Protocol** dialog using the new **Edit Channel Protocol > Compound Addition > Aspirate > Barcode Position** field.

If the selected channel protocol has a barcode position set to anything other than NONE for one or more of the compounds, the software prompts to scan barcodes at the beginning of the experiment.

## Export Barcode Information With Trace Data Export

If used in an experiment, the barcode information is now also exported with Trace Data exports using the **Advanced Export Traces > Include header** option. When barcodes are not included in an experiment, placeholder text is added in the header.



---

**Note:** The additional exported barcode information rows and columns can impact your existing Excel macros that may have been created to read the data from these previously exported files. Adjustments may be required.

---

## End of Day Flush and Rinse Monitoring

To maintain the system properly, you must run End of Day Flush and Rinse within six days of system usage. Otherwise, on the seventh day, the system stops running experiments until you run Utilities > Plenum > End of Day Flush and Rinse.

## Modified

### Include Barcode in Experiment Name

The **Set > Set Experiment Name > Include Barcode** option now only controls the inclusion of the barcode information in the experiment name and is tied to the Channel Protocol Barcode Position field setting. The barcode information only appears in your experiment name when the **dit Channel Protocol > Compound Addition > Aspirate > Barcode Position** field is set to anything other than NONE.

### Continue Experiment Without Barcode Entry

When prompted to enter a barcode, you can now skip the barcode number entry, and run your experiment.

### Add Perforation Agent Duration Change

The fixed duration for **Edit Setup Protocol > Obtain Cell Access > Add Perforation Agent** to complete is now 115 seconds.

## Resolved Issues

### Traces for Scans with Multiple Sweeps are Mapped Incorrectly

Defect ID:10218

Exported trace data switches the Sweep and Well numbers in the file header. For example, data from Sweep1 Well1, Sweep1 Well2 display as Sweep1 Well1, Sweep2 Well1.

### Export Metrics Column Label has wrong sweep numbers

Defect ID:10793

Exported metrics files display incorrect column headers. For example, Scan 7 Sweep10 Metric1 sometimes appears as S7\_SW1\_M1 instead of S7\_SW10\_M1.

### Post Compound Scans Export with Incorrect Names

Defect ID:9534

When multiple post compound scans are selected, not all of the traces are exported as expected.

## Known Issues

### Metrics Export Incorrectly When a Multiple Channel Protocol with More Than one Protocol is set up for Automatic Export

Defect ID:12714

#### Impact

When assays are performed in Multiple Protocol Mode or in Assay Development Mode using Automated Metrics Export, only the metrics for the last channel protocol of the experiment exports automatically.

#### Workaround

Review data and export metrics from any of the channel protocols in offline mode.

## Multiple Channel Protocol Mode does not let users scan barcodes for all channel protocols

Defect ID:12775

### Impact

When running experiments in Multiple Protocol Mode or in Assay Development Mode, only the first channel protocol requiring barcode entry is handled, all subsequent channel protocols requiring barcode entry are ignored.

### Workaround

None.

### Planned Solution

The issue will be fixed in a future version of the software.

## Occasionally the Data Acquisition Starts Before the Amphotericin Filling Completes

Defect ID:12777

### Impact

Under extremely rare circumstances, it has been reported that the data acquisition starts before the amphotericin filling is complete.

### Workaround

Insert a wait duration of 115 seconds or more following the Fill Perforation Agent operation to avoid the risk of acquiring data while the pumps are running.

# IonWorks Barracuda Software Version 2.3 Software Release Notes

## 8

This chapter provides release notes for version 2.3 of the IonWorks Barracuda® Software. Read this entire document before you use the software. The following is a summary of the changes incorporated in this update as compared to version 2.1.1, the last general release of the IonWorks Barracuda® Software.

- [New in IonWorks Barracuda Software Version 2.3](#)
- [Modified](#)
- [Known Issues](#)

## New in IonWorks Barracuda Software Version 2.3

### New Preferences > General > Protocol Editor

The **Protocol Editor > Enable Charge Calculator** check box enables and disables the **Coulomb Meter**, and **Estimated Seal Resistance** field in the **Edit Setup Protocol** dialog and the **Edit Channel Protocol** dialog.

### Estimate Experiment Total Charge

The ability to calculate and display the estimated charge for the experiment is available in the **Edit Setup Protocol** dialog and **Edit Channel Protocol** dialog. Use the new **Estimated Seal Resistance** field and **Coulomb Meter Index** display.

## Modified

### Edit Setup Protocol Dialog

Optional **Wash before cell addition** in **Add Cells** section and **Wash after cell addition** in **Seal Test** section functionality added. Used for top and bottom Plenum buffer washes through external and internal buffer exchanges before cell addition, and external and amphotericin exchanges after cell addition.

## Resolved Issues

### Software Crashes When Some Experiments Run in the Same Order Multiple Times

Defect ID:9060

An identified error will be logged in the Plate.log.txt file. The data files will be flagged and the experiment will continue. When errors are logged and data files are flagged, a warning will display in the **Experiment Name** section. For additional information and details, user access is available to the log files and plate acquisition details.

### Post Compound Scans Not Exported as Expected

Defect ID:9327

Exported trace data for post compound scans are not following the naming convention as others scans. Exported trace data is being written to the same file when more than one compound is used. SW is not picking up the entire name (picks up only the second half of the scan name) of the scan while exporting the data.

### Recovering from Core.Config Corruptions Results in Automatic Restoration of the Proper Mode Selection (Instrument/ Simulator)

Defect ID: 9348

In cases of core.config file corruptions, at the software start up a message appears stating that core.config file has a problem and will be restored using the factory default settings. After closing the message, the software will automatically restore either the Instrument mode or Simulator mode depending on the instrument connection status.



## Known Issues

### Resistance values reported during and after acquisition are not the same

Defect ID:03680

#### Impact

Not all measurements are affected by this problem. Primarily affected measurements include Channel Protocols where the channel pre-signal is a small percentage of the total number of samples collected.

This defect can lead to some confusion and concern about the accuracy of the resistance calculations reported during and after data acquisition. Testing and validation efforts conclude that the resistance measurements reported after data acquisition are correct. Since data analysis is performed after data acquisition, we do not anticipate a significant impact to the users because of this defect.

#### Workaround

None.

### Software fails to load data intermittently with error message: “Some Scans Cannot be Loaded”.

Defect ID:03628

#### Impact

This bug is intermittent. When changing default data folder settings, an error message displays below the **Experiment Name** field that says “Error: Some scans cannot be loaded from file. The data may be corrupt.”

The error message gives the impression that there is something wrong with the data. In reality it means that the software was unable to load the data from the file.

### **Workaround**

Migrate to a different plate. After the display is updated with that newly selected plate, re-select the original plate. If the current data folder contains only one plate, switch to a different data folder. After the display is updated, switch back to the original data folder.

### **Planned Solution**

The issue will be fixed in a future version of the software.

# IonWorks Barracuda Software Version 2.1.1 Software Release Notes

## 9

This chapter provides release notes for version 2.1.1 of the IonWorks Barracuda® Software. Read this entire document before you use the software. The following is a summary of the changes incorporated in this update as compared to version 2.1, the last general release of the IonWorks Barracuda® Software.

- [New in IonWorks Barracuda Software Version 2.1.1](#)
- [Modified](#)

### New in IonWorks Barracuda Software Version 2.1.1

#### Plenum Soak

The **Soak Plenum** functionality has been added to the Plenum Utilities for the purpose of manually starting and ending the plenum soaking with external buffer.

#### Plenum Wash

The **Wash Plenum** functionality has been added to the Plenum Utilities for the purpose of setting automated plenum soaking with external buffer. Settings include **Soak Duration** and **Number of cycles**.

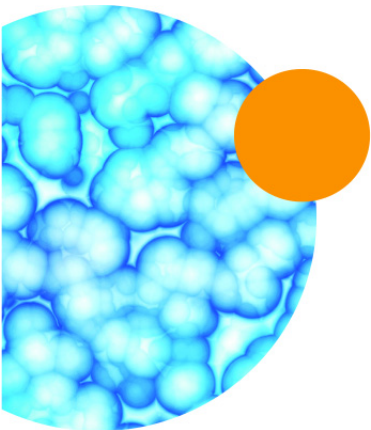
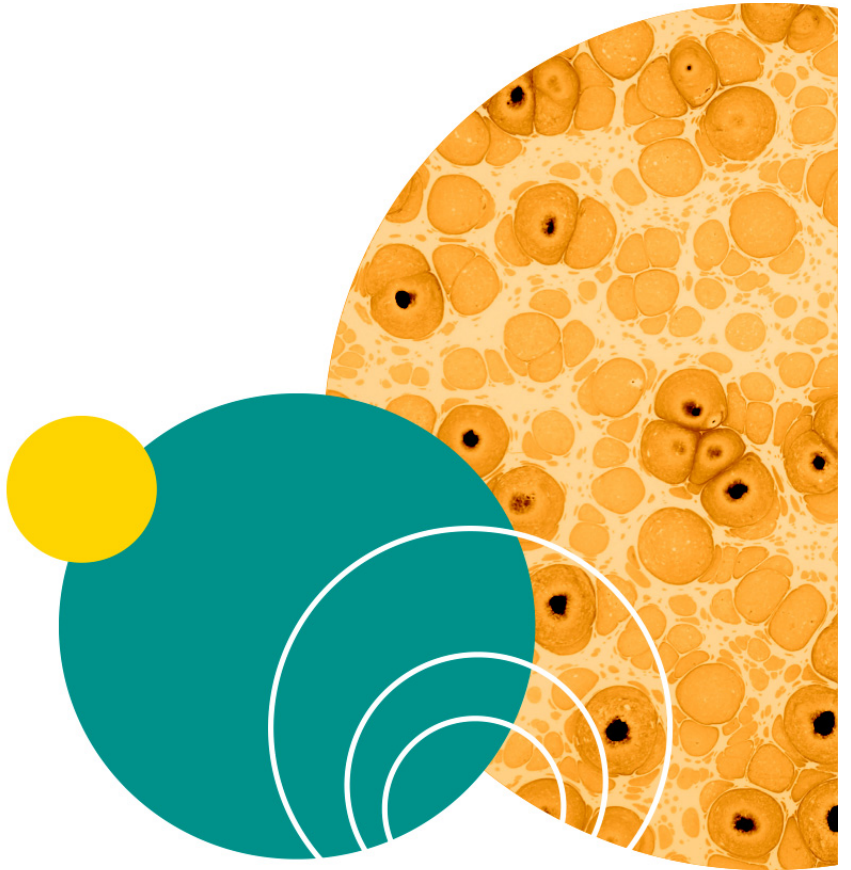
### Modified

#### Utilities Dialog

New Manual Instrument Control buttons added to support the new **Soak Plenum** and **Wash Plenum** functionality.

#### Edit Cleanup Protocol Dialog

Optional **Wash plenum** settings added. Settings include **Soak Duration** and **Number of cycles**.



# IonWorks Barracuda Software Version 2.1 Software Release Notes

10

This chapter provides release notes for Version 2.1 of the IonWorks Barracuda® Software. Read this entire document before you use the software. The following is a summary of the changes incorporated in this update as compared to version 2.0, the last general release of the IonWorks Barracuda® Software.

- [New in IonWorks Barracuda Software Version 2.1](#)
- [Modified on page 29](#)
- [Resolved Issues on page 30](#)
- [Known Issues on page 31](#)

## New in IonWorks Barracuda Software Version 2.1

### iPad and Web Browser Monitor Application Support

View completed experiment data acquired on the IonWorks Barracuda Automated Patch Clamp System remotely through an Intranet connection.



---

**Note:** IonWorks Barracuda Monitor is an optional, and free application for use on a Web browser or an iPad. For the Web application installer, contact Molecular Devices Technical Support. The iPad application can be downloaded from Apple's App Store.

---

## Modified

### Preferences Dialog

New **Preferences > General Options** check box is available for disabling the capturing of plate diagnostic scan screenshots used by the IonWorks Barracuda Monitor applications.

## Resolved Issues

### **When a PatchPlate is not in position, quitting the routine during Start of Day and End of Day Flush and Rinse disables utilities dialog**

Defect ID:03678/5598

The Utilities dialog functions as expected after clicking **Quit**.

### **After changing preferences values in the Instrument Options dialog, the Experiment Mode start button can become disabled**

Defect ID: 03737/5600

The Experiment Mode start button functions as expected after changing preference values in the **Preferences > Instrument Options** dialog.

### **After Renaming A Plate, an Unhandled Exception may Occur in Analysis Summary View**

Defect ID:5601

After renaming a plate, accessing the summary information under Analysis Summary View no longer intermittently crashes the application.

### **Metrics Applied to the Entire Signal are Exported Incompletely**

Defect ID:5604

Exporting a metric to be applied to the entire signal, now contains the entire signal.

### **In a Fresh Installation on the Instrument, the Installer 2.0.0.335 Installs the Simulator Configuration File**

Defect ID:5606

The software now installs the instrument configuration file as expected.

## **Barracuda Trace Export with Leak Correction On, Ends too Soon and Misses Some Trace Exports**

Defect ID:5782

The software now exports all of the traces as expected.

## **Seal Test Quality Control Checks on the First Scan and Aborts the Experiment if it Fails**

Defect ID:5807

The software now performs the Quality Control (QC) check on the last scan.

## **The Default Buffer Volume Setting in the Channel Protocol Editor Displays the Incorrect Volume (5)**

Defect ID:5808

The default buffer volume setting in the Channel Protocol Editor now displays the correct volume (11).

## **The Software Should Show a Warning Message When the Selected Protocol will Exceed the Supported Number of Samples per Scan**

Defect ID:6176

The software now detects the threshold and shows a warning message when the selected protocol will exceed the supported number of samples per scan.

## **Known Issues**

### **Resistance values reported during and after acquisition are not the same**

Defect ID:03680

#### **Impact**

Not all measurements are affected by this problem. Primarily affected measurements include Channel Protocols where the channel pre-signal is a small percentage of the total number of samples collected.

This defect can lead to some confusion and concern about the accuracy of the resistance calculations reported during and after data acquisition. Testing and validation efforts conclude that the resistance measurements reported after data acquisition are correct. Since data analysis is performed after data acquisition, we do not anticipate a significant impact to the users because of this defect.

### **Workaround**

None.

## **Software fails to load data intermittently with error message: “Some Scans Cannot be Loaded”.**

Defect ID:03628

### **Impact**

This bug is intermittent. When changing default data folder settings, an error message displays below the **Experiment Name** field that says “Error: Some scans cannot be loaded from file. The data may be corrupt.”

The error message gives the impression that there is something wrong with the data. In reality it means that the software was unable to load the data from the file.

### **Workaround**

Migrate to a different plate. After the display is updated with that newly selected plate, re-select the original plate. If the current data folder contains only one plate, switch to a different data folder. After the display is updated, switch back to the original data folder.

### **Planned Solution**

The issue will be fixed in a future version of the software.



This chapter provides release notes for version 2.0 of the IonWorks Barracuda® Software. Read this entire document before you use the software. The following is a summary of the changes incorporated in this update as compared to version 1.0.1, the last general release of the IonWorks Barracuda® Software.

- [New in IonWorks Barracuda Software Version 2.0](#)
- [Resolved Issues on page 35](#)
- [Known Issues on page 37](#)

## New in IonWorks Barracuda Software Version 2.0

### Windows 7, 64-bit Operating System Support

IonWorks Barracuda Software v2.0 runs only on a Windows 7, 64-bit operating system.

### Computer Replacement

To accommodate the newly supported Windows 7, 64-bit operating system, the instrument computer gets replaced. A Molecular Devices Field Services Engineer replaces the instrument computer. The computer replacement process includes the v2.0 software update.

### Support for Running Multiple Instances of Analysis Only Mode

New functionality. Multiple instances of the Analysis only software can run simultaneously, because it does not control the instrument, while one instance of the Barracuda data acquisition software runs controlling the instrument.

### Rename an Experiment

New functionality in the Data Acquisition and Review Workspace enables renaming completed experiments. Select a completed experiment from the **Experiment Name** field and then click **Rename**.

## Edit Native Filter Quality Control in Setup Protocol

New functionality enables specifying when a PatchPlate fails by setting the percentage of wells with holes or no seals threshold. Automatically stop the experiment by selecting the **Stop the experiment if the PatchPlate has failed** check box. When the Stop check box is clear, the experiment continues to run and a quality control fail message appears in the status bar.

## Data Review in Assay Development mode

New functionality enables users to pause assay development experiments for reviewing Channel Protocol data in the Data Acquisition and Review workspace. The plate view, Fixed Well view and Floating Well view functions are all active for data review.

## Multiple Protocol Mode

New functionality enables running multiple channel protocols consecutively. This allows users to chain together protocols with different sampling frequencies and combine them in a single experiment.

## Preferences Dialog

New dialog box. From the software toolbar, click **Edit > Preferences** to open the dialog. The Preferences dialog allows you to customize some general settings and some instrument settings. The general options involve the appearance of the software exit message and the plate diagnostic scans. The instrument options involve customizing parameters for the plenum vacuum, compound addition, and mixing.

## Well Statistics in Data Analysis Mode

New interface pane in the Data Analysis Workspace. View metrics for individual wells.

## Experiment Notes

New dialog. You are automatically given the opportunity to add Experiment Notes before running an experiment. You can view the added Experiment Notes after the experiment runs. After the experiment runs, the Experiment Notes can be viewed in the Data Acquisition and Review Workspace in the plate view by clicking the View Experiment Notes icon tab. Experiment Notes in review cannot be edited.

## Set Default Folders

New field. View the last ten saved file paths for the **\Protocols** and **\Data** directories.

## Instrument Utilities Shortcut Toolbar

New icon button to appear on the Shortcut toolbar. Same functionality as selecting the **Instrument > Utility** menu.

## Resolved Issues

### The system ran out of memory when data points exceeded 100K.

Defect ID: 028176

Operating system has changed to a Windows 7, 64-bit. Data points limit has been increased to 600K per scan with or without multiple sweeps.

### Data analysis was disabled for aborted data.

Defect ID: 027604

Software users can analyze all available collected data regardless of an aborted experiments.

### No warning message for Bypass Cellpettor.

Defect IDs: 027469

There is now a warning message at the beginning of the experiment reminding the user to manually place a cellboat with cells on the process deck. Keep in mind that if the cell boat is present and has not been manually replaced, the experiment will still continue irrespective of whether or not there are cells in the boat.

### Pause message appeared even after the bottle was filled.

Defect ID: 027987

The pause message no longer intermittently appears twice.

### **Export family of traces exported only the last trace.**

Defect ID: 028302

The software now exports multiple sweeps of the dataset.

### **Error on closing IonWorks Barracuda Software software application for non-admin users.**

Defect ID: 028346

Administrative files that used to be written in `\ProgramFiles` are now written in `\ProgramData`.

### **Uninstall or reinstall did not start because IonWorks Barracuda Software was running.**

Defect ID: 028716

The message stating that the software is currently running is now enhanced to clearly inform users how to terminate a process before proceeding.

### **If waste bottles were not emptied before an assay, and source bottles were filled, waste bottles might have overflowed undetected during the assay.**

Defect ID: 31777

The software now detects the overflow condition and pauses the assay.

### **Instrument entered fault state when experiment was terminated after a pause.**

Defect ID: 31778

The software no longer goes into a fault state and the system continues to operate properly.

### **CellPettor dimensions were incorrect in the Instrument Overview section of the Help for IonWorks Barracuda Software.**

Defect ID: 31819

The dimensions are now listed correctly in the Help file.

## Known Issues

### Software fails to load data intermittently with error message: “Some Scans Cannot be Loaded”.

Defect ID:03628

#### Impact

This bug is intermittent. When changing default data folder settings, an error message displays below the **Experiment Name** field that says “Error: Some scans could not be loaded from file. The data may be corrupt.”

The error message gives the impression that there is something wrong with the data. In reality it means that the software was unable to load the data from the file.

#### Workaround

Migrate to a different plate. After the display is updated with that newly selected plate, re-select the original plate. If the current data folder contains only one plate, switch to a different data folder. After the display is updated, switch back to the original data folder.

#### Planned Solution

The issue will be fixed in a future version of the software.

### When a PatchPlate is not in position, quitting the routine during Start of Day and End of Day Flush and Rinse disables utilities dialog

Defect ID:03678

#### Impact

At the beginning of the **Start of Day Flush and Rinse** and **End of Day Flush and Rinse** operations, instrument sensors check to make sure that the PatchPlate is in the proper position. When the sensors detect a problem with the PatchPlate placement, the software displays the **User Intervention Required** dialog and unlocks the process deck door to enable users to properly position the PatchPlate.

If the user fixes the problem and clicks **Continue**, the operation functions as expected. If the user clicks **Quit**, it has been noticed that all the

options on the **Utilities** dialog are disabled and the software does not allow users to perform any other operations.

### **Workaround**

Exit and restart the application to continue.

### **Planned Solution**

This issue will be fixed in the next version of the software.

## **Resistance values reported during and after acquisition are not the same**

Defect ID:03680

### **Impact**

Not all measurements are affected by this problem. Primarily affected measurements include Channel Protocols where the channel pre-signal is a small percentage of the total number of samples collected.

This defect may lead to some confusion and concern about the accuracy of the resistance calculations reported during and after data acquisition.

Testing and validation efforts conclude that the resistance measurements reported after data acquisition are correct. Since data analysis is performed after data acquisition, we do not anticipate a significant impact to the users because of this defect.

### **Workaround**

None.

### **Planned Solution**

The issue will be fixed in the next version of the software.

## **After changing preferences values in the Instrument Options dialog, the Experiment Mode start button may become disabled**

Defect ID: 03737

### **Impact**

This is an intermittent problem. When in the **Edit > Preferences >Instrument Options** dialog, changing and saving values may result in

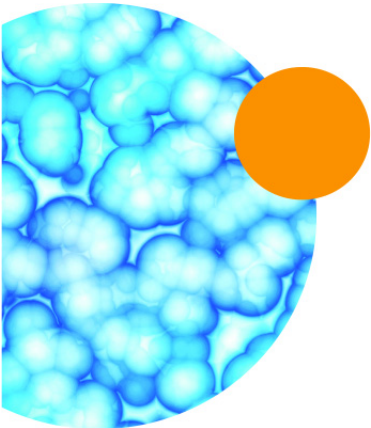
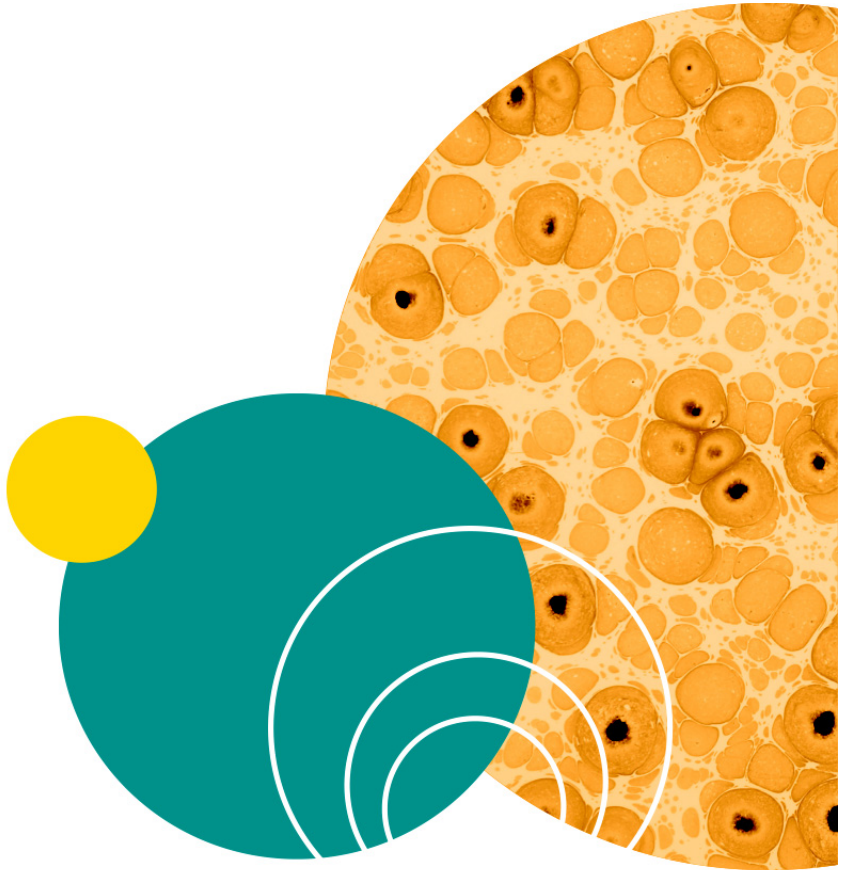
the **Experiment Mode** start button being disabled and the **Running** message appearing in the **Instrument Status** bar.

### **Workaround**

Exit and restart the application to continue.

### **Planned Solution**

The issue will be fixed in the next version of the software.





## Contact Us

Phone: [800.635.5577](tel:800.635.5577)  
Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
Email: [info@moldev.com](mailto:info@moldev.com)

Check our website for a current listing  
of worldwide distributors

## Regional Offices

USA and Canada <a href="tel:+1.800.635.5577">+1.800.635.5577</a>	China (Beijing) <a href="tel:+86.10.6410.8669">+86.10.6410.8669</a>	Japan (Osaka) <a href="tel:+81.6.7174.8831">+81.6.7174.8831</a>	Brazil <a href="tel:+55.11.3616.6607">+55.11.3616.6607</a>
United Kingdom <a href="tel:+44.118.944.8000">+44.118.944.8000</a>	China (Shanghai) <a href="tel:+86.21.3372.1088">+86.21.3372.1088</a>	Japan (Tokyo) <a href="tel:+81.3.6362.5260">+81.3.6362.5260</a>	
Europe* <a href="tel:00800.665.32860">00800.665.32860</a>	Hong Kong <a href="tel:+852.2248.6000">+852.2248.6000</a>	South Korea <a href="tel:+82.2.3471.9534">+82.2.3471.9534</a>	

*\*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden, and Switzerland*