



IonWorks® product update

IonWorks Software Version 2.0.5 Release Notes

5 November, 2013

IonWorks Software Version 2.0.5 has been developed to address known issues in IonWorks Version 2.0.3. Included in this software upgrade is a firmware upgrade from Version 2.0.3 to 2.0.5. This upgrade is for customers currently using IonWorks Software Version 2.0 and higher only. If you are running a version earlier than IonWorks Software Version 2.0, you should contact your local Molecular Device Representative to get upgraded to IonWorks Software Version 2.0. A summary of the changes incorporated into IonWorks Software Version 2.0.5 are:

- **Fix – When interleaving was used, incorrect compound mapping occurred.**
When interleaving is used, compound mapping is now correct.
- **Fix – Conditioning Train used in combination with Read Bank Interleaving resulted in incorrect command-voltage response measurements.**
When Conditioning Train is used with Read Bank Interleaving, the command-voltage responses measure correctly.
- **Fix – Two Help menu hyperlinks opened incorrect Support site webpages.**
The Help menu hyperlinks now open the correct Support site webpages.

Upgrading the Instrument Firmware

The Firmware Upgrade utility allows you to install new versions of the instrument firmware without requiring a field service call. In most cases firmware updates will accompany updates of the Host application as well. This procedure covers updating both software components.

Note: This utility can only be used to update an instrument that is currently running version 2.0 Firmware. Upgrading to v2.0 from an earlier version requires a call to Molecular Devices Field Service.

How to install software updates:

| Step | Action |
|------|---|
| 1 | Uninstall the previous version of IonWorks V2 application. Make sure you are logged in to the instrument using an administrator's account. If you go to the Control Panel and select Add/Remove Programs and then click on IonWorks V2, you should see buttons labeled Change and Remove. If not, you don't have the proper rights. If so, select Remove. The uninstaller will appear. Click the Next button if necessary to start the uninstaller. Click Close if necessary to close the uninstaller when it is finished. |
| 2 | Install the new version. Put in the CD and wait for the drive folder to open. Once you see the Setup2.msi file, double-click on it to run the installer. Select Yes to agree to the license, then select the option to install for Everyone. Keep clicking Next until the installer starts copying files. Click the Close button when the installer is finished. |
| 3 | If the unit is not on, power it up. |
| 4 | Once the unit has fully booted up and the motion stages have homed (the D-axis of the pipettor will go through the homing sequence), start the Application. The splash screen should show the new version. |
| 5 | Wait for the green connect light to come on. |
| 6 | When the Host has connected, select Utilities / Upgrade Firmware . Wait for the dialog to appear, it may take a few moments. |
| 7 | When the Firmware Upgrade Utility dialog is visible, it should show the currently installed instrument Firmware version. The Firmware version we are installing is shown in the dialog also, labeled as 'Source File Version'. The network connection section should show that the connection to the instrument has been established. |
| 8 | Click the Terminate Firmware button to shut down the Firmware. A timer will count the number of seconds. It typically takes about 10-12 seconds for the process to complete. |
| 9 | When the dialog indicates the Firmware has terminated, click the Update Firmware button. A message will quickly appear indicating the Firmware has been updated. Power down the unit as instructed, wait about a minute and then power the unit back up. |
| 10 | Close the Firmware Upgrade Utility dialog. The green connect light will be dark. |
| 11 | When the unit has finished rebooting, the green connect light will come on again. The unit is ready for use. |