



ClonePix™ 2 Clone Picking Software

Version 1.4.14

Software Release Notes

August 2013



www.moleculardevices.com

ClonePix 2 Software Release Notes

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.

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

Product manufactured by Molecular Devices, LLC.
1311 Orleans Drive, Sunnyvale, California, United States of America 94089.
Molecular Devices, LLC is ISO 9001 registered.
© 2013 Molecular Devices, LLC.
All rights reserved.

Contents

Chapter 1: ClonePix 2 Software General Information	5
Starting the System and the Software	5
Obtaining Support	6
Chapter 2: ClonePix 2 Software Version 1.4.14 Software Release Notes	7
Issue Addressed in ClonePix 2 Software v1.4.14	7
Chapter 3: ClonePix 2 Software Version 1.4 Software Release Notes	9
Modifications Made to ClonePix 2 Software v1.4	9
Issues Addressed in ClonePix 2 Software v1.4	10

Chapter 1: ClonePix 2 Software General Information

The ClonePix™ 2 Clone Picking Software controls the ClonePix™ 2 Clone Picking System.

The following topics are included in this chapter:

- [Starting the System and the Software on page 5](#)
- [Obtaining Support on page 6](#)

Starting the System and the Software

Before starting the system and the software confirm the following conditions:

- The **Emergency Stop** button on the front panel of the instrument is pulled out. The instrument will not start if this button is pushed in.
- The instrument bed is clear of obstructions and loose items.
- All motor tracks are free of obstruction.
- There are no obstructions to movement of the head.
- The main robot axes, XY, are roughly in the center of their respective travel positions. The axes need to be away from their respective end-stops on start up so that they can reference themselves correctly to their home positions.
- The instrument door is fully closed.
- No plates or cassettes are in the stackers.

Power-Up Procedure

1. Turn on the power supply to the compressor.
 2. Push the **Start** button on the front panel of the instrument.
The **Power On** light illuminates on the front indicator panel. If the power to the system does not turn on, it is possible that the door is open or the Emergency Stop button is pushed in.
The instrument cycles through various start-up processes indicated on the front indicator panel.
 3. Check that the **Air Pressure OK** icon illuminates on the front indicator panel.
 4. Switch on the computer and wait for it to finish initializing.
 5. From the computer desktop, double-click the ClonePix 2 Clone Picking Software icon.
- Every time the instrument is used, the three axes sequentially run through their “Initialize drives” routine. This enables the drives to find their respective home positions. The system must be allowed to complete this routine without interference to ensure that there is no damage to the instrument or its ancillary equipment.

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, 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.



WARNING! BIOHAZARD. It is your responsibility to decontaminate components of the instrument before requesting service by a service engineer or returning parts to Molecular Devices for repair. Molecular Devices will not accept any items which have not been decontaminated where it is appropriate to do so. If any parts are returned, they must be enclosed in a sealed plastic bag stating that the contents are safe to handle and are not contaminated.

Chapter 2: ClonePix 2 Software Version 1.4.14 Software Release Notes

The ClonePix™ 2 Clone Picking Software version 1.4.14 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 1.4, the last general release of the software.

Issue Addressed in ClonePix 2 Software v1.4.14

The following issue was addressed in ClonePix 2 Clone Picking Software version 1.4.14.

Camera Substage Misalignment After Restart

Tracking ID: 4398

After restarting the instrument and the software, the camera substage sometimes does not align properly during initialization. This intermittent issue can effect the accuracy of picking.

Resolution:

The initialization routine properly aligns the camera substage for consistent picking accuracy.

Impact of fix:

This fix has no impact on current workflow or data.

Chapter 3: ClonePix 2 Software Version 1.4 Software Release Notes

The ClonePix™ 2 Clone Picking Software version 1.4 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 1.2.94, the last general release of the software.

- [Modifications Made to ClonePix 2 Software v1.4 on page 9](#)
- [Issues Addressed in ClonePix 2 Software v1.4 on page 10](#)

Modifications Made to ClonePix 2 Software v1.4

The following modifications were made to ClonePix 2 Clone Picking Software version 1.4.

Subsystem Firmware Update

A critical subsystem had its a firmware updated, and the software was modified to work with the new firmware. Before running this version of the ClonePix 2 Clone Picking Software, you need to run a firmware update for this subsystem.

The firmware update and installer are contained in a file named **Molecular Devices Firmware Loader.zip**. If you did not receive this file along with the software upgrade, contact Molecular Devices Technical Support. See [Obtaining Support on page 6](#).

To update the subsystem firmware:

1. Download the **Molecular Devices Firmware Loader.zip** file to the system computer.
2. Power on the system.
3. Extract the contents of the **Molecular Devices Firmware Loader.zip** file.
4. Open the folder where you extracted the contents of the **Molecular Devices Firmware Loader.zip** file and then double-click the **MolecularDevicesFirmwareLoader.exe** file.
5. In the **Molecular Devices Firmware Loader** utility, click **Update Firmware**.

If the firmware is up-to-date, the utility will inform you of this and discontinue the firmware update process.



CAUTION! After starting the firmware update, do not close the utility until after the Status message indicates that the firmware update is complete.

6. Click **Close**.



Note: A system firmware update is also required for this release. See [The Firmware Times Out Giving a Transport Layer Error on page 10](#).

Issues Addressed in ClonePix 2 Software v1.4

The following issues were addressed in ClonePix 2 Clone Picking Software version 1.4.

The Firmware Times Out Giving a Transport Layer Error

Tracking ID: 4046

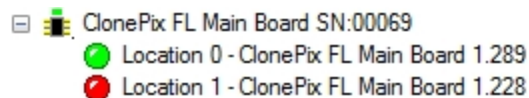
The instrument firmware times out causing the instrument to periodically freeze at various times during use.

Resolution:

A thread management issue was resolved in firmware version 1.289 requiring a firmware update for existing systems.

To view the current firmware version:

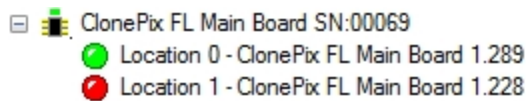
1. In the ClonePix 2 Software window, click **Tools > Configuration**.
2. In the **Edit Configuration** window, double-click **Cold Fire Devices**.
3. Under **ClonePix Main Board**, view the firmware version for the listing next to the green dot.



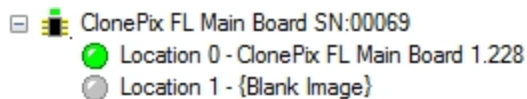
The firmware should be version 1.289 or later. If you have an older version of the firmware, contact Molecular Devices technical support to obtain a file containing the latest version of the firmware. See [Obtaining Support on page 6](#).

To update the firmware:

1. Save the firmware-update file on the system computer.
2. In the ClonePix 2 Software window, click **Tools > Configuration**.
3. In the **Edit Configuration** window, double-click **Cold Fire Devices**.
4. Under **ClonePix Main Board**, right-click the listing next to the red dot and then click **Update Firmware**.



If there is no red dot, then right-click the listing next to the gray dot.

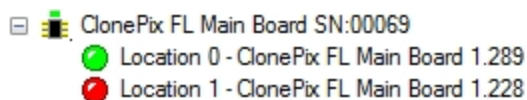


CAUTION! Do not update the firmware for the listing next to the green dot, since this is the version that is currently running on the system.

5. In the **Open** dialog, navigate to and select the firmware-update file you saved in step 1.
6. Click **Open** to start the firmware update.
An hourglass replaces the icon next to **ClonePix Main Board**.
7. After the icon reappears next to **ClonePix Main Board**, click **Close**.
8. Exit the software.
9. Turn off and restart the instrument.
10. Restart the software.

To confirm the updated firmware version:

1. In the ClonePix 2 Software window, click **Tools > Configuration**.
2. In the **Edit Configuration** window, double-click **Cold Fire Devices**.
3. Under **ClonePix Main Board**, view the firmware version for the listing next to the green dot.



The firmware should be version 1.289 or later.

Impact of fix:

This fix has no impact on current workflow or data.

Picking Pins are Misaligned

Tracking ID: 4108

The picking pins alignment is incorrect reducing the picking accuracy.

Resolution:

The picking pins now align properly.



Note: This issue has also been resolved for side-by-side installations of the ClonePix 2 Software and the ClonePix FL Software.

Impact of fix:

This fix has no impact on current workflow or data.

Picking Pins Sometimes Miss the Target Area if Fluorescence is Used as the Prime Configuration

Tracking ID: 4166

If fluorescence is used as the Prime Configuration, the picking pins sometimes miss the target area and can strike the edges of the source tray.

Resolution:

The picking pins now align properly when fluorescence is used as the Prime Configuration.

Impact of fix:

This fix has no impact on current workflow or data.

Contact Us

Phone: +1-800-635-5577
Web: www.moleculardevices.com
Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada +1-800-635-5577
Latin America +55 11 3616 6607
China (Beijing) +86 10 6410 8669
China (Shanghai) +86 21 3372 1088
China (Hong Kong) +852 3971 3520

Japan (Tokyo) +81 3 6362 5260
Japan (Osaka) +81 6 7174 8831
South Korea +82 2 3471 9531
United Kingdom +44 (0) 118 944 8000
Germany/Austria +49 (0) 89 9605 880

