



# ClonePix™ 2 Clone Picking Software

Version 1.4.14

## Software Upgrade Guide

5029604 A  
September 2013



[www.moleculardevices.com](http://www.moleculardevices.com)

## ClonePix 2 Software Upgrade 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 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.

## Chapter 1: Upgrading ClonePix 2 Software

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

This document describes how to upgrade the ClonePix 2 Clone Picking Software from a previous version of the software. If you need to install the software on a computer that does not already have a version of the software installed, see the *ClonePix 2 Software Installation Guide*.

The following topics are included in this document:

- [Computer System Requirements on page 3](#)
- [Backing Up the Previous Version Configuration on page 4](#)
- [Uninstalling the Previous Version of the Software on page 4](#)
- [Installing the ClonePix 2 Software Upgrade on page 4](#)
- [Subsystem Firmware Update on page 5](#)
- [Obtaining a License for the Software on page 5](#)
- [Starting the System and the Software on page 6](#)
- [Re-Installing the Previous Version of the Software on page 7](#)
- [Obtaining Support on page 7](#)

### Computer System Requirements

The ClonePix 2 software upgrade can be installed on a computer with the a previous version of the software already installed and the following system specifications:

**Table 1-1: Minimum Computer System Requirements**

Item	Description
Operating system	Windows 7, 32-bit (x86) The software has not been validated on the Windows 7, 64-bit (x-64) operating system.
Memory	2 GB RAM
Data Connection	10/100 Ethernet port
Camera Connection	USB 2.0 port

## Backing Up the Previous Version Configuration

Before removing the previous installation of the software, you must back up the configuration in case you need to reinstall it later.

1. Click **Start > All Programs > Molecular Devices > ClonePix 2 > Configuration Manager**.
2. In the **Fusion Configuration Manager**, click **Backup To File**.
3. In the **Backup Fusion Configuration** dialog, type 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, type a **File name**, and then click **Save**.

## Uninstalling the Previous Version of the Software

1. Click **Start > Control Panel**.
2. Under **Programs**, click **Uninstall a program**.
3. From the list of programs, click **ClonePix 2**.
4. Click **Uninstall**.
5. Follow the on-screen instructions to finish uninstalling the software.

## Installing the ClonePix 2 Software Upgrade

1. Download the upgrade installation file to the system computer.
2. Double click the **ClonePix 2 v.1.4.14.msi** file to start the installation program.
3. Follow the on-screen instructions to finish installing the software upgrade.



**Note:** Before starting the software, update the subsystem firmware. See [Subsystem Firmware Update on page 5](#).

---

## Subsystem Firmware Update

A critical subsystem had its 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 7](#).

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**.

## Obtaining a License for the Software

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

To license the software:

1. Contact Molecular Devices customer support to acquire a valid license file. See [Obtaining Support on page 7](#).
2. From the computer desktop, double-click the ClonePix 2 Clone Picking Software icon.
3. Follow the on-screen instructions to license the software.

## 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.

## Re-Installing the Previous Version of the Software

If for any reason the newly installed version of software is not functioning correctly you can re-install the previous version.

1. Uninstall the newly installed software as described in [Uninstalling the Previous Version of the Software on page 4](#).  
You do not need to back up the configuration for this version.
2. Insert the Recovery DVD or locate the downloaded installation file for the previous version of software.
3. Start the installation program and follow the on-screen instructions to re-install the software.
4. Click **Start > All Programs > Molecular Devices > ClonePix 2 > Configuration Manager**.
5. In the **Fusion Configuration Manager**, click **Restore From File**.
6. Locate and select the backup file that you saved in [Backing Up the Previous Version Configuration on page 4](#).
7. Click **Open**.
8. After the configuration is restored, close the **Fusion Configuration Manager**.
9. Start and license the software. See [Obtaining a License for the Software on page 5](#).

## 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.



**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.

---

#### Contact Us

Phone: +1-800-635-5577  
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 +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

