



# ClonePix 2

Mammalian Colony Picker

Software Release Notes

## 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 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.

©2023 Molecular Devices, LLC.

All rights reserved.



# Contents

<b>Chapter 1: ClonePix 2 Software General Information</b> .....	<b>5</b>
Computer System Requirements .....	5
Backing Up the Previous Version Configuration .....	6
Uninstalling the Previous Version of the Software .....	6
Subsystem Firmware Update .....	7
Installing the ClonePix 2 Software Upgrade .....	7
Updating the System Firmware .....	8
Obtaining a License for the Software .....	9
Re-Installing the Previous Version of the Software .....	12
Obtaining Support .....	13
<b>Chapter 2: ClonePix 2 Software Version 1.6.16 Software Release Notes</b> .....	<b>15</b>
New in ClonePix 2 Software v1.6.16 .....	15
Issue Addressed in ClonePix 2 Software v1.6.16 .....	15
<b>Chapter 3: ClonePix 2 Software Version 1.6 Software Release Notes</b> .....	<b>17</b>
Modifications Made to ClonePix 2 Software v1.6 .....	17
<b>Chapter 4: ClonePix 2 Software Version 1.5.52 Software Release Notes</b> .....	<b>19</b>
Issue Addressed in ClonePix 2 Software v1.5.52 .....	19
<b>Chapter 5: ClonePix 2 Software Version 1.5.49 Software Release Notes</b> .....	<b>21</b>
Issue Addressed in ClonePix 2 Software v1.5.49 .....	21
<b>Chapter 6: ClonePix 2 Software Version 1.5.46 Software Release Notes</b> .....	<b>23</b>
Issue Addressed in ClonePix 2 Software v1.5.46 .....	23
<b>Chapter 7: ClonePix 2 Software Version 1.5.45 Software Release Notes</b> .....	<b>25</b>
New in ClonePix 2 Software v1.5.45 .....	25
Modifications Made to ClonePix 2 Software v1.5.45 .....	25
<b>Chapter 8: ClonePix 2 Software Version 1.5 Software Release Notes</b> .....	<b>27</b>
Issue Addressed in ClonePix 2 Software v1.5 .....	27



# Chapter 1: ClonePix 2 Software General Information



The ClonePix<sup>®</sup> 2 Clone Picking Software controls the ClonePix<sup>®</sup> 2 ClonePix 2 Mammalian Colony Picker.

The following topics are included in this chapter:

- [Computer System Requirements on page 5](#)
- [Backing Up the Previous Version Configuration on page 6](#)
- [Uninstalling the Previous Version of the Software on page 6](#)
- [Subsystem Firmware Update on page 7](#)
- [Installing the ClonePix 2 Software Upgrade on page 7](#)
- [Installing the ClonePix 2 Software Upgrade on page 7](#)
- [Obtaining a License for the Software on page 9](#)
- [Re-Installing the Previous Version of the Software on page 12](#)
- [Obtaining Support on page 13](#)

## Computer System Requirements

The ClonePix 2 software requires the following computer specifications:

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

Item	Description
Operating system	Windows 11 Windows 10, 64-bit Windows 7, 32-bit
Memory	8 GB RAM or more
Data Connection	10/100 Ethernet port
Camera Connection	SBIG: USB 2.0 port 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.

The recommended process is two steps:

1. [Add a Windows System Restore Point on page 6](#)
2. [Back up ClonePix 2 Software Configuration File on page 6](#)

### 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 message **The restore point was created successfully** appears, click **Close**.
6. In any other related open dialogs, click **Close** to finish.

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

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

## Subsystem Firmware Update

The most recent version of the software is modified to work with new firmware for a critical subsystem. If you are updating from a Clone Picking Software version older than version 1.5.45, you must run this firmware update before updating the ClonePix 2 Clone Picking Software. If the firmware is up-to-date, the utility informs you of this and ends the firmware update process. After this firmware version verification, you can continue installing the ClonePix 2 Clone Picking Software.

The firmware update and installer are contained in a file named **FirmwareLoader-5.5.0.zip**. If you did not receive this file along with the software upgrade, contact Technical Support. See [Obtaining Support on page 13](#).

To run the subsystem firmware update utility:

1. Download the **FirmwareLoader-5.5.0.zip** file to the system computer.
2. Power on the ClonePix 2 system computer and instrument.
3. Extract the contents of the **FirmwareLoader-5.5.0.zip** file.
4. Open the folder where you extracted the contents of the **FirmwareLoader-5.5.0.zip** file and then double-click the **MolecularDevicesFirmwareLoader.exe** file.
5. In the **FirmwareLoader-5.5.0** utility, click **Update Firmware**.



**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**.

## Installing the ClonePix 2 Software Upgrade

1. Download the upgrade installation file to the system computer.
2. Double click the **ClonePix 2 v.1.6.16.msi** file to start the installation program.
3. Follow the on-screen instructions to finish installing the software upgrade.



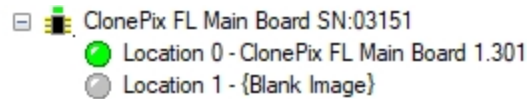
**Note:** Before starting the software, update the subsystem firmware. See [Subsystem Firmware Update on page 7](#).

---

## Updating the System Firmware

To view the current firmware version:

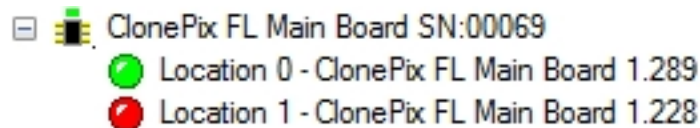
1. In the ClonePix 2 Software window, click **Tools > Configuration**.
2. In the **Edit Configuration** window, double-click **Cold Fire Devices**.
3. Under **ClonePix FL Main Board**, view the firmware version for the listing next to the green dot.



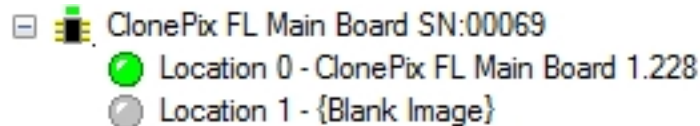
The firmware should be version 1.301 or higher. If you have an older version of the firmware, contact technical support to obtain a file containing the latest version of the firmware. See [Obtaining Support on page 13](#).

To update the firmware:

1. Power on the ClonePix 2 system computer and instrument.
2. Save the firmware-update file on the system computer.
3. In the ClonePix 2 Software window, click **Tools > Configuration**.
4. In the **Edit Configuration** window, double-click **Cold Fire Devices**.
5. Under **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.

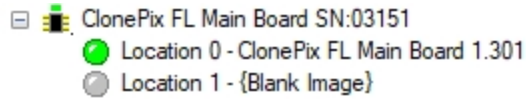
---

6. In the **Open** dialog, navigate to and select the firmware-update file you saved in step 1.
7. Click **Open** to start the firmware update.  
An hourglass replaces the icon next to **Main Board**.
8. After the icon is restored next to **Main Board**, click **Close**.
9. Exit the software.
10. Turn off and restart the instrument.
11. 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 **Main Board**, view the firmware version for the listing next to the green dot.



The firmware should be version 1.301 or later.

## 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.

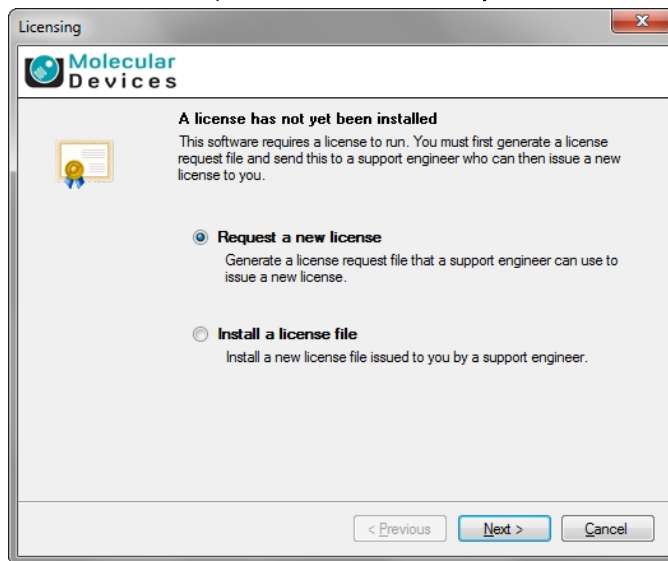
Licensing is a two step process:

1. [Requesting a Software License on page 9](#)
2. [Installing a Software License on page 11](#)

### 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 text "User", "Company/Institute Name\*" with the text "Company", and "Instrument Serial Number (optional):". There is also a large text area for "Notes/Comments:". At the bottom, it says "\*These fields are required" and has three buttons: "< 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 bottom buttons remain: "< Previous", "Next >", and "Cancel".

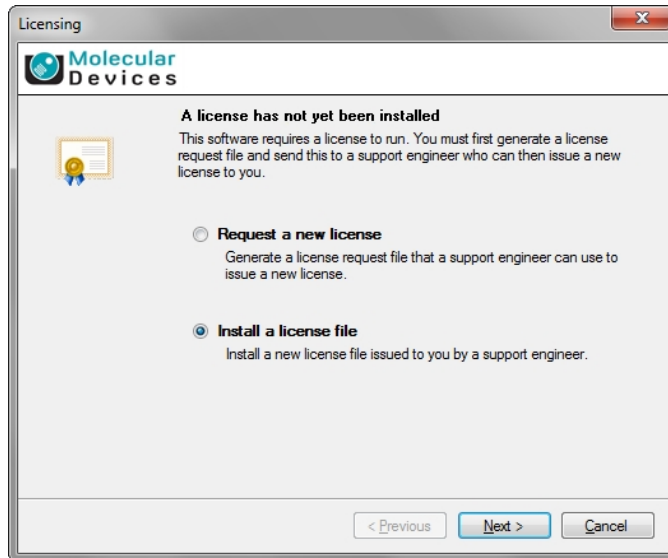
5. In the Save the Request to a File dialog, save the file in a location where you can find it.
6. To close the wizard, click **Finish**.
7. Change the **LicenseRequest.req** file name to a **LicenseRequest.txt** file, and then email it to Molecular Devices Technical Support. See [support.moleculardevices.com/](http://support.moleculardevices.com/).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

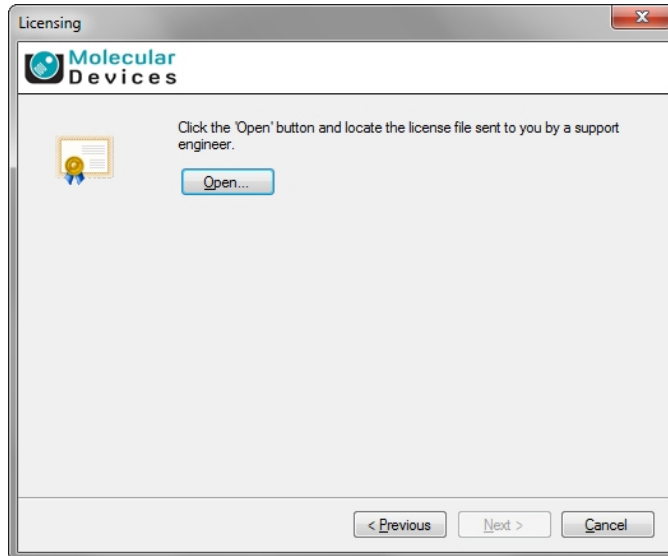
After you receive the license from Molecular Devices Technical Support, do the following:

To install a software license:

1. From the computer desktop, double-click the CloneSelect ImagerClonePix 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**.

## Re-Installing the Previous Version of the Software

If for some 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 6](#).  
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 >> 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 6](#).
  7. Click **Open**.
  8. After the configuration is restored, close the **Fusion Configuration Manager**.
  9. If you are downgrading to Clone Picking Software version 1.5.44 or older, downgrade the firmware:
    - a. Download the **FirmwareLoader-4.1.2.zip** file to the system computer.
    - b. Power on the ClonePix 2 system computer and instrument.
    - c. Extract the contents of the **FirmwareLoader-4.1.2.zip** file.
    - d. Open the folder where you extracted the contents of the **FirmwareLoader-4.1.2.zip** file and then double-click the **MolecularDevicesFirmwareDowngrade.exe** file.
    - e. In the **Firmware Loader** utility, click **Update Firmware**.  
If the firmware is up-to-date, the utility informs you of this and discontinues the firmware update process.
- 
-  **CAUTION!** After starting the firmware installation, do not close the utility until after the Status message indicates that the firmware installation is complete.
- 
- f. Click **Close**.
10. Start and license the software. See [Obtaining a License for the Software on page 9](#).

## 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/](https://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.

---



## Chapter 2: ClonePix 2 Software Version 1.6.16 Software Release Notes



The version 1.6.16 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 1.6, the last general release of the software.

### New in ClonePix 2 Software v1.6.16

The following new features are included in version 1.6.16.

#### Windows 11

Added support for the Windows 11 operating system.

### Issue Addressed in ClonePix 2 Software v1.6.16

The following issue was addressed in version 1.6.16.

#### Pin Datuming Out of Focus

Tracking ID: 5363

When running the source plate datuming and pin offset datuming, the pin tip image view is sometimes blurry, and is worse using the Basler camera.

#### Resolution:

The source plate datuming and pin offset datuming pin tip image view is now clearly in focus.

#### Impact of fix:

This fix has no impact on current workflow or data.

#### Navitar Board Ports Not Closing Consistently

Tracking ID: 5341

The Navitar hardware board ports are not closing properly intermittently, which causes an error when trying to launch the software the next time.

#### Resolution:

The Navitar hardware board ports now close consistently as expected.

#### Impact of fix:

This fix has no impact on current workflow or data.





## Chapter 3: ClonePix 2 Software Version 1.6 Software Release Notes



# 3

The version 1.6 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 1.5.52, the last general release of the software.

### Modifications Made to ClonePix 2 Software v1.6

The following modifications were made to ClonePix 2 Clone Picking Software version 1.6.

#### **New Navitar PCBA Support**

The current PCBA board is obsoleted by the vendor and replaced by a new Navitar PCBA board. This ClonePix Software update supports both versions of the board.

#### **New Basler Camera Support**

A Basler camera replaces the SBIG camera in new instruments. This ClonePix Software update supports both new and existing cameras.



## Chapter 4: ClonePix 2 Software Version 1.5.52 Software Release Notes



The version 1.5.52 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 1.5.49, the last general release of the software.

### Issue Addressed in ClonePix 2 Software v1.5.52

The following issue was addressed in version 1.5.52.

#### Intermittent Software Crash During FITC Imaging

Tracking ID: 5308

When the FITC acquisition option is selected, the software intermittently crashes during a picking run.

#### Resolution:

The software now retries capturing correct data from the camera.

#### Impact of fix:

This fix has no impact on current workflow or data.



## Chapter 5: ClonePix 2 Software Version 1.5.49 Software Release Notes



The version 1.5.49 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 1.5.46, the last general release of the software.

### Issue Addressed in ClonePix 2 Software v1.5.49

The following issue was addressed in version 1.5.49.

#### Intermittent Instrument Connection Error Using the Newest Digi RealPort Driver

Tracking ID: 5282

Using the newest Digi RealPort driver (version 4.7.410), there is an intermittent error when the software tries to connect to the ThorFilterWheel of the ClonePix 2 ClonePix 2 instrument.

#### Resolution:

The software now tries to connect a few times.

#### Impact of fix:

This fix has no impact on current workflow or data.



## Chapter 6: ClonePix 2 Software Version 1.5.46 Software Release Notes



The version 1.5.46 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 1.5.45, the last general release of the software.

### Issue Addressed in ClonePix 2 Software v1.5.46

The following issue was addressed in version 1.5.46.

#### Incorrect Order of Source and Destination Plates in Results Overview and Exported Reports

Tracking ID: 5260

After a batch run where several plates are used as source and destination, the results overview and exported reports list the plates in the wrong order in which they were picked and deposited.

#### Resolution:

The results overview, the exported statistics, and the HTML report now all list the plates in the correct order.

#### Impact of fix:

This fix has no impact on current workflow or data.





## Chapter 7: ClonePix 2 Software Version 1.5.45 Software Release Notes



The version 1.5.45 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 1.5, the last general release of the software.

### **New in ClonePix 2 Software v1.5.45**

The following new features are included in version 1.5.45.

#### **Windows 10 Operating System Support**

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

### **Modifications Made to ClonePix 2 Software v1.5.45**

The following modifications were made to ClonePix 2 Clone Picking Software version 1.5.45.

#### **ClonePix FL Software End of Life**

ClonePix FL Software is no longer supported.



## Chapter 8: ClonePix 2 Software Version 1.5 Software Release Notes



The version 1.5 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 1.4.14, the last general release of the software.

### Issue Addressed in ClonePix 2 Software v1.5

The following issue was addressed in version 1.5.

#### Error in Barcode Reader Causes Plate Skipping

Tracking ID: 5131

During a batch picking, if a source plate barcode cannot be read, the plate gets skipped, which affects the order of plate deposit.

#### Resolution:

Now the source plate barcode reads only once, resulting in a complete and accurate picking and deposit.

#### Impact of fix:

No data changes, however the workflow is changed. The instrument reads the barcodes only during the read, and only once.

#### Picking Order is Opposite of Plate Read Preview Order

Tracking ID: 5195

The picking read preview shows the destination plate template wells in the order of the source plates, however, the picking actually picks in the reverse source plate order.

#### Resolution:

The picking read preview and the actual picking order are now the same.

#### Impact of fix:

This fix has no impact on current workflow or data.



## 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.