

# Quick Start Guide

Thank you for purchasing the ImageXpress™ Pico Automated Cell Imaging System and the CellReporterXpress™ Image Acquisition and Analysis Software from Molecular Devices®. Review this guide to get started.

## ImageXpress Pico Automated Cell Imaging System

The ImageXpress Pico Automated Cell Imaging System is an affordable all-in-one platform for automatically acquiring and analyzing images from fluorescently labeled biological samples in plates and slides. It enables you to increase the throughput of your image acquisition and analysis, allowing you to gain insights in minutes. With the addition of a modular option, the system provides environmental control for live cell imaging.

The core hardware component of the imaging system is a custom-designed, fully automated, epi-illumination fluorescence microscope. The rapid autofocus and precision sample movement features of the microscope allow large numbers of high-resolution images to be acquired in the shortest possible time. All key optical and mechanical elements are motorized, which allows complete control of the instrument configuration.

When used in combination with the CellReporterXpress Image Acquisition and Analysis Software, the ImageXpress Pico System becomes an extremely flexible device, ideally suited for user-defined, automated assays.

## CellReporterXpress Image Acquisition and Analysis Software

The CellReporterXpress Image Acquisition and Analysis Software is the user interface for the ImageXpress Pico System.

The CellReporterXpress Software integrates image acquisition and analysis into a unified workflow. Along with the ImageXpress Pico System, the CellReporterXpress Software streamlines automated imaging to offer a simplified solution for scaling up microscopy. Its features include:

- A web-based interface that runs on many browsers, including those found on iPads and Android tablets.
- Over 25 available predefined experimental protocols.
- High-powered analysis tools equivalent to those found in desktop applications.
- Easy-to-manage data with no requirement to configure a database.
- A simplified user interface that is easy to learn and easy to use.

## 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, [www.moleculardevices.com/service-support](http://www.moleculardevices.com/service-support), has a link to the Knowledge Base, which contains technical notes, software upgrades, safety data sheets, and other resources. If you still need assistance after consulting the Knowledge Base, you can submit a request to Molecular Devices Technical Support.

## Technical Support

You can contact Molecular Devices Technical Support by submitting a support request through the Knowledge Base or by phone. To find regional support contact information, visit [www.moleculardevices.com/contact](http://www.moleculardevices.com/contact).

You will need the instrument serial number and the software system ID.

## Documentation

Review the product documentation on the Knowledge Base, including installation guides and user guides. In addition, online Help is available within the CellReporterXpress Software. Press **F1** to access Help for the active page.

## Additional Resources

Web-based microscopy courses:

- [www.leica-microsystems.com/science-lab](http://www.leica-microsystems.com/science-lab)
- [www.ibiology.org/ibioeducation/taking-courses/ibiology-microscopy-short-course.html](http://www.ibiology.org/ibioeducation/taking-courses/ibiology-microscopy-short-course.html)

The Molecular Probes Handbook offers advice on fluorescent probes and can help you determine if there are better stains available for your analysis:

- [www.lifetechnologies.com/us/en/home/references/molecular-probes-the-handbook.html](http://www.lifetechnologies.com/us/en/home/references/molecular-probes-the-handbook.html)

## Product Documentation

The following guides are available on the Molecular Devices Knowledge Base at [mdc.custhelp.com](http://mdc.custhelp.com):

- *CellReporterXpress Installation Guide*
- *CellReporterXpress Release Notes*
- *CellReporterXpress User Guide*
- *ImageXpress Pico Pre-Installation Guide*
- *ImageXpress Pico Installation Guide*
- *ImageXpress Pico User Guide*
- *ImageXpress Pico Calibration Kit Guide*

In addition, the CellReporterXpress Software includes context-sensitive Help that you can access from within the software. Just press the **F1** key from within the software to view Help for the current page.



**Tip:** Molecular Devices recommends that you review the documentation before installing or using the ImageXpress Pico System or the CellReporterXpress Software.

## Items Required to Install the Software

You will need the following items to install the CellReporterXpress Software:

- **Host Computer:** Unless you purchased the ImageXpress Pico System with the optional computer, you must provide the computer that will be the host computer.
- **Dongle:** Molecular Devices provides the SafeNet Sentinel USB hardware key (called a "dongle"), which contains the license required to run the software.



- **Installer:** Download the CellReporterXpress Installation and Configuration Utility from the Molecular Devices Knowledge Base. Go to [www.meta.moleculardevices.com/pico](http://www.meta.moleculardevices.com/pico) and click the **CellReporterXpress Software (CRX)** link under **Downloads**.

CellReporterXpress Imaging Acquisition and Analysis Software

## Next Steps

Review the *ImageXpress Pico Installation Guide* for details on unpacking and setting up your ImageXpress Pico System.

Review the *CellReporterXpress Installation Guide* for details on installing and configuring your CellReporterXpress Software.

All documentation is available in the Knowledge Base on the Molecular Devices Support website at [www.moleculardevices.com/service-support](http://www.moleculardevices.com/service-support).