

CellReporterXpress Imaging Software Quick Start

Items Required to Complete the Software Installation

- The **CellReporterXpress Image Acquisition and Analysis Software Installer** flash drive contains the software components and product documentation.
- The **SafeNet Sentinel** CellReporterXpress Software USB hardware key (called a dongle) contains licensing information required to run the CellReporterXpress Software.



Software Components on the Flash Drive

- **CellReporterXpress Image Acquisition and Analysis Software** is the primary tool for acquiring and analyzing images.
- **CellReporterXpress Software Help** file is available on the flash drive and also within the software.

Before You Install the Software Components

Do the following tasks before installing any of the components that are on the CellReporterXpress Software flash drive:

- Make sure your computer has at minimum a 2.4 GHz CPU, 10 logical processors, and at least 12 GB of RAM to support the CellReporterXpress Software and 4 concurrent analysis processes.
- Make sure the computer is running the 64-bit Windows 10 Professional or Enterprise operating system.
- Make sure that you are logged in as a user with administrative privileges on the computer where you are installing the software.
- Make sure the computer can connect to the imaging device either through the local network or through a direct Ethernet connection.

For help, visit www.moleculardevices.com/support.

CellReporterXpress Imaging Software Quick Start

Software Installation and Start

1. Plug in the CellReporterXpress Software Installation flash drive.
If the CellReporterXpress Software Installation window does not appear after you plug in the flash drive, then browse to the flash drive folder on your computer, right-click the **setup.exe** file, and click **Run as administrator**.
2. Plug in the **SafeNet Sentinel** USB hardware key.
3. Click **Display End-User License Agreement**. Review the license agreement, select **I accept the End-User License Agreement**, and then click **Continue**.
4. Confirm or install the required components.
5. Confirm or install the drivers for the **SafeNet Sentinel** USB hardware key.
6. Confirm or install the license file that matches the **System ID** of the **SafeNet Sentinel** USB hardware key.
7. On the **Install the CellReporterXpress Software** page, next to **CellReporterXpress : Complete**, click **Install**.
8. To start the software, double-click on the **MD.CellReporterXpress** icon on your desktop.
9. In the log in screen, type your Windows credentials, and then click **LOG IN**.

For information about connecting to the instrument and more advanced software configurations, including firewall and anti-virus setting, see the *CellReporterXpress Software Installation and Configuration User Guide*.

Documentation Available on the Flash Drive

The CellReporterXpress Software flash drive contains installation and user guides for the supported software and hardware.

The following documents are available in the **Manuals** folder on the flash drive:

- *CellReporterXpress Software Installation and Configuration User Guide*
- *ImageXpress® Nano Automated Imaging System With CellReporterXpress Software User Guide*

Contact Us

Phone: [+1-800-635-5577](tel:+1-800-635-5577)
Web: moleculardevices.com
Email: info@moldev.com

Visit our website for a current listing of worldwide distributors.

The trademarks used herein are the property of Molecular Devices, LLC or their respective owners.
Specifications subject to change without notice. Patents: www.moleculardevices.com/productpatents
FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

© 2017 Molecular Devices, LLC.
All rights reserved.
5058431 B