

MetaXpress Imaging Software Suite

Items Required to Complete the Software Installation

- The **MetaXpress® Imaging Software Suite Installer** flash drive contains the software components and product documentation.



- The **SafeNet Sentinel** MetaXpress Software USB hardware key (sometimes called a dongle) contains licensing information required to run the MetaXpress Software. This is either a single-license hardware key or a multi-license network key. You can install a multi-license network key on a server, so that multiple client computers can run the analysis software from a single hardware key. Other hardware keys might be required depending on your installation configuration. All required hardware keys are provided in the package, or can be obtained from your Molecular Devices representative.



Note: The MetaXpress Software and related components must be installed on a 64-bit Windows operating system. For more information about supported operating systems, see the *Molecular Devices High-Content Computer and Server Specifications for Offline Analysis* guide located on the Knowledge Base.

Read this entire document before installing the software components.

- [Software Components on the Flash Drive, see page 2](#)
- [Required Reading, see page 2](#)
- [Before You Install the Software Components, see page 3](#)
- [Documentation Available on the Flash Drive, see page 4](#)

Software Components on the Flash Drive

- **MetaXpress® High-Content Image Acquisition and Analysis Software** is the primary tool for acquiring and analyzing images.
- **MetaXpress® PowerCore™ High-Content Distributed Image Analysis Software** is an optional component that accelerates image analysis by distributing image-processing tasks across multiple computer processes.
- **MDCStore™ High Content Data Management Solution** is a database schema that can be used with the Microsoft SQL Server or Oracle platforms. It supports multi-server and multi-user configurations.
- **MDCStore Tools™ Data Management Utility** is a stand-alone tool for setting up and maintaining an MDCStore database schema, managing user accounts and data, and optimizing the supporting SQL and Oracle servers.
- **MDCStore™ Xchange Data Conversion Service** is an optional service that imports and exports images to and from the MDCStore database.
- **MDC File Server** is an optional network service used to manage file servers to store MetaXpress Software images outside the MDCStore database.
- **ImageXpress® System USB Driver** is the driver installation program required for the ImageXpress® Micro System.
- **Andor SDK 3** is the driver installation program required for an ImageXpress® Micro System that uses an Andor camera.
- **Microsoft SQL Server 2014 Express** is a small database server that is useful for demonstration or evaluation purposes. Its size is limited to 10 GB, making it impractical for production environments.

Required Reading

Before you install the MetaXpress Software and its related components, read the information in the following documents:

- *Molecular Devices High Content Screening Computer Specifications*
- *Molecular Devices Complete Solution HCS IT Requirements*

These documents are available on the flash drive or in the knowledge base on the Molecular Devices support site: www.moleculardevices.com/support. Click **Visit the Knowledge Base**.

Before You Install the Software Components



CAUTION! The regional settings on your computer must be set to **English (United States)** even if you are located outside of the United States. If this setting is changed to anything other than **English (United States)**, the MetaXpress Software will not work correctly. This setting is found in the **Control Panel** in **Region and Language**.

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

- 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 C: drive has at least 1 GB of available space after all other requirements have been met to install the software.

To install all the required prerequisites and all the software on the flash drive, you need at least 4 GB of available space.

- Make sure that the appropriate database server software is installed:
 - For databases larger than 10 GB (most databases), Microsoft SQL Server 2012 or 2014, or Oracle 11g or 12c, must be installed and accessible. Microsoft SQL Server can be pre-installed on instrument workstations.



Note: Molecular Devices recommends installing the database on a separate server. See [Required Reading, see page 2](#).

These products are not available on the MetaXpress Software Suite flash drive. If you need assistance determining which product is most appropriate, contact your System Administrator or a Molecular Devices representative.

- For databases smaller than 10 GB (in general, databases that contain data from fewer than 10 plates), you can use Microsoft SQL Server 2014 Express, which is suitable for desktop operation. The 64-bit version of SQL Server 2014 Express is provided on the MetaXpress Software Suite flash drive.
If SQL Server is already installed, then SQL Server 2014 Express cannot be installed.
- If the MetaXpress Software Suite installation window does not appear after you insert the flash drive, then browse to the flash drive folder on your computer, right-click the **autorun.exe** file, and click **Run as administrator**.

MetaXpress Imaging Software Suite

Documentation Available on the Flash Drive

The MetaXpress Software Suite flash drive contains user guides and installation guides for the software and hardware supported by this release of the MetaXpress Software.

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

- *Molecular Devices High Content Screening Computer and Server Specifications for Offline Analysis*
- *Molecular Devices Complete Imaging Solution Network Requirements*
- *MDCStore™ High Content Data Management Solution Database Schema Installation and Upgrade Guide*
- *MDCStore Tools™ Data Management Utility User Guide*
- *MDCStore™ Xchange Data Conversion Service Quick Start Guide*
- *MetaXpress® PowerCore™ High-Content Distributed Image Analysis Software User Guide*
- *MetaXpress® PowerCore™ High-Content Distributed Image Analysis Software Release Notes*
- *MetaXpress® High-Content Image Acquisition and Analysis Software Installation and Update Guide*
- *MetaXpress® High-Content Image Acquisition and Analysis Software Analysis Guide*
- *ImageXpress® Micro Widefield High Content Imaging System User Guide*
- *ImageXpress® Micro Confocal High Content Imaging System User Guide*

In addition, the MetaXpress Software application help is available when running the software.

Contact Us

Phone: +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

© 2016 Molecular Devices, LLC.
All rights reserved.
5039824 C