

MetaXpress Offline Software with IN Carta Software

Read Me First

Do the following to get started with the MetaXpress® High-Content Image Acquisition and Analysis Software (offline) and the IN Carta® Image Analysis Software:

- Connect the MetaXpress software USB hardware key to the analysis computer.
- Install the IN Carta software.
- License the IN Carta software.
- Connect the preinstalled MetaXpress software to a MDCStore data source.

Connecting the USB Hardware Key

The **SafeNet Sentinel** MetaXpress Software USB hardware key (sometimes called a dongle) contains licensing information required to run the MetaXpress offline software on the analysis computer. Insert the dongle in a USB port on the analysis computer to get started.



Installing the IN Carta Software

The analysis computer that ships from Molecular Devices is set up to run both the MetaXpress software and the IN Carta software. Your IT Department may have changed some settings to comply with your organization's policy. In that case, before you begin, confirm that the following is true:

- The Regional Format setting in Windows is set to **English (United States)**.
- The user name for the Windows account on the analysis computer contains only alphanumeric characters. No special characters (including the period or full stop character) are allowed. You may need to work with your IT Department to get an account that meets this requirement.

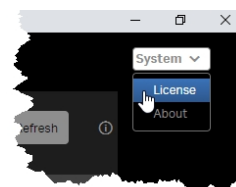
To install the IN Carta software on the analysis computer:

1. In Windows, open File Explorer and browse to the C:\INCarta_Installers folder.
2. Right-click **INCarta_1.x.x.exe** and select **Run as administrator**.
3. Review the license agreement.
4. If you accept the license agreement, select the **I agree to the license terms and conditions** check box.
5. Click **Install**.
6. If a message displays about monitor resolution, click **Yes** to continue installing.
7. When the installation completes, click **Close**.

Licensing the IN Carta Software

To license the IN Carta software on the analysis computer:

1. On the Windows Desktop, double-click the **INCARTA** icon.
2. If a Windows Security Alert displays, click **Allow Access**.
3. In the upper right corner of the IN Carta window, click **System > License**.
4. In the License Information window, confirm the licenses you purchased are selected.
5. Click **Close**.



Connecting to the MDCStore Data Source

After installing and licensing the IN Carta software, connect the MetaXpress offline software on the analysis computer to the MDCStore data source.

To connect to the MDCStore data source:

1. On the Windows Desktop, right-click the **MetaXpress** icon, and select **Run as administrator**.
2. In the Welcome to MetaXpress dialog, click **New Data Source**.
3. Select the **Microsoft SQL Server or MSDE** option and click **OK**.
If you are using Oracle, see the *MetaXpress Installation and Update Guide*.
4. In the **Computer\Instance** field, enter the database server name and the instance of the MDCStore data source.



Note: If you do not know the database server name and the database name, log in to another computer that connects to the MDCStore data source (for example, the acquisition computer) and select **Help > About the Database**.

5. In the **Login Name** and **Password** fields, enter the system administrator credentials for the MDCStore data source.
6. Click **Connect**.
7. In the **Database Name** field, confirm the database.
8. In the **New Data Source** field, use the suggested name.
9. Click the **ODBC Driver** drop-down, and select the latest driver.
10. Click **Create**.

Obtaining Support

To contact Molecular Devices Technical Support, submit a support request through the Molecular Devices Knowledge Base at support.moleculardevices.com.

You can also submit a support request by phone. For regional support contact information, go to www.moleculardevices.com/contact.

Contact Us

Phone: +1-800-635-5577
Web: moleculardevices.com
Email: info@moldev.com

Visit our website for a current listing of worldwide distributors.