



PRE-INSTALLATION GUIDE

IN Carta

Image Analysis Software

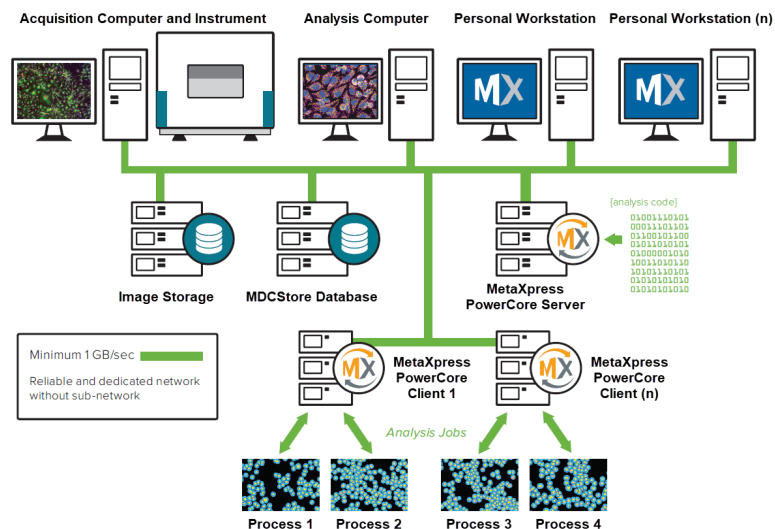
Welcome to the family of Molecular Devices software users. This guide provides you with the information required to prepare your site for your new software.

The IN Carta™ Image Analysis Software combines powerful analytics with a modern intuitive interface to simplify workflows for advanced phenotypic classification and 3D image analysis.

When used in combination with the MetaXpress® High-Content Image Acquisition and Analysis Software, the IN Carta software makes it easy to embrace the complexity in an image. You can get meaningful data quickly and reliably.

To help you take advantage of all that the software has to offer, the IN Carta software package also includes:

- A GPU-enabled, artificial intelligence workstation—the Analysis Computer—that becomes an integral part of your ImageXpress® High-Content Imaging System.
- An offline version of the MetaXpress software.



Package Contents

The IN Carta software package includes the following:

- Analysis Computer: Dell Precision 5280 Tower
- 22" or 27" Monitor (optional)
- Keyboard and Mouse
- Power Supply
- Monitor Cables
- MetaXpress Dongle Case with USB Hardware Key

The Analysis Computer that ships from Molecular Devices is configured to run both the IN Carta software and the MetaXpress offline software. After you set up the computer, you are just a few quick steps away from getting started.



Optional 27" Monitor



Analysis Computer: Dell Precision Tower 5280

Space and Table Requirements

The Analysis Computer requires a table suitable for the size and weight of the components, which are as follows:

Component	Height	Width	Depth	Weight
Computer	41.8 cm (16.5 in.)	17.7 cm (7.0 in.)	51.8 cm (20.4 in.)	15.4 kg (34 lb)
Monitor 27" (with stand)	39.5 cm (15.6 in.)	61.2 cm (24.1 in.)	18.5 cm (7.3 in.)	6.9 kg (15.2 lb)
Monitor 22" (with stand)	35.3 cm (13.9 in.)	48.7 cm (19.2 in.)	16.6 cm (6.5 in.)	4.7 kg (10.4 lb)

To minimize vibration, the Analysis Computer should not be placed on the same table as the ImageXpress instrument. The Analysis Computer can be used with multiple monitors, so ample desk space is needed for monitor placement.



IN Carta Image Analysis Software

Power Requirements

Consider the following power requirements for the Analysis Computer:

- Direct connections to all international supply voltages are available.
- Use the included power supply to connect the Analysis Computer to a GROUNDED power receptacle.
- Fluctuations must be within 10% of the nominal voltage.
- When using a power strip, connect the Analysis Computer to a different power strip than the ImageXpress Micro Systems Power and Options Controller (which connects to the instrument and light source).

To limit the risk of interruption during power loss, use an uninterruptible power supply (UPS) to provide backup power and power line conditioning for the Analysis Computer. We recommend that the UPS capacity is 20% higher than the power requirement for your system.

Component	Watts	VA	Power Cables
Analysis Computer	120	133	1
27" Monitor (each)	40	44	1
22" Monitor (each)	20	22	1

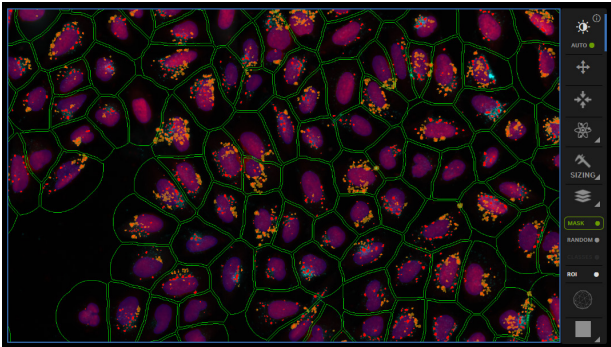
To determine the power consumption (watts) or apparent power (volt-amperes or VA) for the Analysis Computer, add the power requirements of all applicable components. For example, the power consumption of a computer with two large monitors is 200 watts (120 + 40 + 40).

Facility Receiving Requirements

The IN Carta software package arrives in multiple boxes. The size and weight of the boxes are as follows:

Box	Height	Width	Depth	Weight
Computer and Accessories	59.7 cm (23.5 in.)	67.9 cm (26.8 in.)	30.5 cm (12 in.)	20.4 kg (45 lb)
Monitor 27"	31.2 cm (12.3 in.)	61.3 cm (24.1 in.)	21.5 cm (8.5 in.)	9.6 kg (21.16 lb)
Monitor 22"	37.9 cm (14.9 in.)	49.9 cm (19.7 in.)	23.3 cm (9.18 in.)	7.5 kg (16.62 lb)

In addition, the USB Hardware Key arrives in its own box along with a Read Me First sheet that will help you get started with the Analysis Computer and the software.



IN Carta Image Viewer: Cell and Organelle Segmentation Masks

Installation Requirements

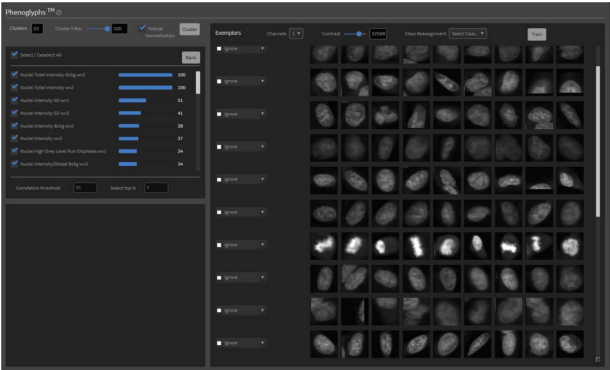
You will need to provide the following to set up the Analysis Computer:

- Sturdy table appropriate for the size and weight of the Analysis Computer with the keyboard, mouse, and at least one monitor
- Network cable and 1 gigabit network connection
- Online power conditioning UPS (optional)
- Power strip (optional)

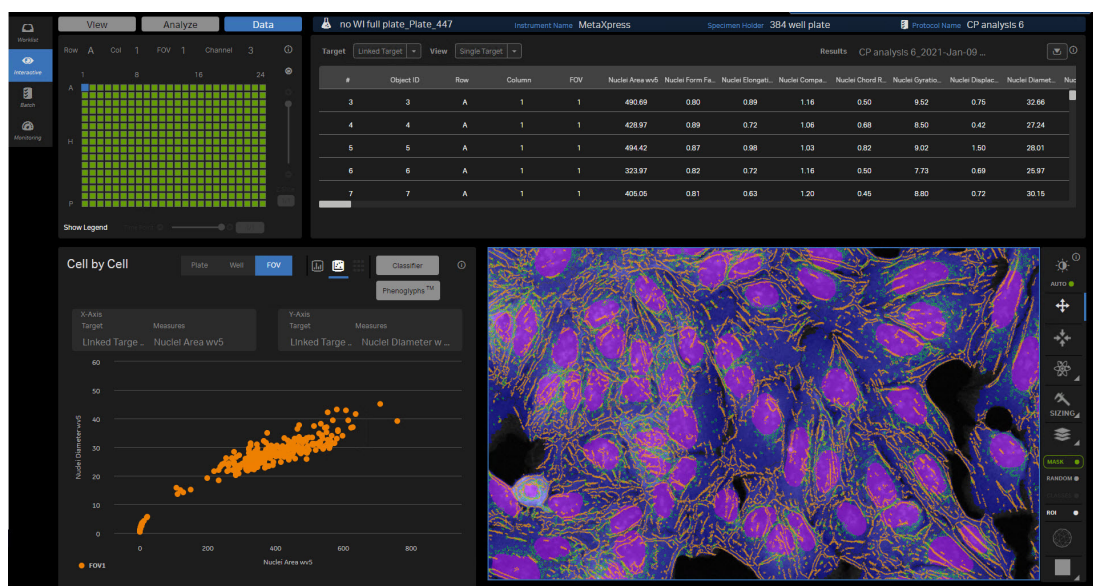
After setting up the Analysis Computer, you will need to do the following:

- Connect the USB Hardware Key to the Analysis Computer.
- Install and license the IN Carta software.

See the Read Me First sheet in the package with the USB Hardware Key for details.



IN Carta Software: Phenoglyphs™ Machine Learning Classification Module



IN Carta Software: Intuitive User Interface

IT Requirements

The Analysis Computer is installed on your network, so you may need to involve your IT professional during installation. In addition, consider the following IT requirements:

- The MetaXpress offline software must be connected to an MDCStore data source, which may reside on the Acquisition Computer or (more likely) as a Microsoft SQL Server database on your network. See the Read Me First sheet in the package with USB Hardware Key for details.
- Windows Defender is pre-installed on the Analysis Computer with all the required settings and exclusions to run the IN Carta software and the MetaXpress offline software. If you are not using Window Defender, you must create these exclusions on your own. See the Molecular Devices High-Content Screening IT Requirements for details.
- The user name for the Windows account on the Analysis Computer cannot contain any special characters (including the period or full stop character). Only alphanumeric characters are allowed.

Documentation

The Molecular Devices Knowledge Base includes complete documentation for the IN Carta software, the MetaXpress software, and the ImageXpress system. Along with user guides, you will find technical notes, software upgrades, safety data sheets, and other resources. To access the Knowledge Base, go to support.moleculardevices.com.

Support and Service

Molecular Devices provides comprehensive support and service solutions for the ImageXpress Confocal HT.ai system supported by a global network of factory trained engineers. To contact us, go to:

Web: www.moleculardevices.com/service-support
www.moldev.com/support
 Email: support@moldev.com

Contact Us

Web: www.moleculardevices.com
 Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada
 UK and Europe*

+1.800.635.5577
 00800.665.32860

China +86.400.821.3787
 Taiwan +886.2.2656.7585

Japan +81.3.6362.9109
 Korea +82.2.3471.9531

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Luxembourg, Netherlands, Sweden, and Switzerland