



IN Carta

Image Analysis Software

Version 2.1

Installation Guide



IN Carta Image Analysis Software Installation Guide

This document is provided to customers who have purchased Molecular Devices equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a non-transferrable license. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Each product is shipped with documentation stating specifications and other technical information. Molecular Devices products are warranted to meet the stated specifications. Molecular Devices makes no other warranties or representations express or implied, including but not limited to, the fitness of this product for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom. The sole obligation of Molecular Devices and the customer's sole remedy are limited to repair or replacement of the product in the event that the product fails to do as warranted.

For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

The IN Carta software is sold under use license from Cellomics Inc.

Patents: www.moleculardevices.com/patents

Product manufactured by Molecular Devices, LLC.
3860 N. First Street, San Jose, California, 95134, United States of America.
Molecular Devices, LLC is ISO 9001 registered.
©2023 Molecular Devices, LLC.
All rights reserved.



Contents

Chapter 1: IN Carta Image Analysis Software	5
Obtaining Support	6
About This Guide	6
Chapter 2: IN Carta Software Requirements	7
Chapter 3: Before Installing the IN Carta Software	9
Setting the Windows Region Formats	9
Removing an Earlier Version of the Software	13
Checking the Computer	13
Chapter 4: Installing the IN Carta Software	15
Downloading the Software Installers	16
Installing the Software	17
Verifying the GPU Driver for the SINAP Module and IN Carta 3D Viewer	18
Installing the SINAP Deep-Learning Models	19
Activating the Software License	20
Updating Image Stacks, Results Data, and Protocols	22
Chapter 5: After Installing the IN Carta Software	27
Granting Folder Permissions	27
Granting Software Permissions	29
Chapter 6: Setting Up Network Licensing	31
Setting Up a Network Licensing Server	31
Setting Up a Network Licensing Client	32



Chapter 1: IN Carta Image Analysis Software



1

The IN Carta® Image Analysis Software provides powerful analytics for advanced phenotypic classification and 3D image analysis. It delivers robust, quantitative results from complex biological images and datasets using advanced AI technology.

Faster Data

- Intuitive design makes complex analysis accessible with minimal training.
- Shorten analysis time with true parallel processing.

Reliable Data

- Sophisticated algorithms generate reliable data with minimal user input.
- Improved segmentation algorithms represent cellular structures more accurately.

Results That Matter

- See real results quickly—from populations to single cells—using integrated data visualization tools.
- User-friendly interface guides you through your discoveries with continual updates that grow with your needs.

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—describes the support options offered by Molecular Devices, including service plans and professional services. It also has a link to the Molecular Devices Knowledge Base, which contains documentation, technical notes, software upgrades, safety data sheets, and other resources. If you still need assistance, you can submit a request to Molecular Devices Technical Support.

Technical 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.

To expedite support, be prepared to provide the software version and your activation ID. To

display this information, at the bottom left of the IN Carta window, click  **Settings**.

Documentation

Review the product documentation on the Molecular Devices Knowledge Base at support.moleculardevices.com. In addition, online Help is available within the IN Carta software.

Additional Resources

Web-based microscopy courses:

- www.microscopyu.com
- 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.thermofisher.com/us/en/home/references/molecular-probes-the-handbook.html

The *Assay Guidance Manual* details state-of-the-art approaches to high-content screening (HCS) and discusses challenges specific to HCS. It serves as a good introduction for new HCS practitioners.

- www.ncbi.nlm.nih.gov/books/NBK100913

About This Guide

This guide is intended for the scientist or IT professional who will be installing and configuring the IN Carta software. It contains critical information that helps you successfully install the software.

The information in this guide is valid for IN Carta software version 2.1 and is subject to change without notice. We recommend that you review the guide on the Molecular Devices Knowledge Base at support.moleculardevices.com for the most up-to-date information.



CAUTION! We recommend installing the IN Carta software on a computer that meets the minimum requirements. If you install on a computer that does not meet the minimum requirements, Molecular Devices cannot guarantee system functionality and performance.

The following table describes the minimum and recommended requirements for the IN Carta computer:

Item	Requirement
Operating System	One of the following: <ul style="list-style-type: none"> • Windows 10 Pro, 64-bit • Windows 10 Enterprise 64-bit • Windows 11 Pro, 64-bit • Windows 11 Enterprise 64-bit
Windows Region	English (US)
Processor Frequency	Minimum: 3.2 GHz Recommended: 4.6 GHz
Processor Cores	Minimum: 4 Recommended: 18
RAM	Minimum: 16 GB Recommended: 128 GB
Storage	Minimum: 1 TB (50 GB free) Recommended: Both a 2 TB SSD (50 GB free) and a 4 TB HDD
Nvidia GPU Memory (optional, see below)	Minimum: 4 GB Recommended: 24 GB
CUDA Compute Capability of Nvidia GPU (optional, see below)	3.5
Monitor Size	Optimized for a 24-inch screen
Monitor Resolution	Optimized for 1920 × 1080



Note: An Nvidia graphics processing unit (GPU) is required only for the optional SINAP module or the optional IN Carta 3D Viewer.




This section describes the procedures required before running the IN Carta software installer:

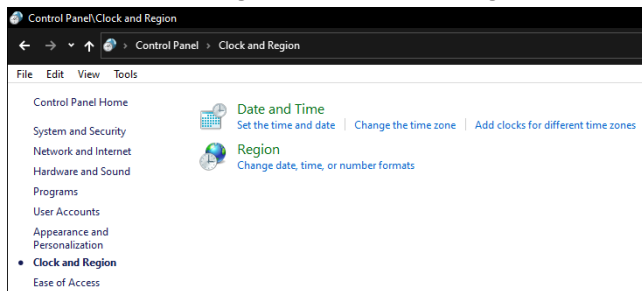
- [Setting the Windows Region Formats, see below](#)
- [Removing an Earlier Version of the Software, see page 13](#)
- [Checking the Computer, see page 13](#)

Setting the Windows Region Formats

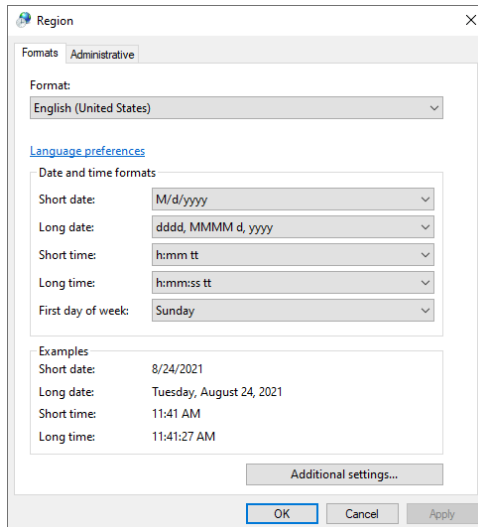
Installing the IN Carta software requires that the Windows Region Formats be set to English (United States).

To set Windows Region Formats for the software:

1. In Windows, click  **Start**, enter `control`, and select **Control Panel**.
2. Click the **View by** drop-down, and select **Category**.
3. Click **Clock and Region**.
4. In the **Clock and Region** window, click **Region**.

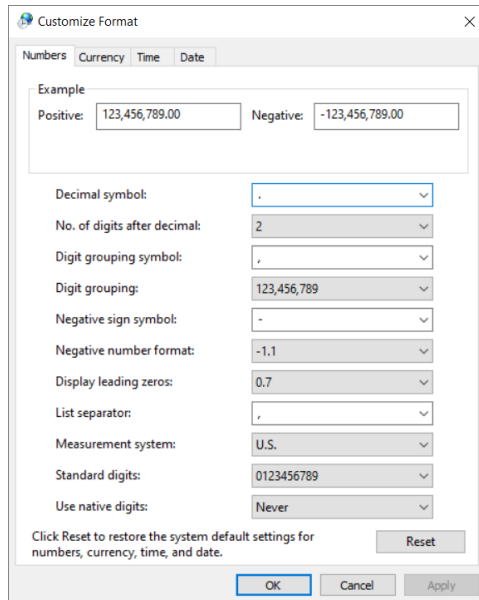


5. In the **Region** window, on the **Formats** tab, set the **Format** field to **English (United States)**.

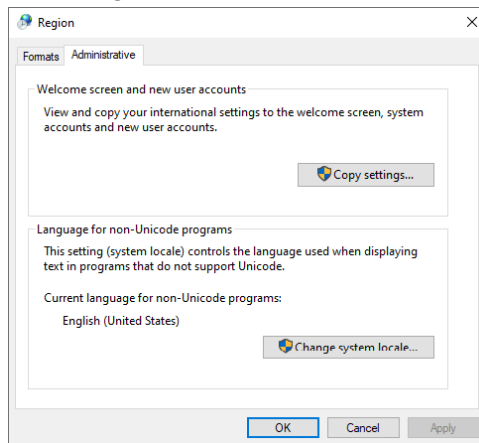


6. Click **Additional settings**.

- In the **Customize Format** window, on the **Numbers** tab, click **Reset**. This ensures that the default English language number settings are used.



- Click **Yes**.
- Click **OK**.
- In the **Region** window, click the **Administrative** tab.



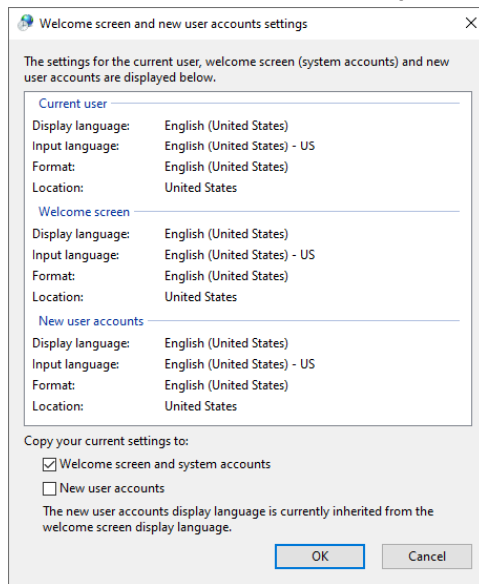
- On the **Administrative** tab, set the **Language for non-Unicode programs** to **English (United States)**.



Note: If **English (United States)** is not already set, click **Change system locale** to set it.

- In the **Welcome screen and new user accounts** section, click **Copy settings**.
- If prompted, click **Apply**.

14. Select the **Welcome screen and system accounts** check box.



Note: This check box may not be available in all Windows versions.

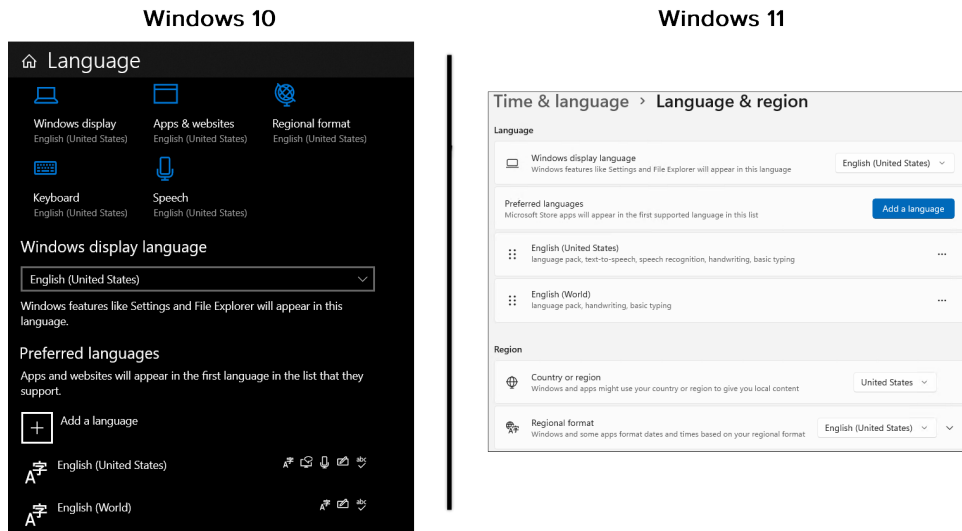
15. Click **OK**.
16. Click the **Formats** tab.
17. Click **Language preferences**.

18. On the **Language** window (with Windows 10) or the **Language & Region** window (with Windows 11), set the **Windows display language** to **English (United States)**.



Note: If **English (United States)** is not available in the drop-down, click **Add a language** to install it.

19. Under **Preferred languages**, set **English (United States)** as the default language. The default language appears at the top of the list of languages.



Note: If **English (United States)** is not already set, click and drag it to the top of the list.

20. If you changed any settings, restart the computer.

Removing an Earlier Version of the Software

If you are upgrading to version 2.1 of the IN Carta software from version 1.17 or version 1.16, you must uninstall the earlier version of the software before running the installer. This includes the following:

- INCarta
- INCarta_Base
- INCarta_Models



Tip: We recommend restarting the computer BOTH before and after uninstalling an earlier version of the software.



Note: If you are upgrading the IN Carta software from a version earlier than version 1.16, contact Molecular Devices Technical Support. See [Obtaining Support on page 6](#) for details.

Checking the Computer

To run the IN Carta software, confirm that the following is true:

- The computer meets the minimum requirements described in [IN Carta Software Requirements on page 7](#).
- The **User Account Control Setting** in Windows settings (**Control Panel > System and Security > Security and Maintenance > Change User Account Control**) is set to **Never Notify** during installation.



Note: This setting is only required during installation. You can revert the setting after the software is installed.

- You are logged in to the computer with a Windows administrator account.
- The user name for the Windows administrator account on the IN Carta computer contains only alphanumeric characters. No special characters (including spaces, periods, or underscores) are allowed. You may need to work with your IT professional to get an account that meets this requirement.
- No other installation of the IN Carta software exists on the computer. See [Removing an Earlier Version of the Software on this page](#) for details.
- The IN Carta folder (C:\Program Files\INCarta) does not exist on the computer. If it does exist, delete it.
- All files, folders, and programs are closed.
- The Windows Region Formats are set to **English (United States)** as described in [Setting the Windows Region Formats on page 9](#).



This section describes the procedures required to install version 2.1 of the IN Carta software:

- [Downloading the Software Installers](#), see page 16
- [Installing the Software](#), see page 17
- [Verifying the GPU Driver for the SINAP Module and IN Carta 3D Viewer](#), see page 18 (optional)
- [Installing the SINAP Deep-Learning Models](#), see page 19 (optional)
- [Activating the Software License](#), see page 20
- [Updating Image Stacks, Results Data, and Protocols](#), see page 22



Note:

- If you are upgrading to version 2.1 of the IN Carta software from version 1.17 or version 1.16, you must uninstall the earlier version of the software before running the installer.
 - If you are upgrading the IN Carta software from a version earlier than version 1.16, contact Molecular Devices Technical Support. See [Obtaining Support on page 6](#) for details.
-

Downloading the Software Installers

You can download the IN Carta software installers from the Molecular Devices Knowledge Base.




Note: If the IN Carta computer is not connected to the Internet, use another computer that has an Internet connection to download the software installers, and then transfer it to the IN Carta computer.

To download the software installers:

1. In a browser, log in to the IN Carta page of the Molecular Devices Knowledge Base at www.meta.moleculardevices.com/links/incarta.



Note:

- If you do not have a Knowledge Base account, click **Create Account**, and follow the on-screen instructions.
- With a computer purchased from Molecular Devices, you can also click  **IN Carta download** on the Windows Desktop to connect to the IN Carta page on the Molecular Devices Knowledge Base.

2. Under **Software Downloads**, click **IN Carta 2.x Software Download**.
3. Download the current version of the IN Carta installer and corresponding checksum file to the **Desktop** folder:
 - INCarta_2.1.xxxxxxx.exe
 - INCarta_2.1.xxxxxxx_checksum.exe
4. Download the current version of the IN Carta Base installer and corresponding checksum file to the **Desktop** folder:
 - INCarta_Base_2.1.xxxxxxx.exe
 - Base_2.1.xxxxxxx_checksum.exe
5. If you are installing the SINAP module, download the IN Carta Models installer and corresponding checksum file to the **Desktop** folder:
 - INCarta_Models_1.4.xxxxxxx.exe
 - Models_1.4.xxxxxxx_checksum.exe
6. Double-click each checksum file to verify the integrity of the installers.



Note: If a security warning appears, click **Run**.

- If the checksum test passes, a success message displays.
- If the checksum test fails, a file corruption error displays. Repeat these steps to download that installer again.

Installing the Software



CAUTION! You must complete the steps described in [Before Installing the IN Carta Software on page 9](#) in order to run the IN Carta software installer. Otherwise, system functionality and performance may be affected.

To install the IN Carta software:

1. Do the following to install INCarta_Base:
 - a. On the Windows Desktop, right-click **INCarta_Base_2.1.xxxxxx.exe** and select **Run as administrator**.
-



Note: If a security warning appears, click **Run**.

- b. Click **Yes**.
 - c. Click **Install**.
 - d. When the installation completes, click **Close**.
 2. Do the following to install IN Carta software.
 - a. On the Windows Desktop, right-click **INCarta_2.1.xxxxxx.exe** and select **Run as administrator**.
-



Note: If a security warning appears, click **Run**.

- b. Review the license agreement.
 - c. If you accept the license agreement, select the **I agree to the license terms and conditions** check box.
 - d. Click **Install**.
 - e. If a message displays about monitor resolution, click **Yes** to continue installing.
 - f. When the installation completes, click **Close**.


Verifying the GPU Driver for the SINAP Module and IN Carta 3D Viewer

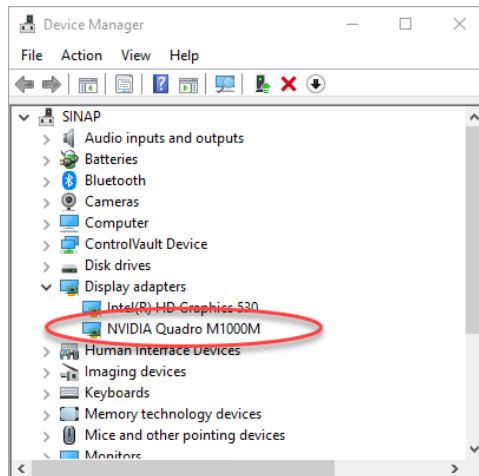


Note: If you did not purchase the SINAP module or the IN Carta 3D Viewer, you can skip this procedure.

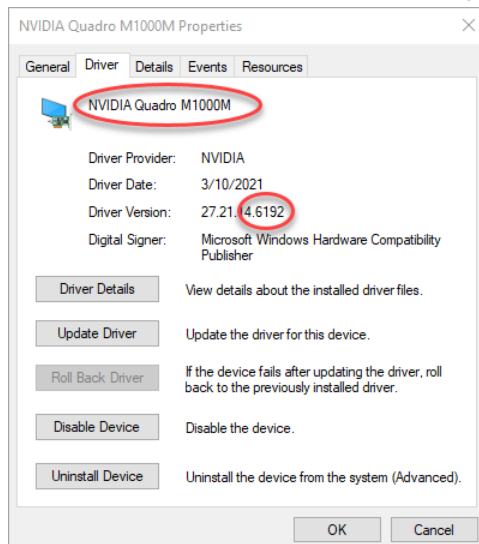
The optional IN Carta SINAP module requires a supported Nvidia graphics processing unit (GPU) with the latest driver and a CUDA compute capability of at least 3.5. For details on determining the CUDA compute capability of your Nvidia GPU, go to developer.nvidia.com/cuda-gpus.

To verify the GPU driver installed in the IN Carta computer:

1. Right-click  **Start**, and select **Device Manager**.
2. Under **Display Adapters**, verify that an Nvidia GPU is listed.



3. Double-click the Nvidia GPU and click the **Driver** tab.
4. Note the GPU model and the last five digits of the driver version.



5. In a browser, go to www.nvidia.com/download/index.aspx.

6. On the **NVIDIA Download Drivers** page, select your GPU model and operating system.

NVIDIA Driver Downloads

Select from the dropdown list below to identify the appropriate driver for your NVIDIA product.

Product Type:

Product Series:

Product:

Operating System:

Download Type: ?

Language:



Note:

- If a **Download Type** drop-down appears, select **Production Branch/Studio** or **Studio Driver (SD)**.
- With a Tesla series GPU, in the **CUDA Toolkit** drop-down, select **11.0**.

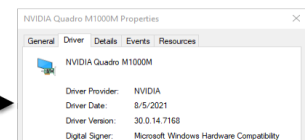
7. Click **Search**.
8. Check the version of the driver that displays and do one of the following:
- If the displayed driver is the same as your installed driver, no action is required. You can skip the rest of this procedure.
 - If the displayed driver is a later version than your installed driver, continue this procedure to download and install the latest driver.



Note: The driver version number appears differently on the Nvidia website and in the Device Manager Properties dialog. The decimal point is shown in different place. For example, the driver shown on the website as version 471.68 appears in the Device Manager Properties dialog as version 4.7168.



Version 471.68
is the same driver as
Version 4.7168



9. Click **Download** and follow the on-screen prompts to download the driver.
10. In Windows File Explorer, browse to the location of the downloaded driver.
11. Double-click the driver and follow the on-screen prompts to install it.

Installing the SINAP Deep-Learning Models

The optional SINAP module requires the installation of the deep-learning models.



Note: If you did not purchase the SINAP module, you can skip this procedure.

To install the deep-learning models for the SINAP module:

1. On the Windows Desktop, right-click **INCarta_Models_1.4.xxxx.exe** and select **Run as administrator**.



Note: If a security warning appears, click **Run**.

2. Click **Install**.
3. When the installation completes, click **Close**.

Activating the Software License

To activate the IN Carta software license, you need the following:

- IN Carta License Activation utility, which is installed on your computer along with the IN Carta software.
- Activation Code, which you received from Molecular Devices.



Note:

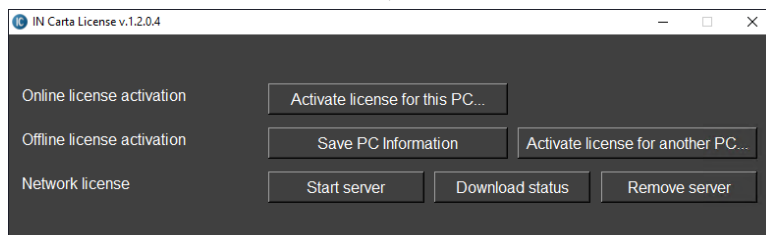
- If you do not have your activation code, contact Molecular Devices Technical Support. See [Obtaining Support on page 6](#) for details.
- An activation code can only be used to activate a license one time. If you need to re-use a code, it must be reset by Molecular Devices Technical Support.

Activating the software license requires a computer with an Internet connection to download a license file. If the IN Carta computer is not connected to the Internet, you can use another computer that has an Internet connection to download a license file, and then transfer it to the IN Carta computer.

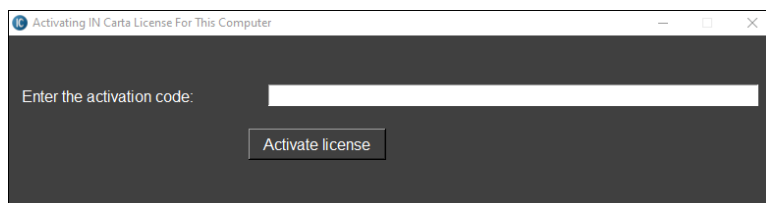
Activating on an IN Carta Computer Connected to the Internet

To activate the software license on an IN Carta computer with an Internet connection:

1. On the IN Carta computer, in Windows File Explorer, browse to the `C:\Program Files (x86)\INCarta_License` folder.
2. Right-click **INCartaLicensing.exe** and select **Run as administrator**.
3. Next to **Online license activation**, click **Activate license for this PC**.



4. In the **Enter the activation code** field, enter the activation code you received from Molecular Devices.

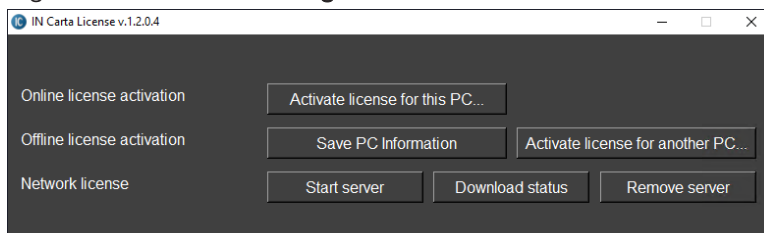


5. Click **Activate license**.

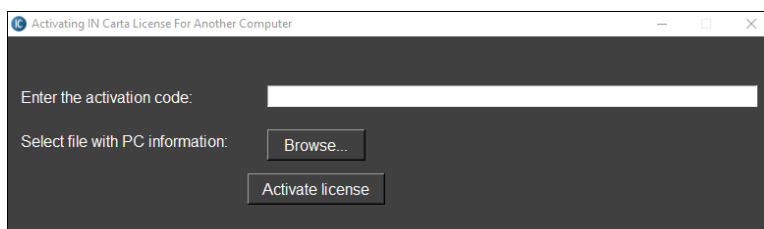
Activating on an IN Carta Computer Not Connected to the Internet

To activate the software license with the IN Carta computer without an Internet Connection:

1. On the IN Carta computer, in Windows File Explorer, browse to the `C:\Program Files (x86)\INCarta_License` folder.
2. Right-click **INCartaLicensing.exe** and select **Run as administrator**.



3. Next to **Offline license activation**, click **Save PC Information**.
The IN Carta License Activation utility creates an **INCartaLicenseActivation** folder on the Windows Desktop.
4. Transfer this folder to the **Desktop** folder of a computer connected to the Internet.
5. On the computer connected to the Internet, right-click **INCartaLicensing.exe** and select **Run as administrator**.
6. Next to **Offline license activation**, click **Activate license for another PC**.
7. In the **Enter the activation code** field, enter the activation code you received from Molecular Devices.



8. Next to **Select file with PC information**, click **Browse**.
9. Navigate to the **INCartaLicenseActivation** folder on the **Desktop**.
10. Select **PCInfo.lcf**.
11. Click **Open**.
12. Click **Activate license**.
The IN Carta License Activation utility creates a License file (**license.lic**) the **INCartaLicenseActivation** folder on the **Desktop**.
13. Transfer the License folder back to the `C:\Program Files (x86)\INCarta_License` folder on the IN Carta computer.

Updating Image Stacks, Results Data, and Protocols

After upgrading from version 1.17 or version 1.16 of the IN Carta software, you must use the IN Carta Converter Tool to update existing results data and protocols to the new format. You are not required to converting existing image stacks. The software will convert existing image stacks to the new format on-the-fly when you open them, but you can also update them all at once with the IN Carta Converter Tool.

With some exceptions (noted below), the IN Carta Converter Tool does the following:

- Converts image stacks from XDCE format (for version 1.x) to JDCE format (for version 2.x).
- Converts protocols created for the **Mono-nucleated Cells**, **Time-lapse Analysis**, and **2D-Analysis** applications to the new format for use with the new **Flexi-Protocol** application.
- Converts results data to the new format.



Note:

- Converted protocols can no longer be opened in version 1.x of the IN Carta software.
 - You must import all converted protocols in version 2.1 of the IN Carta software before they display in the Protocol Editor Table. See the *IN Carta Help* for details on importing protocols.
 - Protocols created in version 1.x for the **Custom Module Editor 2D** and **Custom Module Editor 3D** applications do not require conversion. However, you must import them in version 2.1 of the IN Carta software. See the *IN Carta Help* for details on importing protocols.
 - Protocols created in version 1.x for the **Volumetric Analysis** application are no longer supported in version 2.1 of the IN Carta software and cannot be converted.
 - Measures in protocols created in version 1.x that require the calculation of more than one target (for example, intensity ratio of nucleus to cytoplasm) cannot be converted and are excluded from the converted protocol.
 - Classifiers in protocols created in version 1.x that include a measure that cannot be converted are excluded from the converted protocol.
-


Downloading the IN Carta Converter Tool

To download the IN Carta Converter Tool:

1. In a browser, log in to the IN Carta page of the Molecular Devices Knowledge Base at www.meta.moleculardevices.com/links/incarta.



Note:

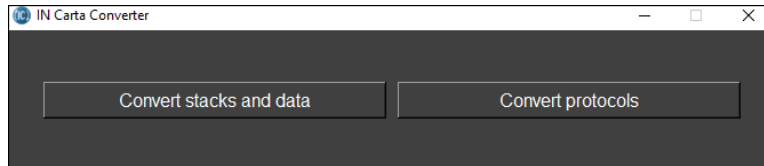
- If you do not have a Knowledge Base account, click **Create Account**, and follow the on-screen instructions.
 - With a computer purchased from Molecular Devices, you can also click  **IN Carta download** on the Windows Desktop to connect to the IN Carta page on the Molecular Devices Knowledge Base.
-

2. Under **Software Downloads**, click **IN Carta 2.x Software Download**.
3. Download `INCartaConverter.exe` to the **Desktop** folder.

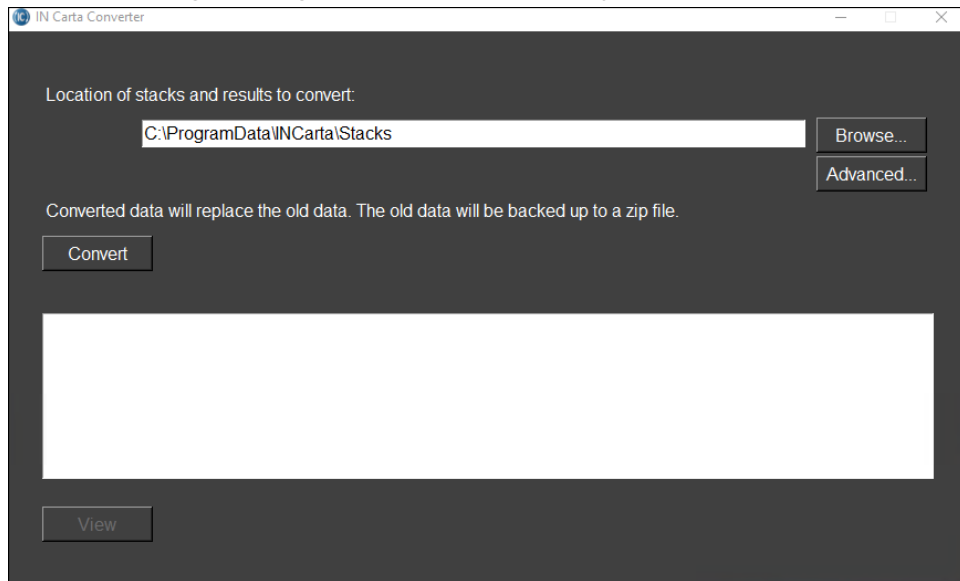
Updating Image Stacks and Results Data

To update image stacks and results data to the current format:

1. On the Windows Desktop, double-click **INCartaConverter.exe**.
2. Click **Convert stacks and data**.



3. In the **Location of stacks and results to convert** field, click **Browse** and browse to select the folder containing the image stacks and results data you want to convert.



4. Click **Select Folder**.
5. If needed, click **Advanced** to set the following options:
 - Select the **Archive all images** check box to create a ZIP file for each stack. The ZIP file contains the XDCE file and all images for the stack. This enables you to easily open the original stack (in version 1.x), if needed. Note that archiving all images can significantly slow down the conversion process and substantially increase a disk space usage.
 - Select the **Backup the results folder before conversion** check box to create a ZIP file for each Data folder. This enables you to easily open the original results data (in version 1.x), if needed. The update will take noticeably longer with this option selected.



Note: If a ZIP file of the Data folder exists, selecting this check box restores the version 1.x Data folder before conversion. This can be useful if the previous conversion was interrupted, resulting in a corrupted converted Data folder.



6. Click **Convert**.
In the selected folder, all image stacks and results data are converted to the current format.
7. If needed, click **View** to display details of the conversion.

Updating Protocols

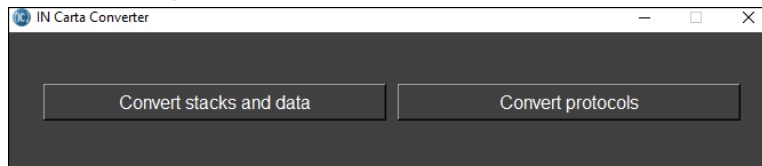
To update protocols to the current format:

1. On the Windows Desktop, double-click **INCartaConverter.exe**.



Note: If a security warning appears, click **Run**.

2. Click **Convert protocols**.

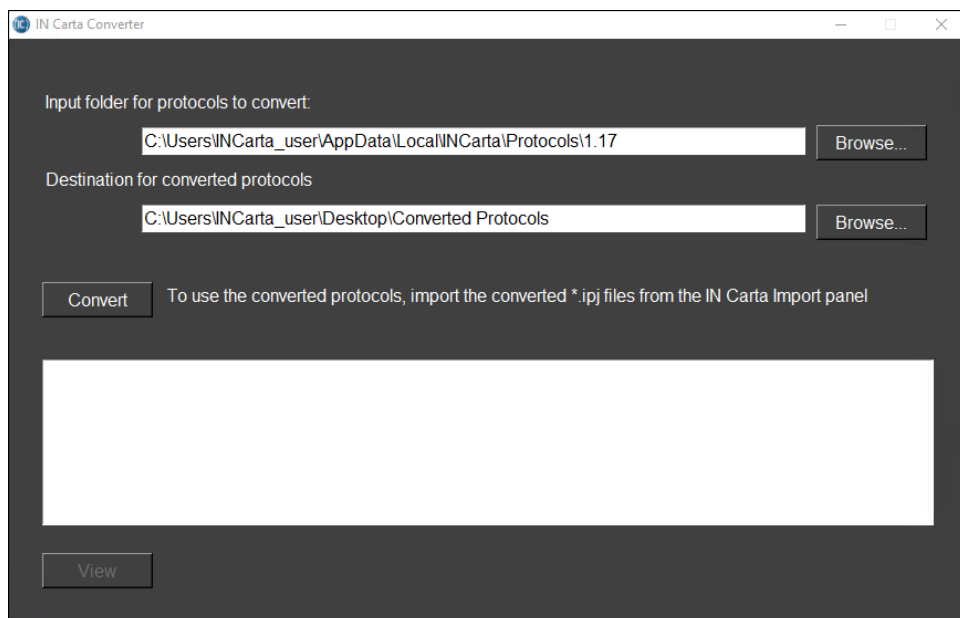


3. In the **Input folder for protocols to convert** field, click **Browse** and browse to select the folder containing the protocols you want to convert.



Note: Existing protocols are stored in the following folders:

- %localappdata%\INCarta\Protocols\1.16
- %localappdata%\INCarta\Protocols\1.17



4. Click **Select Folder**.
5. In the **Destination for converted protocols** field, click **Browse** and browse to select the output folder for converted protocols.



Note: The destination folder must not be under %localappdata%\INCarta. We recommend specifying the **Desktop** folder.

6. Click **Select Folder**.
7. Click **Convert**.

In the specified destination folder, protocol files are created in the current format.

You must import updated protocols in the version 2.1 of the IN Carta software before they appear in the Protocol Editor Table. See the *IN Carta Help* for details on importing protocols.



This section describes the procedures required after running the IN Carta software installer:

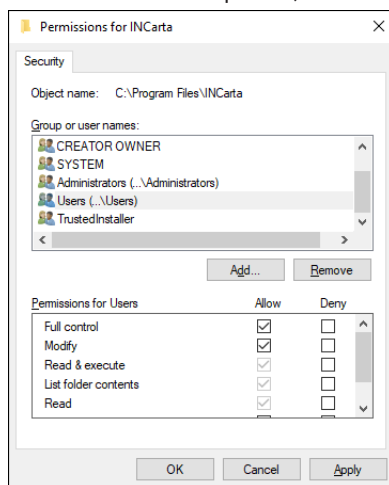
- [Granting Folder Permissions](#), see below
- [Granting Software Permissions](#), see page 29

Granting Folder Permissions

To run the IN Carta software, you must grant the following folder permissions for the user account where the software is installed.

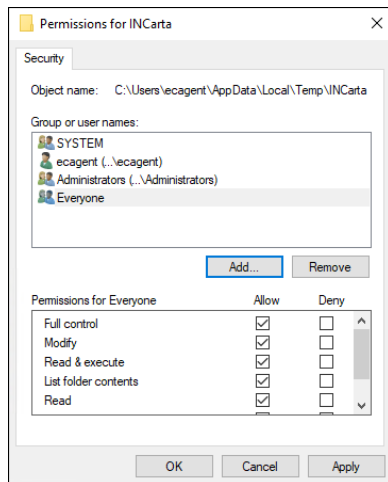
To grant folder permissions:

1. In Windows File Explorer, browse to the `C:\Program Files` folder, and do the following:



- a. Right-click on the **INCarta** folder, and select **Properties**.
- b. Click the **Security** tab.
- c. Click **Edit**.
- d. In the **Group or user names** list, select **Users**.
- e. In the **Permissions for Users** list, in the **Allow** column, select the **Full control** check box.
- f. Click **OK**.
- g. Click **OK**.

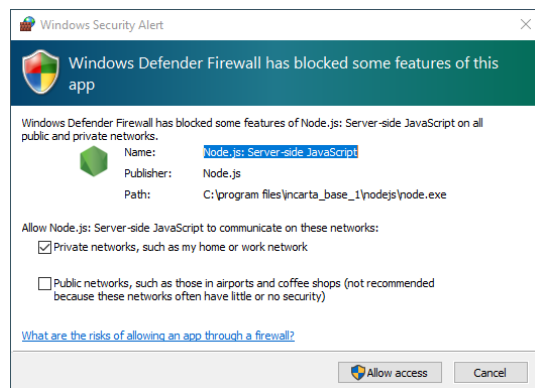
2. In Windows File Explorer, click in the address bar, enter %temp%, and do the following:



- a. Right-click on the **INCarta** folder, and select **Properties**.
- b. Click the **Security** tab.
- c. Click **Edit**.
- d. Click **Add**.
- e. In the **Enter the object names to select** list, enter *Everyone*.
- f. Click **Check Names**.
- g. Click **OK**.
- h. In the **Permissions for Everyone** list, in the **Allow** column, select the **Full control** check box.
- i. Click **OK**.
- j. Click **OK**.

Granting Software Permissions

The first time you run the IN Carta software after installing, one or more Windows Security Alerts may appear.

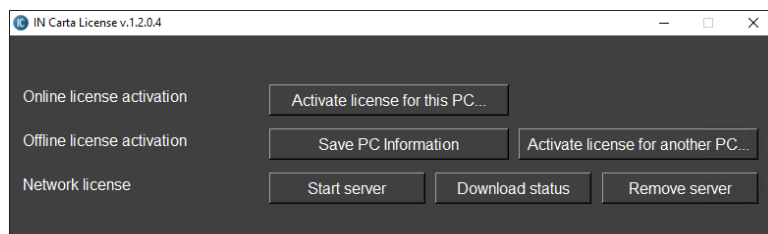


These alerts are requesting that you grant the permissions required to run the software. If an alert appears, click **Allow access**.



The network licensing server and clients must run on the same network.

The server computer must run at all times. To limit the risk of interruption during power loss, we recommend using an uninterruptible power supply (UPS) to provide backup power and power line conditioning for the server computer.



Setting Up a Network Licensing Server

To set up a network licensing server:

1. If you do not plan to install the IN Carta software on the server computer, do the following:
 - a. Install the software on the client computer. See [Installing the Software on page 17](#) for details.
 - b. Create a `C:\Program Files (x86)\INCarta_License` folder on the server computer.
 - c. Copy **INCartaLicensing.exe** from the `C:\Program Files (x86)\INCarta_License` folder on the client computer to the `C:\Program Files (x86)\INCarta_License` folder on the server computer .
2. On the server computer, in Windows File Explorer, browse to the `C:\Program Files (x86)\INCarta_License` folder.
3. Right-click **INCartaLicensing.exe**, and select **Run as administrator**.
4. Activate the (network) license, if not yet activated. See [Activating the Software License on page 20](#) for details.
5. Confirm that the **license.lic** file is present in the `C:\Program Files (x86)\INCarta_License` folder.
6. In the **INCartaLicensing** window, click **Start server**.



Note:

- As part of starting a network licensing server, Java is installed on the server computer.
- If you are unable to start the network licensing server, you may need assistance from your IT professional. Contact Molecular Devices Technical Support to learn more. See [Obtaining Support on page 6](#) for details.

Setting Up a Network Licensing Client



Note: Before you set up a network licensing client, install the IN Carta software on the client computer. See [Installing the Software on page 17](#) for details.

To set up a network licensing client:

1. Copy the **license.lic** file from the `C:\Program Files (x86)\INCarta_License` folder on the server computer to the `C:\Program Files (x86)\INCarta_License` folder on the client computer. Because of permissions, you may have to copy this file first to the desktop of the client computer and then move it to the desktop to the destination folder.
2. Create a `c:\ProgramData\INCarta_License` folder on the client computer.
3. Copy the **nls_address.txt** file from the `c:\ProgramData\INCarta_License` folder on the server computer to the `c:\ProgramData\INCarta_License` folder on the client computer.



Visit us on Spectranet.

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.