



# IN Carta

Image Acquisition and Analysis Software

Version 1.14

## SINAP User Guide

## IN Carta Image Analysis Software SINAP User 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.

Patents: [www.moleculardevices.com/patents](http://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.  
©2021 Molecular Devices, LLC.  
All rights reserved.



# Contents

<b>Chapter 1: IN Carta Image Analysis Software</b>	<b>5</b>
Terminology	5
Obtaining Support	6
About This Guide	6
<b>Chapter 2: IN Carta SINAP</b>	<b>7</b>
Machine Learning	7
AI for Image Analysis	7
IN Carta SINAP	8
<b>Chapter 3: Using IN Carta SINAP</b>	<b>9</b>
Saving Data and Data Storage	9
Launch the Application	10
Create New Protocol with Deep Learning Analysis	11
Refining a Pre-Trained Model	12
AI-Assisted Segmentation	13
Manual Segmentation	14
Using a Training Set to Train a Model	15
<b>Chapter 4: Troubleshooting</b>	<b>17</b>
Update Parameters for Non-Optimal GPUs (if required)	17



# Chapter 1: IN Carta Image Analysis Software

# 1

The IN Carta™ Image Analysis Software provides powerful analytics combined with an 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, you can get meaningful data quickly and reliably.

## 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.

## Terminology

The following table defines the terms and abbreviations used in this guide.

Term	Definition
Annotation	User-created labeling (target/background) of input data to define areas of interest.
Background	Area of an image that is not of interest.
Base Model	One of the pre-trained neural network models provided with IN Carta SINAP for a range of segmentation applications.
Epoch	Single iteration of learning algorithm working through entire training set in an attempt of creating a model.
FOV	Field of view.
Ground Truth	User-annotated input data.
Input data	Images.
Model	Trained deep learning neural network.
ROI	Region of Interest.
Target	Biological structure that is of interest.
Training Set	Set of annotated images used to train a model.

## 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](http://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](http://support.moleculardevices.com).

You can also submit a support request by phone. For regional support contact information, go to [www.moleculardevices.com/contact](http://www.moleculardevices.com/contact).

To expedite support:

- For the IN Carta software, please be prepared to provide the software version and the service tag for the analysis computer. To find the software version, in the IN Carta software, select version in the **System > About** in the top right corner of the window. The service tag is on the label on the top of the analysis computer.
- For the MetaXpress software, please be prepared to provide the system ID number, the software version, and the name of the system owner. To find this information, in the MetaXpress software, select **Help > About MetaXpress**.

### Documentation

Review the product documentation on the Molecular Devices Knowledge Base at [support.moleculardevices.com](http://support.moleculardevices.com). In addition, online Help is available within the IN Carta software.

### Additional Resources

Web-based microscopy courses:

- [www.microscopyu.com](http://www.microscopyu.com)
- [www.ibiology.org/ibioeducation/taking-courses/ibiology-microscopy-short-course.html](http://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](http://www.thermofisher.com/us/en/home/references/molecular-probes-the-handbook.html)

Filter information:

- [www.semrock.com](http://www.semrock.com)
- [www.chroma.com](http://www.chroma.com)
- [www.omegafilters.com](http://www.omegafilters.com)

## About This Guide

This guide is intended for the scientist using the IN Carta software. It provides an overview of IN Carta SINAP and describes how to use IN Carta SINAP to run your methods.

The information in this guide is subject to change without notice. We recommend that you review the guide on the Molecular Devices Knowledge Base at [support.moleculardevices.com](http://support.moleculardevices.com) for the most up-to-date information.

Segmentation is the foundation of the image analysis pipeline, allowing researchers to identify regions of interest in tissue, whole organisms, individual cells, nuclei, and organelles. Through segmentation, researchers can extract information from images to quantitatively compare differences across diverse concentrations, treatments, time, genetics, and so on.

It is crucial to get segmentation right, as it is the foundation for every analysis that follows. Accurate and reliable identification of structures, shapes, and sizes is vital for robust analyses. Errors in segmentation will be propagated and multiplied through the rest of the analysis, effectively lowering the assay robustness (less accurate dose curve, inability to detect small modulations of a phenotype, lower Z' for classification experiments). This leads to false-positives and false-negatives, and it ultimately wastes time, money, and resources.

### Machine Learning

Current segmentation algorithms often struggle to deal with non-optimal images where poor contrast, signal variability, or high complexity of biological structures mean that you must compromise your analyses and work with a solution that accommodates some of their data, but not all.

This results in multiple, specialized tools, each suited for an application or data set, which often require training to be used effectively.

The advent of Artificial Intelligence (AI) has improved the situation. The development of computer systems able to perform tasks through machine- and deep-learning that would normally have required human intelligence—such as visual perception, speech recognition, decision-making, and translation between languages—has addressed some of the intractable problems. These solutions still require a large amount of annotated *ground truth* data to create a reliable, accurate training set.

A learning algorithm capable of being taught to detect the biology in which a researcher is interested is required.

IN Carta SINAP is a trainable segmentation module that utilizes both classic machine learning and deep-learning in an intuitive format that produces robust, reliable segmentation across a wide variety of applications.

### AI for Image Analysis

IN Carta SINAP uses machine learning as an aid for creating annotated ground truth training sets for the Deep Learning component of the software.

Deep Learning is an area of AI that uses multi-layered neural networks to mimic the way that the human brain processes information. A network becomes more accurate as it is provided with more ground truth data from which to learn.

With Deep Learning, there is no requirement to define specific features -of interest. A model learns in a similar way to humans, and, as a result, it can outperform traditional segmentation methods in many applications.

## IN Carta SINAP

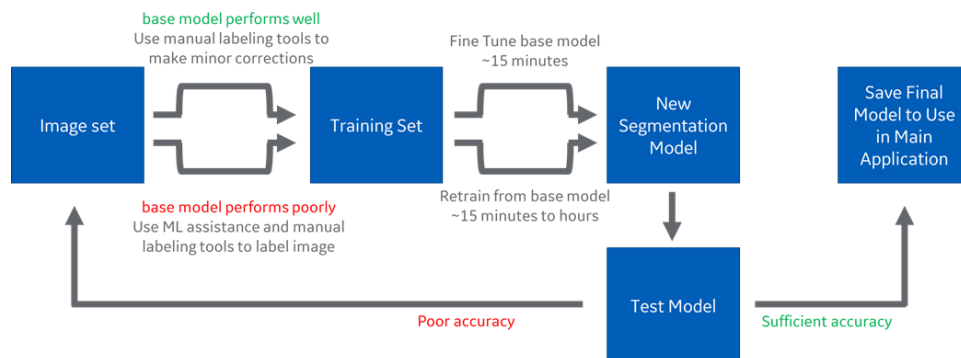
IN Carta SINAP uses pre-trained *base models* to reduce the requirement for large amounts of annotated ground truth data. This saves time and reduces the amount of data required for a model to be accurate. If a model is a good match for the target of interest, the fine tuning will only require a small number of annotated images to be added to a training set.

Where the structure of interest is dissimilar to images used to train a model, the performance of the model may be sub-optimal. In this case, you can re-train an existing model to suit their application.

Re-training a model using IN Carta SINAP adopts an iterative approach, where you add more annotated ground truth data until the model accurately differentiates the target from the background. This takes longer than fine-tuning and requires more data in the training set. Timing and size of training set is highly variable and is dependent on how different the structures are to the base model, and how complex the problem is.

In either case, when you are satisfied with the performance of the model, it can be saved and applied to the full data set.

The saved model will also be available as a pre-trained model for subsequent use in other experiments, as shown in the illustration.



This section contains instructions on how to run your methods using IN Carta SINAP.

- [Saving Data and Data Storage](#), see below
- [Launch the Application](#), see page 10
- [Create New Protocol with Deep Learning Analysis](#), see page 11
- [Refining a Pre-Trained Model](#), see page 12
- [AI-Assisted Segmentation](#), see page 13
- [Manual Segmentation](#), see page 14
- [Using a Training Set to Train a Model](#), see page 15

### **Saving Data and Data Storage**

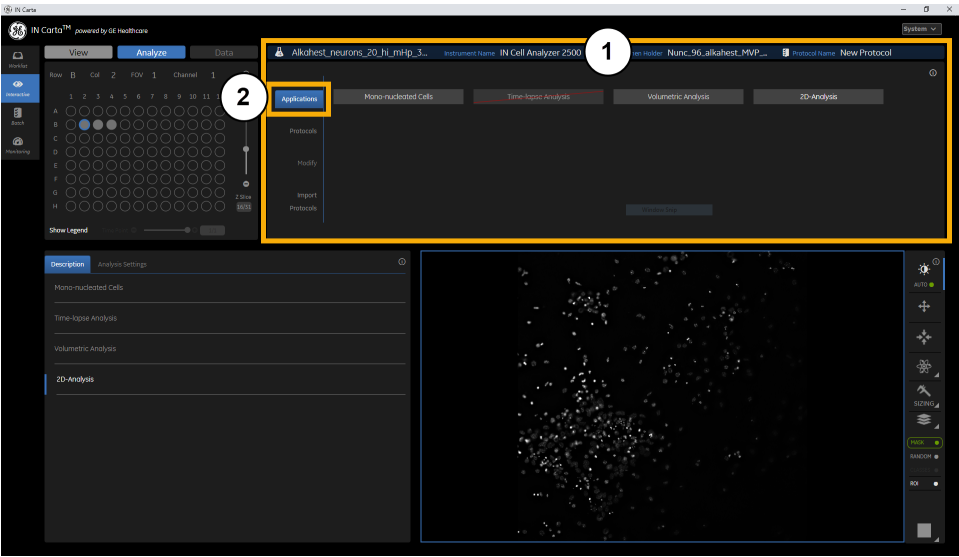
Consider the following when working with IN Carta SINAP:

- It is not possible to delete or overwrite a model once it has been saved.
- Upon editing and saving, a new model is created with a unique identifier.
- Training set data is stored in a temporary folder while the SINAP application is being used.
- Each time the SINAP application is started, all previously stored training-related data is deleted.
- Closing the SINAP application without saving a model means that any changes made to the model will be lost.

## Launch the Application

To launch the application:

From the **Applications Tab** in the **Protocol Editor Table**, click **2D-Analysis** to start the segmentation work flow.



Part	Name
1	Protocol editor table
2	Applications tab

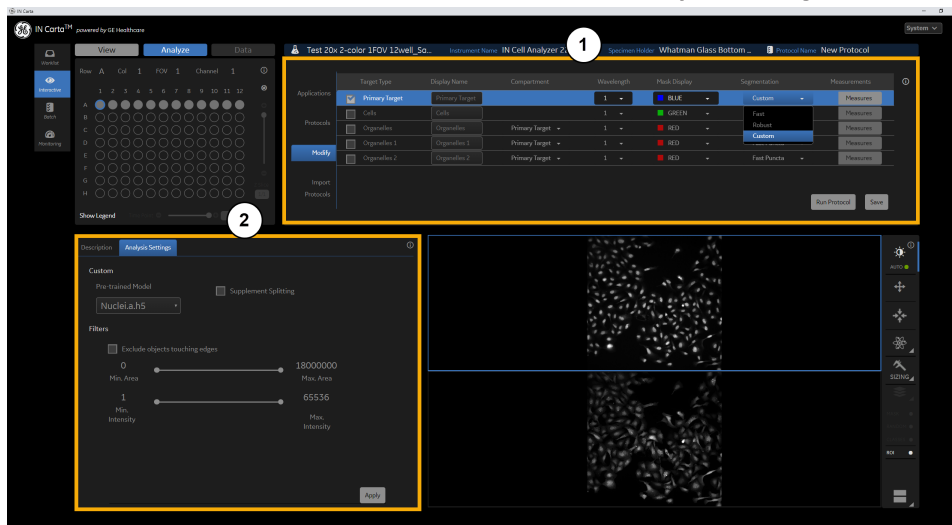
To use a pre-trained model with no further input or training, continue to [Create New Protocol with Deep Learning Analysis on page 11](#).

To refine or create a new model, proceed to [Refining a Pre-Trained Model on page 12](#).


## Create New Protocol with Deep Learning Analysis

To create a new protocol with deep learning analysis:

1. Click **New** to create a new 2D Analysis protocol.
2. Select **Custom** in the **Segmentation** drop-down list.
3. Define the **Wavelength** (channel) of interest.
4. Specify the **Display Name** (for example, nuclei, puncta, mitochondria, other structure).
5. Select a **Pre-trained Model** from the drop-down list in **Analysis Settings**.



Part	Name
1	Protocol editor table
2	Analysis settings

**\* Tip:** More information on filters and 'Supplement Splitting' is accessible through the information panel. Click the  icon in the **Analysis Settings** panel for details.

6. Click **Apply**.
7. If the segmentation is satisfactory, click **Run Protocol** to run the analysis on the data set. Otherwise, proceed to [Refining a Pre-Trained Model on page 12](#).

## Refining a Pre-Trained Model

### General Considerations for Generating a Training Set

Consider the following when generating a training set:

- Different methods of adding images to training set can be used to generate a training set.
- ROI can be used to specify a properly segmented region within an active image. When ROI is active and **Add to Training Set** is clicked, then only the ROI will be added to the training set. The ROI position can be changed by holding **CTRL** on the keyboard and dragging using the left mouse button.
- Multiple ROIs can be added to a training set for any given FOV.
- Consistent labeling of an image is the most important factor. Poor quality ground truth is the biggest obstacle to quickly training a model.



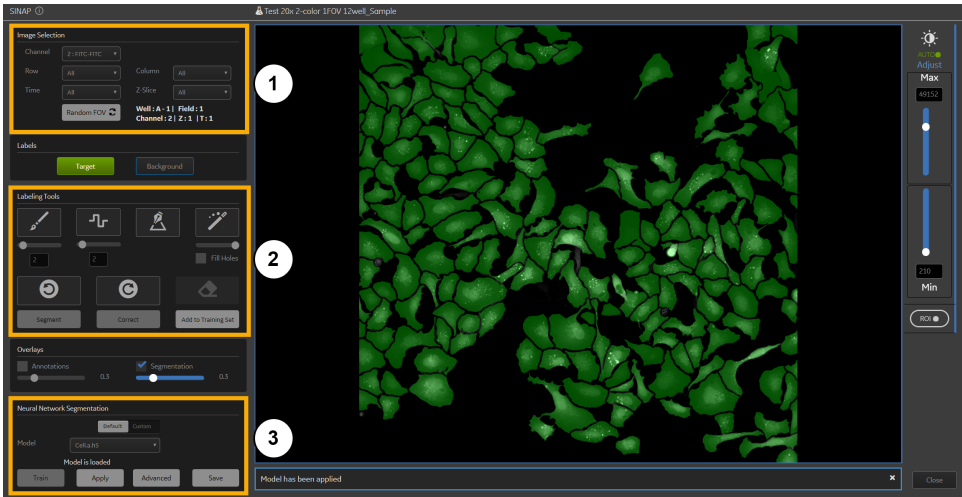
**Tip:** It is more important to label the right pixels than to label many pixels. The most useful information is found at boundaries between target and background. Use the brush tool to precisely annotate these areas.

- Poor image quality, densely packed and overlapping objects, and rare object examples are tougher and typically require more annotated ground truth to be added to a training set.
- The same input images can be annotated differently to train a model to recognize a new target. In this way, models can be trained to recognize multiple targets from the same input images.

### Refining a Pre-Trained Model

To refine a pre-trained model:

1. Click on **SINAP** to launch the IN Carta SINAP application window.
2. Select **Channel**, **Row**, **Column**, **Time**, and **Z-slice** from the drop-down lists in the **Image Selection** panel.



Part	Name
1	Image selection panel
2	Labeling tools panel
3	Neural network segmentation panel

## AI-Assisted Segmentation

### Fine-Tuning

When a pre-trained model produces results that are close to expected and only minor corrections are required, it is simply a matter of fine-tuning the model by correcting areas that have been mislabeled as target or background. This can be done using [Workflow 1: Minor Corrections on page 13](#).

In cases where a model produces sub-optimal segmentation and requires training, refer to [Workflow 2: Major Corrections on page 14](#).

### Workflow 1: Minor Corrections

Do the following if a model produces optimal segmentation and requires only minor corrections:

1. Click **Random FOV** to select a field of view to work on.
2. Adjust brightness/contrast using the control in the right panel.
3. (Optional) Define a region of interest to work on by clicking **ROI** to speed up image segmentation.
4. Select the desired **Model** from the drop-down list in the **Neural Network Segmentation** panel.
5. Click **Apply**.
6. Adjust the opacity of the **Annotation** and **Segmentation** masks using the sliders in the **Overlays** panel.
7. Select the **Label** to be applied (Target or Background) in the **Labeling Tools** panel.
8. Use the **Background** label to train the model to remove false-positive regions.
9. Use the **Target** label to train the model to remove false-negative regions.
10. Select the labeling tool that best fits the objects/areas to be labeled. The following table describes the tool options:

Labeling Tool	Description
Brush	Allows free-hand annotation – the circle size of the cursor corresponds to the size of the brush stroke.
Line	Allows annotating a series of lines using mouse clicks.
Polygon	Allows for annotation of larger regions.
Connected component	Labels image regions with intensities like selected pixels. It can be used for quick annotation of structures of interest. <ol style="list-style-type: none"> <li>Adjust the slider bar to change the sensitivity.</li> <li>Use the <b>Fill Holes</b> check box to include gaps in the annotated area.</li> <li>Add more annotations to correct the segmentation of desired structures.</li> </ol>

11. Click **Correct** to apply the annotations to the segmented image.
12. When the segmentation of the image is satisfactory, the FOV or ROI can be added to the training set by clicking **Add to Training Set** in the **Labeling Tools** panel.

## Workflow 2: Major Corrections

Do the following if a model produces sub-optimal segmentation and requires more ground truth images to be added for training:

1. Click **Random FOV** to select a field of view to work on.
2. Adjust brightness/contrast using the control in the right panel.
3. Using the **Labeling Tools** described in step 9 of [Workflow 1: Minor Corrections on page 13](#), annotate example regions for Target and Background. ROI may be used to speed up segmentation. ROI must include annotations for both Target and Background.
4. Click **Segment** in the **Labeling Tools** panel to segment the current FOV.



**Note:** A machine learning model is used to predict all unlabeled pixels in an image, assigning them as either Target or Background. This model can be re-used for any new FOV that is loaded or when the ROI is moved to a new position.

- a. If the resulting segmentation mask is not satisfactory, more annotations can be added, followed by clicking **Segment**.
- b. (Optional) Corrections can be added to a segmentation mask as described in steps 6 to 10 of [Workflow 1: Minor Corrections on page 13](#).
5. When the segmentation of the image is satisfactory, the FOV or ROI can be added to the training set by clicking **Add to Training Set** in the **Labeling Tools** panel.

## Manual Segmentation

In cases where applying a segmentation model or using the **Segment** button in the **Labeling Tools** panel yields unsatisfactory results, you may want to train a model by annotating only target structures in single or multiple fields of view. In this case, all unlabeled pixels are assigned as background.

Do the following to perform manual segmentation.

1. Click **Random FOV** to select a field of view to work on.
2. Adjust brightness/contrast using the control in the right panel.
3. Using the **Labeling Tools**, annotate only the structures of interest as **Target**. All unlabeled pixels will be assigned as **Background**.
4. Click **Correct**.
5. Repeat steps 3 and 4, if needed.
6. When the segmentation of the image is satisfactory, the FOV or ROI can be added to the training set by clicking **Add to Training Set** in the **Labeling Tools** panel.


## Using a Training Set to Train a Model

After adding images to the training set, they can be used to train a model.

Do the following to train a module:

1. Select either a **Default** model from the drop-down list or a **Custom** model using file browse dialog in the **Neural Network Segmentation** panel.



**Tip:** Consider adjusting **Advanced** parameters using the options seen in the table below. (More information on advanced parameters is accessible through the information panel – to access click the  icon.

Advanced Parameters	Functionality
Fine tune	Use when the base model is closely related to images in the training set.
Retrain	Use when creating a model for a distinct biological structure.
Number of epochs	Increase to allocate more time to obtain a potentially more robust model.

2. Click **Train** to train the model based on the images in the training set.
3. Test the segmentation on the new FOVs to check performance. If it is satisfactory, click **Save** in the **Neural Network Segmentation** panel. Enter the name of the new model and click **OK**.



**Note:**

- If the segmentation results are not robust enough, then more images need to be added to training set.
- Saved models are stored in the c:\ProgramData\TSMModels folder.

The model can now be accessed from **2D Analysis** in the **Analyze** dashboard.



## Update Parameters for Non-Optimal GPUs (if required)

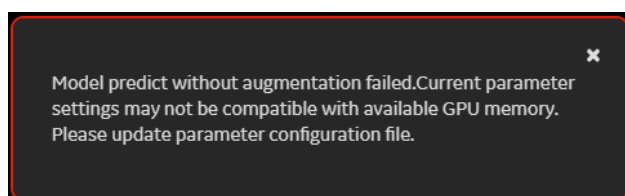
IN Carta SINAP requires an Nvidia GPU card with CUDA compute capability of at least 3.5.

For more information on Nvidia GPUs see [developer.nvidia.com/cuda-gpu](https://developer.nvidia.com/cuda-gpu).



**Note:** Ensure that the latest GPU drivers are installed on the computer.

When using a computer equipped with an Nvidia GPU card with less than 24 GB memory, IN Carta attempts to auto adjust GPU-related settings. However, in some cases, it will be necessary to update the neural net parameters for SINAP to run. The following error message might appear if parameters have not been properly updated.



To update the parameters for deep learning segmentation/training, run the `params_config.bat` file located in `c:\Program Files\INCarta`.

The table below shows default and recommended settings for parameter configuration. Follow the values in the table to adjust parameters for smaller GPU cards. Upon completion, SINAP should now run without displaying an error code.

Parameter	Default Value	Recommended Setting for Smaller GPU
Train Crop Size	1024	512 (if fails, set to 256)
Train Batch Size	2	1
Number of blocks X	1	2 (if fails, set to 4)
Number of blocks Y	1	2 (if fails, set to 4)

## Contact Us

Phone: [+1-800-635-5577](tel:+1-800-635-5577)  
Web: [moleculardevices.com](http://moleculardevices.com)  
Email: [info@moldev.com](mailto: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/patents](http://www.moleculardevices.com/patents)  
FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

©2021 Molecular Devices, LLC.  
All rights reserved.  
5087400 A

