



# IN Carta

Image Acquisition and Analysis Software

Version 1.14

## INPut User Guide

## IN Carta Image Analysis Software INPut 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>  |
| Obtaining Support .....                                   | 6         |
| About This Guide .....                                    | 6         |
| <b>Chapter 2: IN Carta INPut</b> .....                    | <b>7</b>  |
| INPut Main Window .....                                   | 7         |
| <b>Chapter 3: Using INPut to Convert Files</b> .....      | <b>9</b>  |
| Converting DeltaVision/OMX Files .....                    | 9         |
| Converting of PerkinElmer Operetta/Columbus Files .....   | 14        |
| <b>Chapter 4: Using the INPut Annotation Editor</b> ..... | <b>17</b> |
| INPut Annotation Editor Main Window .....                 | 18        |
| Accessing the INPut Annotation Editor .....               | 19        |
| INPut Annotation Editor Options .....                     | 20        |



# 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.

## 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 describes how to use the IN Carta INPut tool to convert image data sets into a format compatible with IN Carta software.

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.

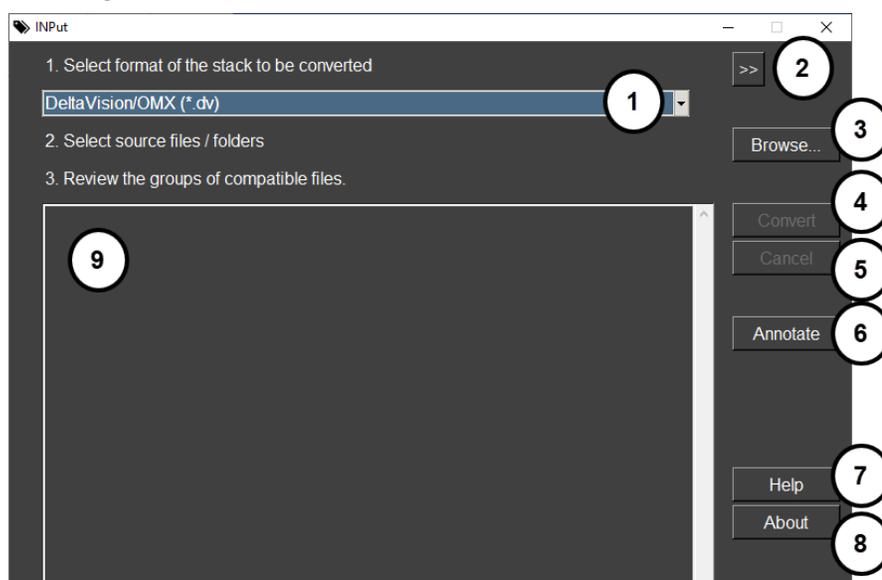
INPut is a tool for converting image data sets into a format compatible with IN Carta analysis software. INPut also allows users to annotate .xdce stacks acquired on IN Cell instrumentation or generated by the INPut converter.

INPut accommodates data sets acquired on or exported from:

- DeltaVision or DeltaVision OMX
- PerkinElmer Operetta/Opera/Columbus

### INPut Main Window

The image below shows the initial view of the INPut main window.



| Part | Name                   | Description                                    |
|------|------------------------|--|
| 1    | Format drop-down list  | Allows selection of a format to convert.       |
| 2    | Expand/Collapse button | Expands/collapses image conversion log window. |
| 3    | <b>Browse</b> button   | Opens file/folder browsing dialog.             |
| 4    | <b>Convert</b> button  | Converts selected files into .xdce stack(s).   |
| 5    | <b>Cancel</b> button   | Cancels current conversion operation.          |
| 6    | <b>Annotate</b> button | Opens INPut: Annotation Editor.                |
| 7    | <b>Help</b> button     | Opens quick guide.                             |
| 8    | <b>About</b> button    | Opens information about software version.      |
| 9    | Review panel           | Displays information during review step.       |



This chapter contains the following instructions for converting files into xdc format:

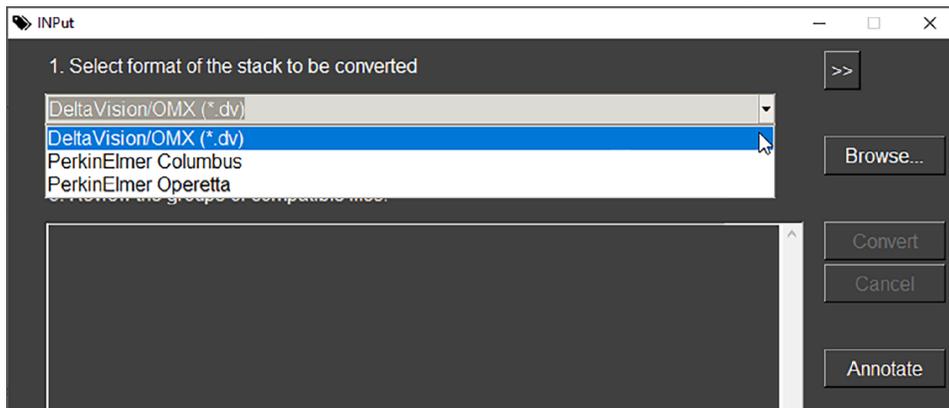
- [Converting DeltaVision/OMX Files, see below](#)
- [Converting of PerkinElmer Operetta/Columbus Files, see page 14](#)

## Converting DeltaVision/OMX Files

### Select Files

To select files for conversion:

1. In the main window, select DeltaVision/OMX from the format drop-down list.



2. Click **Browse**.

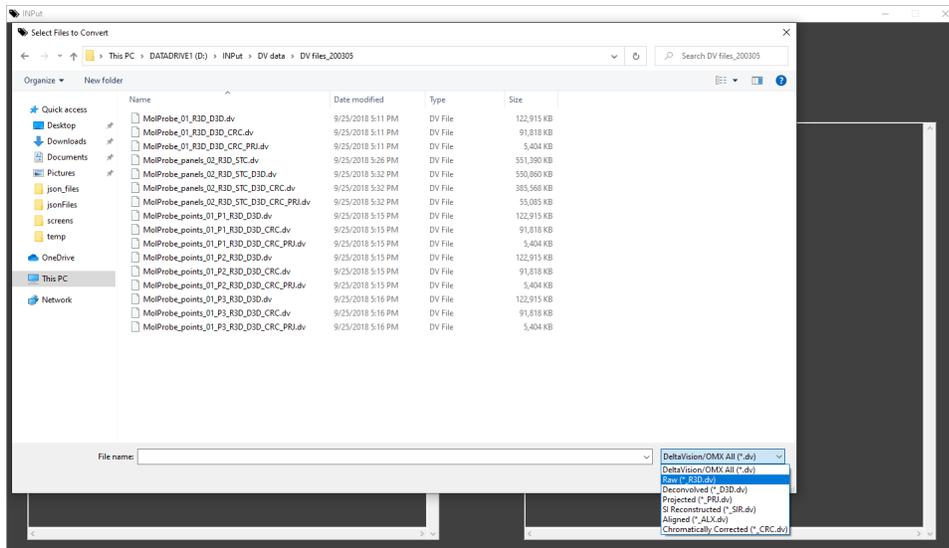
Files with one of the following file extensions appear in the Select Files to Convert dialog.

| File Extension          | Last Post-Processing Step |
|-------------------------|---------------------------|
| DeltaVision/OMX All     | *.dv                      |
| Raw                     | *_R3D.dv                  |
| Deconvolved             | *_D3D.dv                  |
| Projected               | *_PRJ.dv                  |
| SI Reconstructed        | *_SIR.dv                  |
| Aligned                 | *_ALX.dv                  |
| Chromatically Corrected | *_CRC.dv                  |

3. Select one or more files for conversion.



**Note:** Files can be filtered by file extension for display based on last post-processing step.

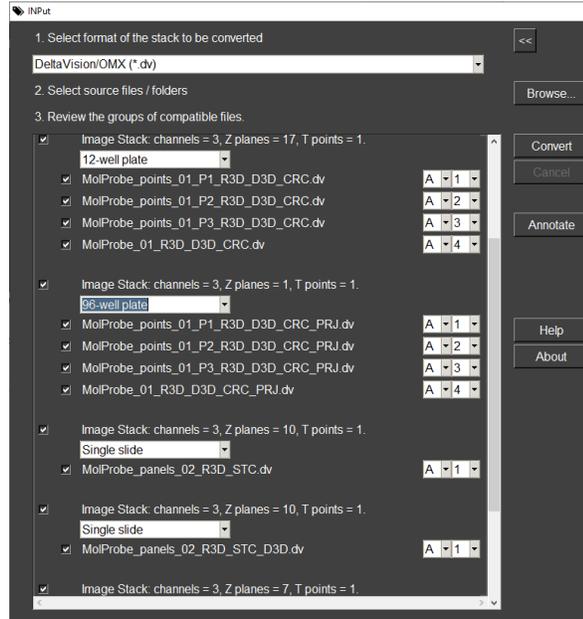


**Tip:** Hold **CTRL** while clicking to select multiple individual files. Hold **SHIFT** while clicking to select a range of files.

4. Click **Open** to import selected files for review.

## Review

During the review step the software displays groups of compatible files.



Files are grouped into plates based on dimensionality. Default plate dimensions depend on the number of files assigned to a plate. All files assigned to one plate have the same:

- magnification/pixel size
- image size
- number of channels
- number of z planes
- number of time-points

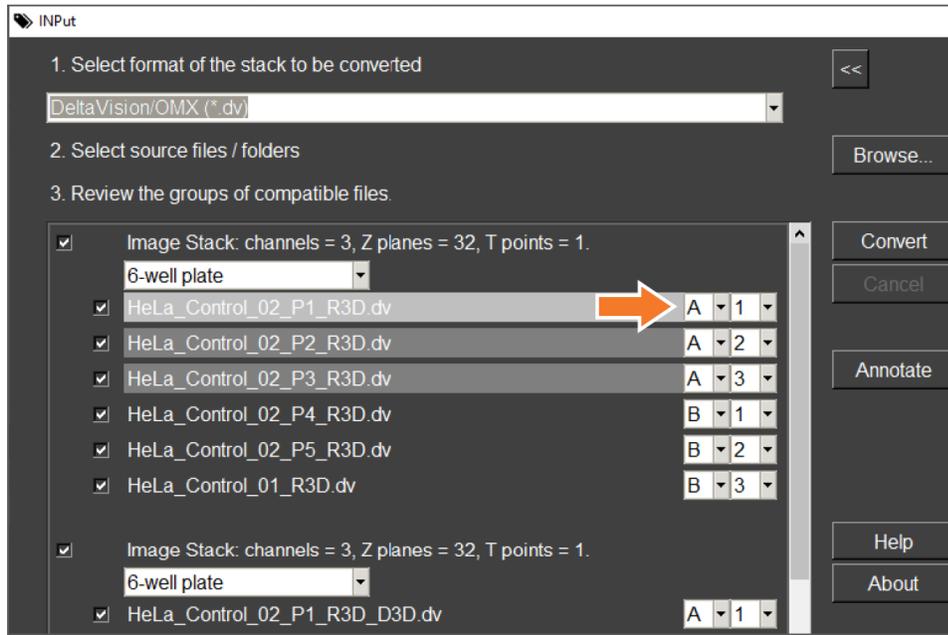
It is possible to modify the plate format or create a virtual plate format for images from a slide. The following table describes the possible actions during the review step.

| Action   | Description   |
|--|---|
| Display ToolTip with <ul style="list-style-type: none"> <li>• stack/data set name</li> <li>• image size/magnification</li> </ul> | Hover the cursor over the plate heading in the review panel.  |
| Change plate format  | Modify using drop-down list under the plate heading. Image files will automatically be assigned well information. |
| Exclude plate from conversion  | Clear the check box for the entire plate to the left of the plate heading.  |
| Exclude individual files from conversion   | Clear the check box for the individual file assigned to a plate.  |
| Change well assignment for individual files  | Modify row/column selection using drop-down list to the right of the file name.                                   |
| Change well assignment for a group of files (files will appear as fields within the selected well)                               | Follow the well-assignment procedure for multiple files, below.   |

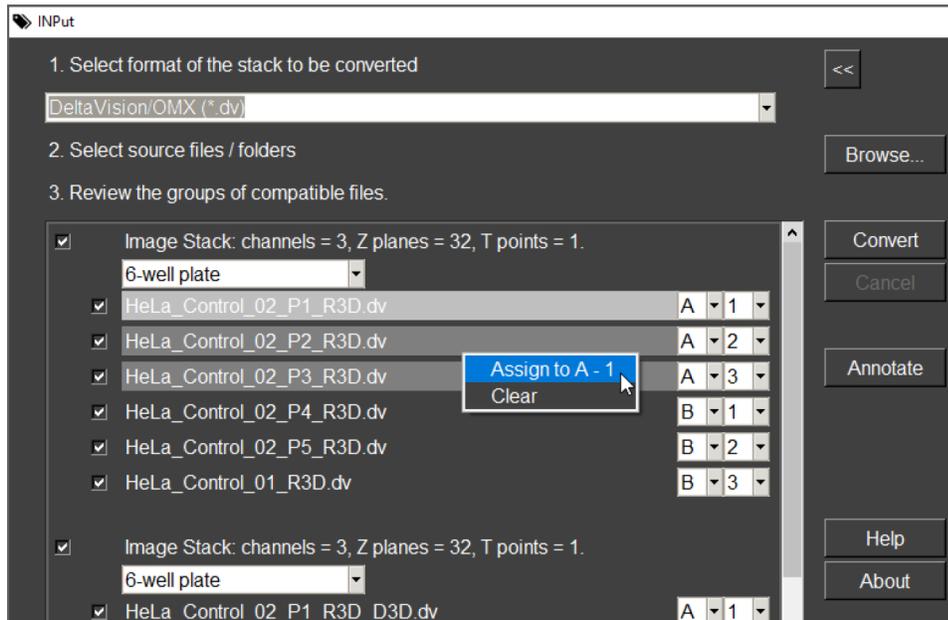
## Modify Well Assignment for Multiple Files

To modify the well assignment for multiple files:

1. Left-click and hold the first file in the target range.
2. Drag the cursor up or down to select a range of files.
3. Release the left mouse button.
4. Change the well assignment of the highlighted file in the selected range manually.



5. Right-click on any file within the selected group.
6. Select **Assign to** to assign all files in the selected group to the well chosen in step 4.



## Convert Selected Files

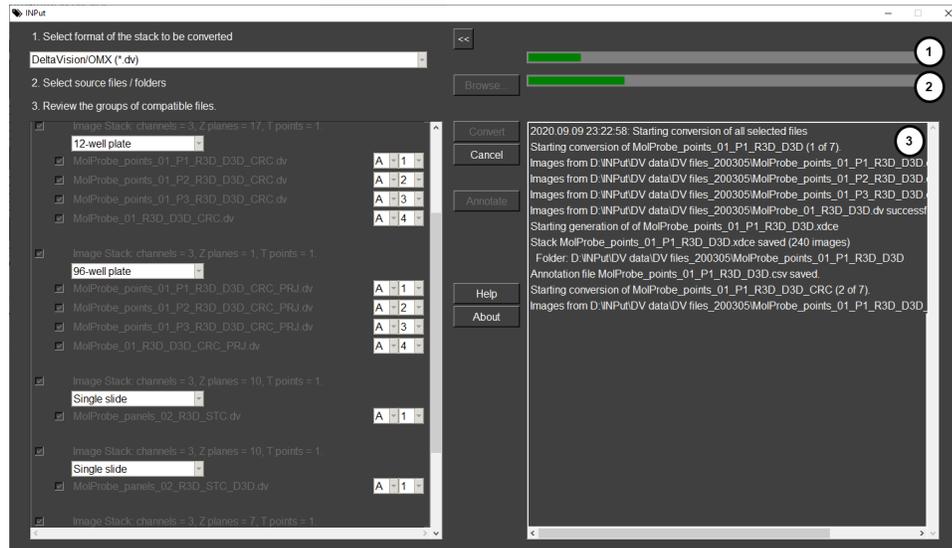
To convert the selected files:

1. Click **Convert**.  
Well assignments for each file are saved in annotation files, which are automatically created for each plate.
2. To cancel the conversion, click **Cancel**.



**Note:** If the conversion is canceled before it is complete, any newly generated files from the incomplete conversion will be automatically deleted.

During the conversion process, the following information is displayed in the image conversion log window.



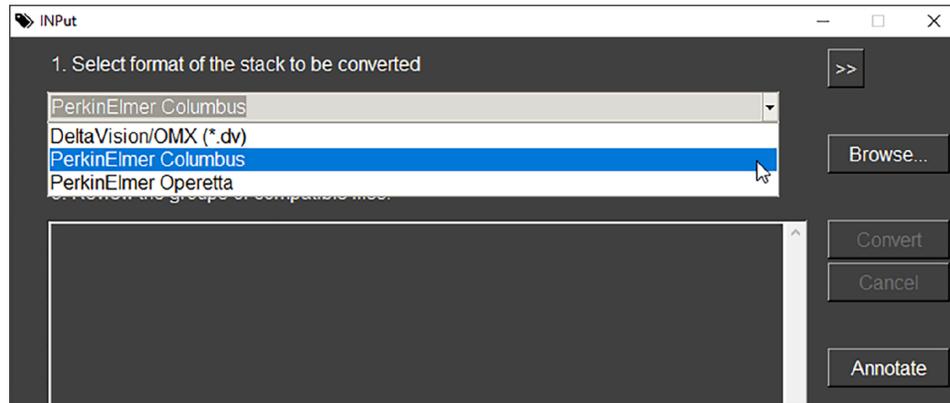
| Part | Name                 | Description   |
|------|----------------------|---|
| 1    | Overall progress bar | Shows overall conversion progress   |
| 2    | Plate progress bar   | Shows progress of current plate conversion  |
| 3    | Log panel            | Shows details about conversion process, including final location path of .xdce stack files. |

## Converting of PerkinElmer Operetta/Columbus Files

### Selecting Files

To select files for conversion:

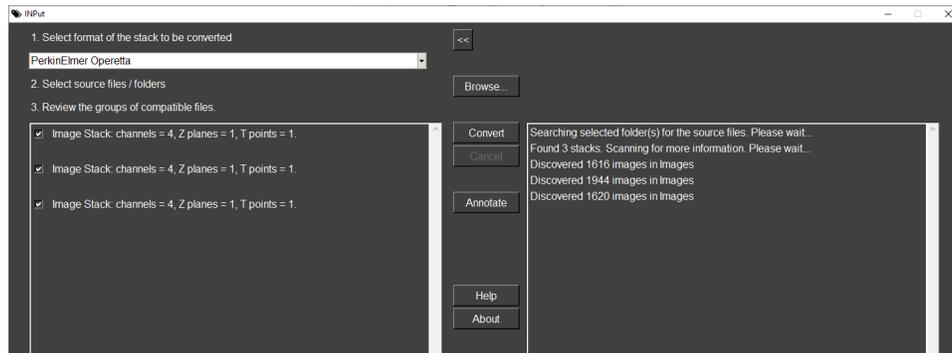
1. In the main window, select PerkinElmer Operetta or PerkinElmer Columbus from the format drop-down list.



2. Click **Browse**.
3. Select a top level folder that contains the image stacks for conversion.

### Review

During the review step the software displays the stacks found in the Review panel.



The following table describes the possible actions during the review step.

| Action   | Description  |
|--|--|
| Display ToolTip with: <ul style="list-style-type: none"> <li>• stack/data set name</li> <li>• image size</li> <li>• magnification or pixel size depending on data available</li> </ul> | Hover the cursor over the plate heading in the review panel                  |
| Exclude image stack from conversion  | Clear the check box for the image stack to the left of the stack description |

## Convert Selected Files

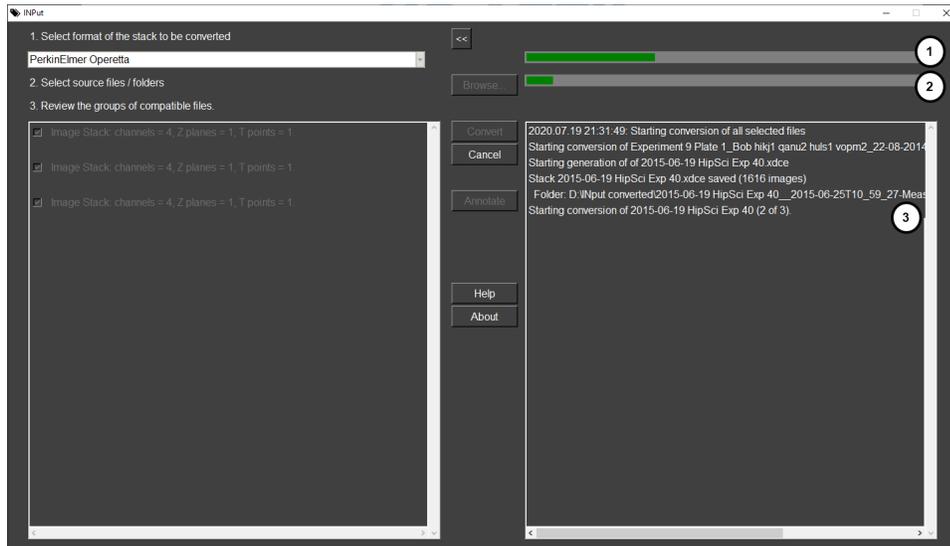
To convert the selected files:

1. Click **Convert**.
2. In the Select Folder dialog, specify the target output folder.
3. Click **Select Folder**.  
INPut initiates conversion of selected data sets into .xdce stacks.
4. To cancel the conversion, click **Cancel**.



**Note:** If the conversion is canceled before it is complete, any newly generated files from the incomplete conversion will be automatically deleted.

During the conversion process, the following information is displayed in the image conversion log window.



| Part | Name                 | Description   |
|------|----------------------|---|
| 1    | Overall progress bar | Shows overall conversion progress   |
| 2    | Plate progress bar   | Shows progress of current plate conversion  |
| 3    | Log panel            | Shows details about conversion process, including final location path of .xdce stack files. |



## Chapter 4: Using the INPut Annotation Editor



# 4

The Annotation Editor tool allows you to create custom plate layouts to store information about each well within a plate. Information about the content of each well can be typed directly into the Annotation Editor window or pasted from a spreadsheet.

This chapter describes the structure and use of the INPut Annotation Editor tool.

## INPut Annotation Editor Main Window

The following image shows the initial view of the INPut: Annotation Editor main window. The following table describes the icons.

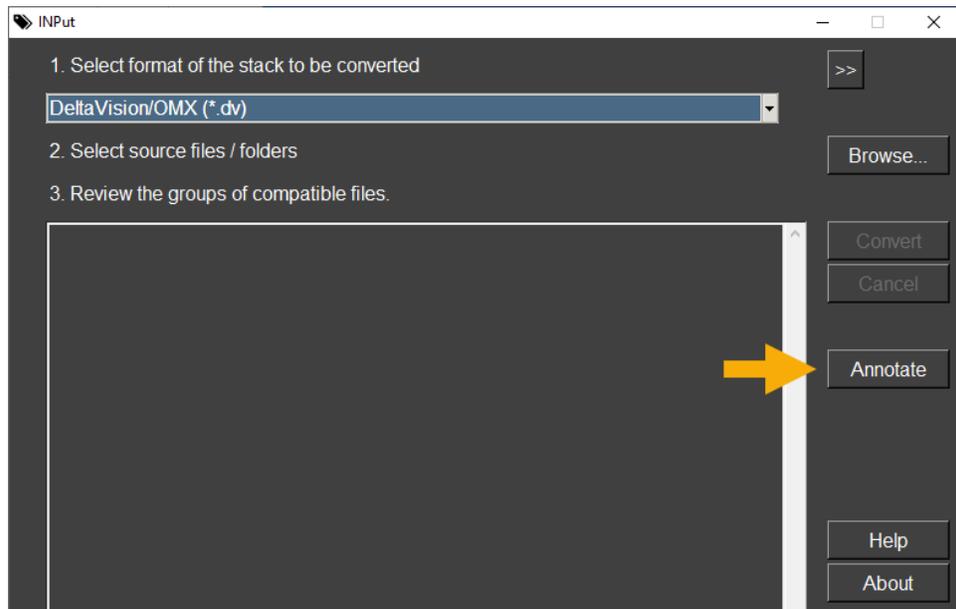


| Icon  | Description   |
|---|---|
|    | Create new annotation file by selecting a predefined sample format or entering a custom format (rows/columns) for the sample.   |
|    | Create new annotation based on an .xdce stack.<br> <b>Note:</b> Plate format will be automatically recognized after the stack is selected. |
|    | Open existing annotation file for review/edit.  |
|    | Save annotation file.   |
|    | Save annotation as a new file.  |
|   | Switch to plate view.<br> <b>Note:</b> Icon visible in list view only.   |
|  | Switch to list view.<br> <b>Note:</b> Icon visible in plate view only.   |
|  | Sort list by row ( <i>default</i> ) or column.  |
|  | Add new annotation type selected from a list of predefined descriptors or custom named.   |
|  | Remove last annotation column added.  |

## Accessing the INPut Annotation Editor

To access the INPut Annotation Editor:

1. In the INPut main window, click **Annotate**.



The INPut: Annotation Editor window opens.

2. Navigate back to the INPut main window by clicking the X in the upper right corner of the INPut: Annotation Editor window.

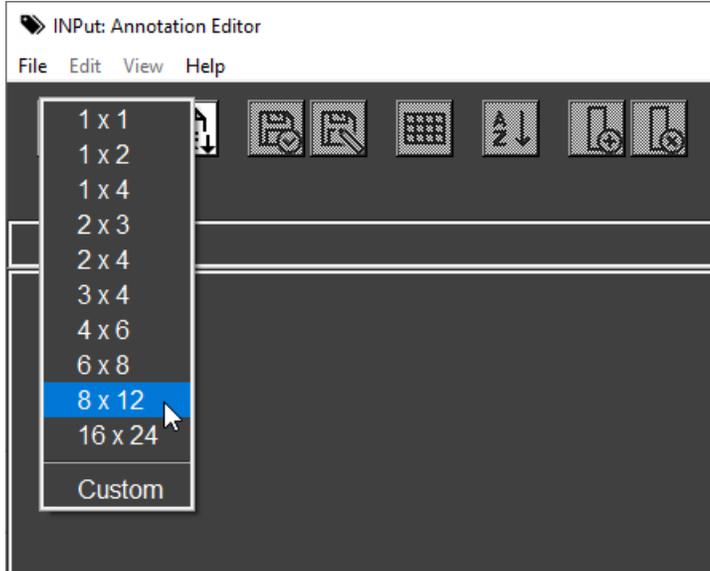


**Note:** Save any open annotation file before closing the Annotation Editor.

## INPut Annotation Editor Options

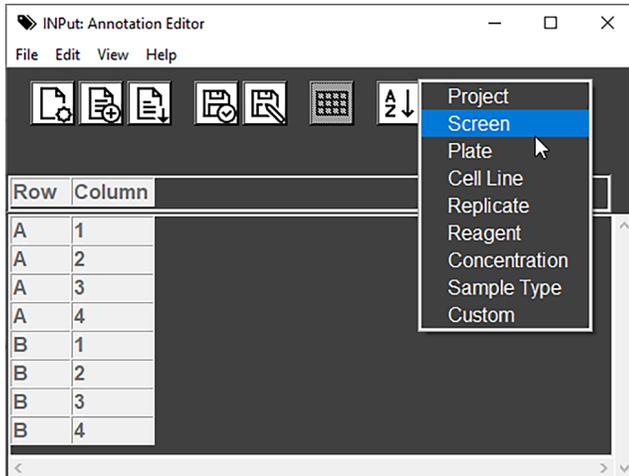
### Predefined Sample Formats

The following images shows the specimen holder sample format choices available for new annotations.



### Annotation Descriptors

The following image shows the annotation descriptors available when adding new columns. Custom columns are named **Custom** by default. You can rename them as needed.



## Editing Content

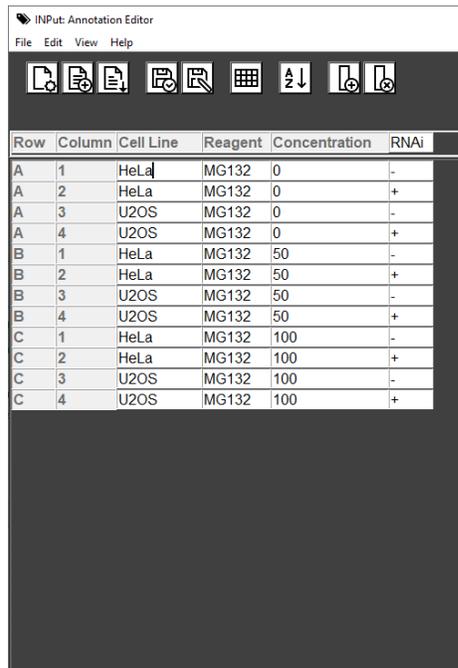
Annotations can be edited by directly typing content in the wells. Edits can also be performed for single or multiple wells according to the following table.

| Edit   | Action  |
|--|---|
| Select multiple wells  | Click and hold while dragging over the target wells.  |
| Repeat information from the top selected well into multiple selected wells below | Right-click selected wells (including the source well) and choose <b>Fill down</b> from the drop-down list. |
| Copy highlighted regions (single or multiple columns)                            | Right-click selected wells and choose <b>Copy region</b> from the drop-down list.                           |
|  | Use the keyboard shortcut <b>CTRL+C</b> .   |
| Paste previously copied regions at cursor position                               | Right-click on one well and choose <b>Paste</b> from the drop-down list.                                    |
|  | Use the keyboard shortcut <b>CTRL+V</b> .   |

## View

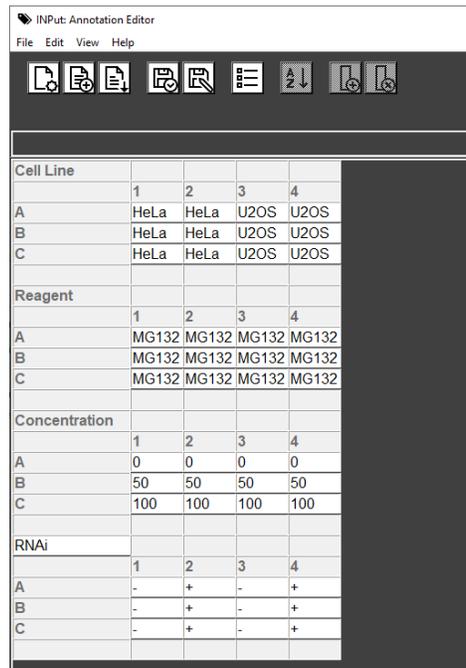
In list view, descriptors are shown as column headings. Sort the data in list view by clicking the sort icon () and choosing column or row.

In plate view, each descriptor is shown in a sample layout.



| Row | Column | Cell Line | Reagent | Concentration | RNAi |
|-----|--------|-----------|---------|---------------|------|
| A   | 1      | HeLa      | MG132   | 0             | -    |
| A   | 2      | HeLa      | MG132   | 0             | +    |
| A   | 3      | U2OS      | MG132   | 0             | -    |
| A   | 4      | U2OS      | MG132   | 0             | +    |
| B   | 1      | HeLa      | MG132   | 50            | -    |
| B   | 2      | HeLa      | MG132   | 50            | +    |
| B   | 3      | U2OS      | MG132   | 50            | -    |
| B   | 4      | U2OS      | MG132   | 50            | +    |
| C   | 1      | HeLa      | MG132   | 100           | -    |
| C   | 2      | HeLa      | MG132   | 100           | +    |
| C   | 3      | U2OS      | MG132   | 100           | -    |
| C   | 4      | U2OS      | MG132   | 100           | +    |

List View (default)



| Cell Line     | 1     | 2     | 3     | 4     |
|---------------|-------|-------|-------|-------|
| A             | HeLa  | HeLa  | U2OS  | U2OS  |
| B             | HeLa  | HeLa  | U2OS  | U2OS  |
| C             | HeLa  | HeLa  | U2OS  | U2OS  |
| Reagent       |       |       |       |       |
| A             | MG132 | MG132 | MG132 | MG132 |
| B             | MG132 | MG132 | MG132 | MG132 |
| C             | MG132 | MG132 | MG132 | MG132 |
| Concentration |       |       |       |       |
| A             | 0     | 0     | 0     | 0     |
| B             | 50    | 50    | 50    | 50    |
| C             | 100   | 100   | 100   | 100   |
| RNAi          |       |       |       |       |
| A             | -     | +     | -     | +     |
| B             | -     | +     | -     | +     |
| C             | -     | +     | -     | +     |

Plate View

## 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.