


# IN Carta Image Analysis Software

## QUICK REFERENCE GUIDE

# Selecting & Viewing Image Data

## Selecting Image Data

1. Click the Worklist  icon.
2. Locate data using **Search** or sort using column headers. To populate the **Worklist** or change directory, click **Browse** and navigate to a parent folder that contains image data and click **OK**.
3. Select image data.

## Navigating and Viewing Image Data

The software includes an image viewer for data exploration. Here is an overview of common visualization tools:

### Setting Up ViewPorts

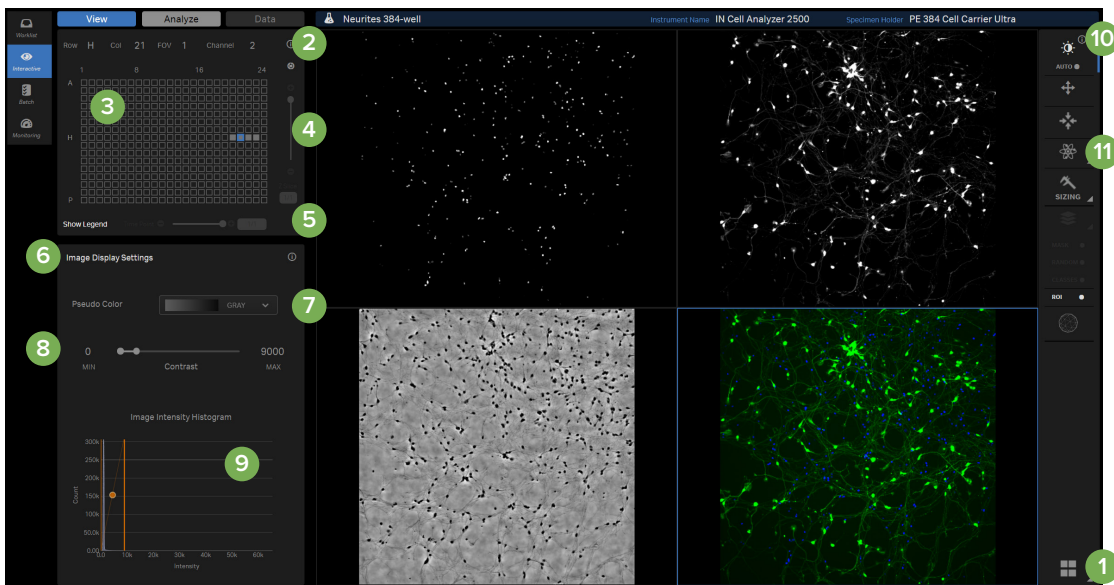
- Click the **ViewPort Layout (1)** icon to change number of images displayed and their orientation. Up to nine channels can be viewed simultaneously.
- To change the image displayed in a ViewPort, click the **ViewPort** to activate it and use drop-downs **(2)** to change field-of-view (FOV) and/or channels.

### Navigating the Dataset

- Select a well in **Plate Map (3)** to display images from that well. Use the mouse wheel to zoom in on the plate map to view and select other FOVs.
- For three-dimensional data, use **Z Slice** slider **(4)** navigate through z.
- For the time lapse data use **Time Point** slider **(5)** to navigate to a time point of interest.

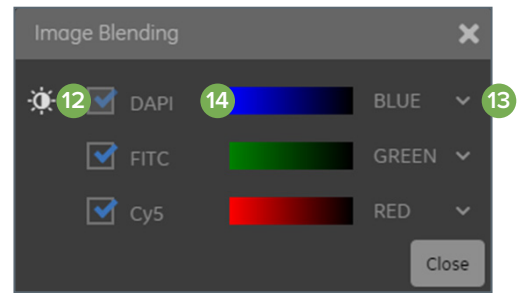
### Adjusting Image Display

- Click **ViewPort** to activate the channels to be adjusted. Use the **Image Display Settings (6)** panel to adjust:
  - Pseudo coloring using the **Pseudo Color (7)** drop-down.
  - Image contrast using either the **Contrast (8)** slider or the **Image Intensity Histogram (9)**.
  - Click the **Auto Contrast (10)** icon to reset contrast to the default min/max values.
  - To manually adjust brightness, click an image, and drag the mouse up and down. To manually adjust contrast, click and drag right and left.



## Creating a Blended Image

1. Click an empty **ViewPort** and click the **Image Blending** (11) icon.
2. Check the box next to each channel (12) to display.
3. Use the **Pseudo Color** drop-down (13) to select a color for each channel.
4. To modify contrast of a blended image, click the name of the wavelength of interest (14). The **Auto Contrast** icon shows which wavelength is selected. Adjust contrast in the **Image Display Settings** (6) panel.



### Contact Us

Phone: +1.800.635.5577  
Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
Email: [info@moldev.com](mailto:info@moldev.com)  
Check our website for a current listing of worldwide distributors.

### Regional Offices

USA and Canada	+1.800.635.5577	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586		

\*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland