

IN Carta software

Select and view image data

Open software and select image data

1. Double-click the IN Carta software  icon.
2. Click the **Worklist**  icon.
3. 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**.
4. Select image data.

Navigate and view image data

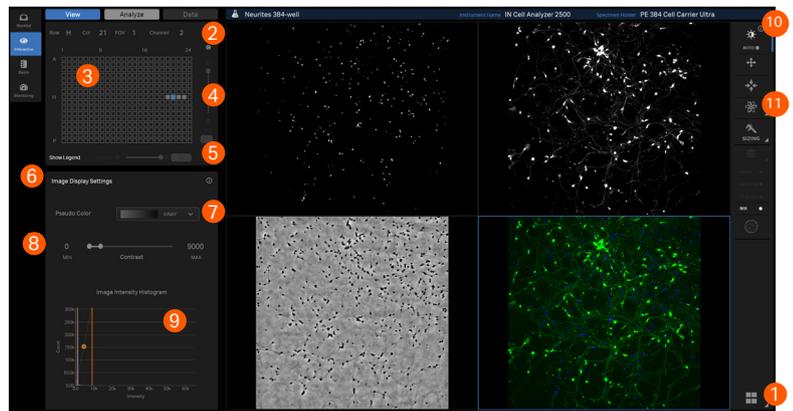
The software includes an image viewer for data exploration. Here is an overview of common visualization tools:

Set up view ports

- Click the **View Port Layout**  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 **View Port**, click the **View Port** to activate it and use drop-downs  to change field-of-view (FOV) and/or channels.

Navigate the dataset

- Select a well in **Plate Map**  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  to navigate through z.
- For the time lapse data use **Time Point** slider  to navigate to a time point of interest.



Adjust image display

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

Create a blended image

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

