

IN Carta Image Analysis Software

QUICK REFERENCE GUIDE

Viewing, Classifying, & Exporting Results

Navigating the Data Table

- Use the **Target (1)** drop-down to select what target data to display.
- Use the **View (2)** drop-down to select the aggregation method:
 - **Single Target:** measures for each segmented object in the selected field-of-view (FOV).
 - **Summary by FOV:** average values for each FOV.
 - **Summary by Well:** average values for each well.
- The **Results (3)** drop-down displays prior results for the current data set. Results are identified by protocol, data set name, and date/time. Select prior results from the drop-down to reload them.

Using the Chart Dashboard Panel

The **Chart Dashboard Panel** is a quick and easy way to display summary or cell by cell data graphically. Here are the most commonly used visualization tools and actions:

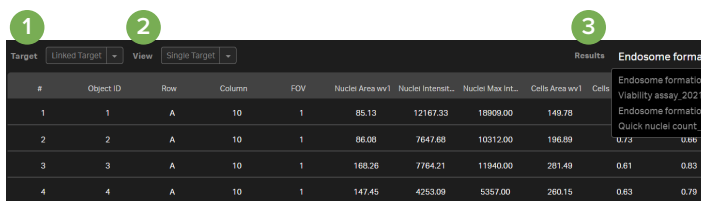
1. Use the **Data Selection (4)** buttons in the **Chart Dashboard** to view **Plate**, **Well** or **FOV** data.




2. Select the **Chart Type (5)** in the Chart Dashboard.

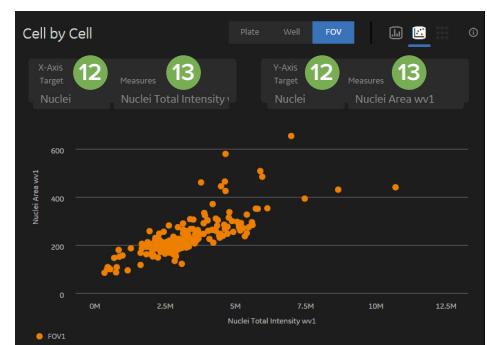
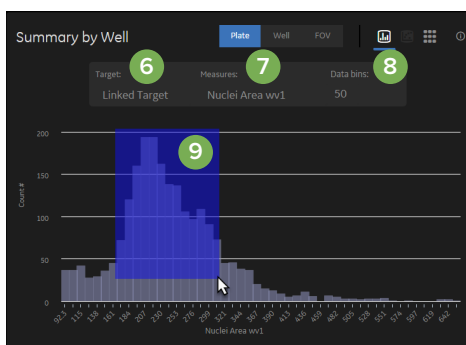
Note: Not all chart types are available with all data selections. See table at bottom for more information.


3. Select chart parameters:

- The **Histogram (5)** displays cell count for a single measure.
 - a. Use **Target (6)** and **Measures (7)** drop-downs to change the data displayed.
 - b. Adjust **Data Bins (8)** as needed. Valid entries are 1 to 99.
 - c. Click and drag over an area of the histogram (9) to zoom in on that portion of the graph (10). Click **Reset zoom (11)** to return to the default display.
- The **Scatter Plot (5)** displays cell by cell data for a single FOV.
 - a. Use **Target (12)** and **Measures (13)** drop-downs to specify data to be plotted on the x- and y-axes.
 - b. Click and drag over an area of the scatter plot to zoom in on a particular area of the graph. Click the **Reset zoom** button to return to the default display.



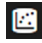
			
Plate	✓		✓
Well	✓		
FOV	✓	✓	



- The **Heat Map**  displays average data for a single measure from each well.
 - Use **Targets/Classes** and **Measures** drop-downs to change the data displayed.
 - Click and drag the arrows at the top and bottom of the **Heat Map Slider** to adjust min/max scaling.

Viewing Linked Data

The IN Carta™ software enables users to interactively connect data points back to a segmented object. This allows you to explore outliers, trends, and perform visual quality control on analyzed data.

1. From the **Chart Dashboard**, set **Data Selection** to **FOV** (14) and **Chart Type** to **Scatter Plot** .
2. Use drop-downs to select **Target** and **Measures** (15) for the x- and y-axes.
3. In the **Data Table**, set **Target** to **Linked Target** (16) and **View** to **Single Target** (17).
4. Verify that the **View Port** is displaying the target identified in the **Scatter Plot** and that **Target Masks** are turned on (18).
5. Click on a point in the **Scatter Plot** (19), cell in the image (20), or row in the **Data Table** (21) to highlight corresponding data for that particular object.


Classifying Data

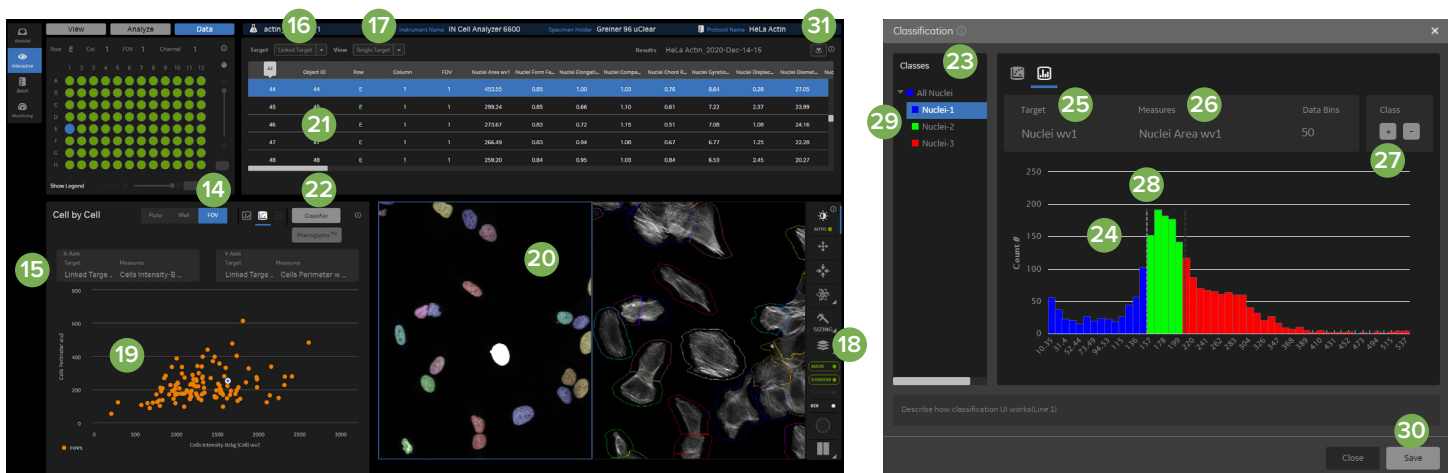
The **Classification** window is composed of the **Classes** navigation panel (23) and the **Histogram** (24). As classes

are added to the **Histogram**, the **Classes** navigation panel will display relationships between parent and daughter classes.

1. From the **Chart Dashboard**, click **Classifier** (22).
2. In the **Classification** window, select **Target** (25) and **Measures** (26) to use for classification.
3. Click the + (27) icon to add a gate to the histogram. Move the gate (28) horizontally to reposition. Up to five gates can be added to a single node.
4. In the **Classes navigation panel**, click on the **Pencil** icon to edit parent and daughter class names.
5. Double-click a node (29) in the **Classes navigation panel** to reload the histogram for that class alone. Use the + (27) icon to add daughter classes for this node.
6. When classification is complete, click **Save** (30) to classify data.

Exporting Measurements

1. Click on **Download Results**  (31) in the upper right hand corner of the data table panel.
2. In the **Download Data** window, choose the **Data Type** to export.
 - **All Data** will export cell by cell measurements AND summary data.
 - **Summary** will export only summary data (FOV and whole well averages).
3. Click **Download**. Results will be saved to the Downloads folder as a single zip file.



Contact Us

Phone: +1.800.635.5577
 Web: www.moleculardevices.com
 Email: info@moldev.com
 Check our website for a current listing of worldwide distributors.

Regional Offices

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

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