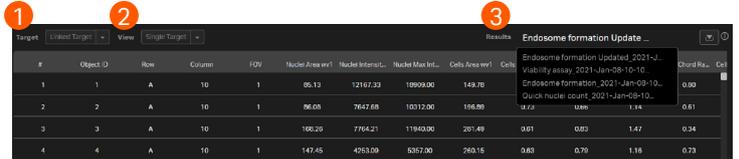


IN Carta software

View, classify and export results

Navigate 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.



#	Object ID	Row	Column	FOV	Nuclei Area ww1	Nuclei Intensity	Nuclei Max Int.	Cells Area ww1	Cells	Endosome formation Updated_2021-1-1	Chord Res.	Cell
1	1	A	10	1	85.13	12167.33	18099.00	149.70		Endosome formation Updated_2021-1-1	0.80	
2	2	A	10	1	86.05	7647.88	10312.00	196.90		Endosome formation_2021-Jan-09-10...	0.61	
3	3	A	10	1	100.29	7764.21	11940.00	201.49	0.61	0.03	1.42	0.34
4	4	A	10	1	147.45	4953.09	6387.00	249.15	0.63	0.79	1.16	0.73

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:

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



- Select the **Chart Type** ② in the Chart Dashboard.

			
Plate	✓		✓
Well	✓		
FOV	✓	✓	

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

- Select chart parameters:

- The **Histogram**  displays cell count for a single measure.
 - Use **Target** ① and **Measures** ② drop-downs to change the data displayed.
 - Adjust **Data Bins** ③ as needed. Valid entries are 1 to 99.
 - Click and drag over an area of the histogram ④ to zoom in on that portion of the graph ⑤. Click **Reset zoom** ⑥ to return to the default display.



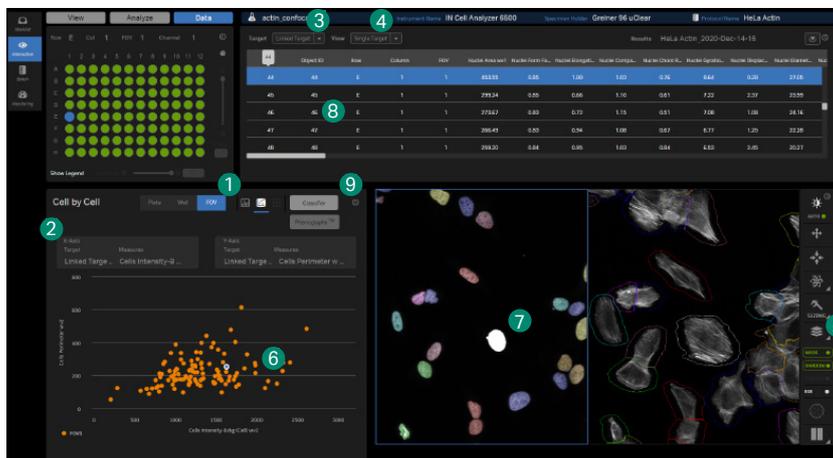
- The **Scatter Plot**  displays cell by cell data for a single FOV.
 - Use **Target** ① and **Measures** ② drop-downs to specify data to be plotted on the x- and y-axes.
 - Click and drag over an area of the scatter plot to zoom in on a particular area of the graph. Click the **Reset zoom** button to return to the default display.
- 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.



View linked data

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** ① and **Chart Type** to **Scatter Plot** .
2. Use drop-downs to select **Target** and **Measures** ② for the x- and y-axes.
3. In the **Data Table**, set **Target** to **Linked Target** ③ and **View** to **Single Target** ④.
4. Verify that the **View Port** is displaying the target identified in the **Scatter Plot** and that **Target Masks** are turned on ⑤.
5. Click on a point in the **Scatter Plot** ⑥, cell in the image ⑦, or row in the **Data Table** ⑧ to highlight corresponding data for that particular object.



Classify data

The **Classification** window is composed of the **Classes** navigation panel ① and the **Histogram** ②. As classes are added to the **Histogram**, the **Classes** navigation panel will display relationships between parent and daughter classes. ⑦

1. From the **Chart Dashboard**, click **Classifier** ⑨.
2. In the **Classification** window, select **Target** ③ and **Measures** ④ to use for classification.
3. Click the **+** ⑤ icon to add a gate to the histogram. Move the gate ⑥ 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 ⑦ in the **Classes navigation panel** to reload the histogram for that class alone. Use the **+** ⑤ icon to add daughter classes for this node.
6. When classification is complete, click **Save** ⑧ to classify data.



Export Measurements

1. Click on **Download Results**  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.

cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

All other third-party trademarks are the property of their respective owners.

Any use of this software is subject to the Cytiva Standard Software End-User License Agreement for Life Sciences Software Products. A copy of this Standard Software End-User License is available on request.

© 2020 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of these terms and conditions is available on request.

Contact your local Cytiva associate for the most current information.

IN Carta Software is for research use only – not for diagnostic use.

For local office contact information, visit [cytiva.com/contact](https://www.cytiva.com/contact)

