

## IN Carta software

# Protocol design (Mono-nucleated Cells application)

### Set up protocol

1. Select the **Analyze** tab.
2. In the **Applications** ① tab, select **Mono-nucleated Cells**.
3. In the **Protocols** ② tab, click the protocol you would like to run and select the **Modify** ③ tab. Alternatively, click **New** to design a new protocol.
4. Check the **Target Type** ⑤ boxes to activate the correct targets.
5. For each **Target Type**:
  - a. Enter a unique **Display Name** ⑥, if desired.
  - b. For organelles, specify the **Compartment** ⑦ in which they should be identified.
  - c. Select **Wavelength** ⑧ and a **Mask Display** ⑨ color.
  - d. Select **Segmentation** ⑩ method. Table at right provides details on the various methods.
  - e. Click the **Measures** ⑪ button and select/deselect measurements for each **Target Type** as needed.  
See "Using measurements menu" for more information.
6. For data sets with multiple z slices, it is possible to define a **Run mode** ⑫. It can be **Multiple Slices**, **Current Slice** or **Best Focus Z**. In order to specify a custom range of z slices, select **Multiple Slices** and click **Controls** button ⑬. Select **Best Focus Z** to allow IN Carta automatically determine the slice to be analyzed via image contrast analysis, for a selected channel. Click **Controls** button ⑬ to specify a channel used for select best focus z-slice.
7. Only one time-point can be analyzed in **Mono-nucleated Cells** application. The first point is selected by default, and another time-point selected for analysis in specimen navigation panel. Switch to **Time-lapse Analysis** application when analysis of multiple time-points is required.
8. Select the **Import Protocols** ④ tab to load previously created analysis protocols or protocols shared by other users.



Target type	Segmentation method	Description
Nuclei	Fast	<ul style="list-style-type: none"> <li>Faster than Robust method, but not as accurate</li> <li>Use for total intensity assays or when morphology information is not critical</li> </ul>
	Robust	<ul style="list-style-type: none"> <li>Slower than Fast method, but will provide more accurate nuclear segmentation results</li> </ul>
Cell	Fast	<ul style="list-style-type: none"> <li>Faster than Robust method, but not as accurate</li> <li>Use for well-separated, reasonably large cells</li> </ul>
	Collar	<ul style="list-style-type: none"> <li>Nuclear dilation method</li> <li>Use when there is no cellular staining, when identifying positive/negative staining, or when precise cell segmentation is not required</li> </ul>
	Robust	<ul style="list-style-type: none"> <li>Watershed algorithm that very accurately defines cell boundary</li> <li>Use when morphology or identification of objects within the cell boundary is important</li> </ul>
	Soma	<ul style="list-style-type: none"> <li>Use with Neurites segmentation</li> <li>Modified Robust Cell segmentation algorithm to segment soma</li> </ul>
Organelle	Fast puncta	<ul style="list-style-type: none"> <li>Multi-scale top hat algorithm</li> <li>Good at identifying puncta of varying size</li> </ul>
	Robust puncta	<ul style="list-style-type: none"> <li>Good at identifying puncta of varying intensity</li> </ul>
	Networks	<ul style="list-style-type: none"> <li>Segments network like organelles within the cell (e.g., endoplasmic reticulum, mitochondria, golgi)</li> </ul>
	Membrane	<ul style="list-style-type: none"> <li>Creates an artificial boundary within the cell.</li> <li>Use to segment a membrane compartment or to create a new compartment within the cell, for example mitochondria</li> </ul>
	Fibers	<ul style="list-style-type: none"> <li>Detects filamentous structures including individual and collapsed fibers</li> </ul>
	Neurites	<ul style="list-style-type: none"> <li>Segments neurite outgrowth and performs its tracing</li> </ul>

### Optimize segmentation parameters

1. Use the **Analysis Settings Panel** to edit available parameters for each **Target Type**.  
**Note:** See the table on the next page or click the **Information** ⓘ icon for more details on parameters.
2. Click **Apply** to see preliminary segmentation results overlaid on the images.
3. Check segmentation results for each target using any or all of the strategies below:
  - Use the **Mask**  tool to toggle mask display on/off and to access the **Target Mask Properties** menu.
  - Measure object length and area to refine size and area estimates using the **Sizing**  tool.
  - Apply an **Area** or **Intensity Filter** as needed.
  - Select the **Results Review** tab to display plots of the measured data to refine estimates for algorithm parameters.
4. Update **Analysis Settings** as necessary and click **Apply**. Repeat steps 2 and 3 until segmentation results are satisfactory.
5. Select another well and repeat steps 2 and 3 to verify that **Analysis Settings** are appropriate across the plate.
6. Click **Save** ⑭.

### Using measurements menu

- **Measures** are divided into **Count**, **Spatial**, **Intensity**, **Position**, **Coordination**, **Texture** and **Colocalization** categories.
- Each category has some default measurements selected and other available measure can be selected as needed.
- A channel selection is located in the **Intensity**, **Texture** and **Colocalization** categories. It allows to perform intensity-relevant measurements on wavelengths other than the one used for segmentation. For example, segment nuclei in the DAPI channel and measure intensity within a nuclear mask in FITC channel.

## Run Protocol

1. Click **Run Protocol** .
2. Click and drag or click **Select All** to select wells for analysis.  
**Note:** Press Ctrl during selection to select non-adjacent wells.
3. Click **Run**.  
**Note:** See Using batch analysis to run more than one protocol at a time.

## Using batch analysis

1. Click **Batch** .
2. Select **New Batch**.
3. Enter a name for the **Batch Container** and click **Ok**.
4. With the **Batch Container** selected in the **Batch Containers panel**, select multiple data sets from the **Acquisition Experiments panel** and select one or multiple analysis protocol. Click **Add Selected**.
5. Click **Submit for Processing**. Results are saved to each image stack location.
6. Click the **Monitoring** icon to see the status of the batch analysis.

Parameter	Targets			Explanation	Tips
	Nuclei	Cells	Organelles		
<b>Area filter</b>	✓	✓	✓	A post segmentation step to remove large or small objects from final data	<ul style="list-style-type: none"> <li>Use to remove objects with an area outside the range selected (i.e., small segmented background noise or large artifacts)</li> </ul>
<b>Background suppression</b>			✓	Advanced parameter for neurite segmentation to correct for low contrast	<ul style="list-style-type: none"> <li>Default value is a good starting point and should be changed only when changing sensitivity and width parameters don't improve segmentation</li> <li>Higher values reduce noise in the segmentation mask</li> </ul>
<b>Cell area</b>		✓		Average cell area ( $\mu\text{m}^2$ )	<ul style="list-style-type: none"> <li>Use size tools to establish appropriate value</li> <li><math>2 \times</math> the nuclear area is a good estimate for monolayer cells</li> <li>After initial segmentation use <b>Results Review</b> plot to refine</li> </ul>
<b>Collar radius</b>		✓		Distance ( $\mu\text{m}$ ) to be dilated from the edges of the nuclear mask	<ul style="list-style-type: none"> <li>Use to generate an artificial cell boundary when there is not a sufficient cell marker</li> </ul>
<b>Connected component segmentation</b>			✓	Advanced parameter for neurite segmentation, which triggers the use of an algorithm to segment broken neurites	<ul style="list-style-type: none"> <li>Turn on by default and turned off only when changing sensitivity and width parameters don't improve segmentation</li> </ul>
<b>Diameter</b>	✓			Typical target diameter ( $\mu\text{m}$ )	<ul style="list-style-type: none"> <li>Use size tools to establish appropriate value.</li> <li>After initial segmentation, use <b>Results Review</b> plot to refine</li> </ul>
<b>Enhance Image</b>			✓	Advanced parameter for neurite target to improve image contrast prior to segmentation	<ul style="list-style-type: none"> <li>Turned off by default and should be turned on only when changing sensitivity and width parameters don't improve segmentation</li> </ul>
<b>Intensity filter</b>	✓	✓	✓	A post segmentation step to remove bright or dim objects from final data	<ul style="list-style-type: none"> <li>Use to exclude objects with a mean intensity outside the selected range</li> </ul>
<b>Min. length</b>			✓	Approximate length ( $\mu\text{m}^2$ ) of the shortest fiber to be segmented	<ul style="list-style-type: none"> <li>Larger values will detect longer continuous objects while smaller values will detect small fragments or break up incorrectly joint objects</li> <li>This value is not a filter so the algorithm may still identify objects smaller than the specified length</li> </ul>
<b>Minimum target area</b>	✓			Approximate area ( $\mu\text{m}^2$ ) of smallest nucleus (Used for optimization of other algorithm functions. This does not mean that smaller nuclei cannot be segmented)	<ul style="list-style-type: none"> <li>Use size tools to establish appropriate value</li> <li>Default value is a good starting point</li> <li>After initial segmentation, use <b>Results Review</b> plot to refine</li> </ul>
<b>Minimum/maximum target area or target diameter</b>			✓	Approximate area ( $\mu\text{m}^2$ ) or diameter ( $\mu\text{m}$ ) of the smallest and largest organelles to be segmented	<ul style="list-style-type: none"> <li>These values adjust the steps of the algorithm to favor objects of a specified size</li> </ul>
<b>Minimum/Maximum width</b>			✓	Approximate thickness ( $\mu\text{m}$ ) of the smallest and largest fibers and neurites to be segmented	<ul style="list-style-type: none"> <li>To segment individual fibers, use very small min value since actin filament width is typically close to the resolution limit</li> <li>Increase Min/Max width for lower magnification images when segmenting individual fibers/neurites isn't possible</li> </ul>
<b>Noise removal</b>	✓	✓	✓	Noise removal filter to improve segmentation results for samples with few cells and/or low contrast	<ul style="list-style-type: none"> <li>Use if background is high, illumination is uneven, or if there is debris in the field-of-view (FOV)</li> </ul>
<b>Ratio</b>		✓		Controls the size of soma required for downstream neurite segmentation	<ul style="list-style-type: none"> <li>Similar values result in larger soma</li> <li>Decrease sensitivity first if soma is over-segmented. If soma mask still expands beyond cell borders then increase ratio value</li> </ul>
<b>Scales</b>			✓	Number of distinct sizes of objects to be identified	<ul style="list-style-type: none"> <li>Increasing the number of scales allows for detection of a larger range of organelle sizes</li> <li>When scales equal                             <ul style="list-style-type: none"> <li><b>1:</b> Only the minimum area is considered</li> <li><b>2:</b> Both minimum and maximum values are used</li> <li><b>3 or more:</b> Values between minimum and maximum are used</li> </ul> </li> </ul>
<b>Sensitivity</b>	✓	✓	✓	Determines whether pixels are assigned to objects or background based on intensity relative to local background	<ul style="list-style-type: none"> <li>Lower contrast images typically require higher sensitivity</li> <li>If object borders are not well defined, adjust sensitivity</li> <li>If too many small false positives are found in background, use the area filter to remove</li> </ul>
<b>Sensitivity threshold</b>			✓	Noise suppression filter	<ul style="list-style-type: none"> <li>Increase if too many organelles are identified with <b>Sensitivity</b> slider at lowest value</li> <li>Decrease if too few organelles are identified with <b>Sensitivity</b> slider at maximum value</li> <li>Increase or decrease by increments of 0.25</li> </ul>
<b>Thickness (membrane)</b>			✓	Distance ( $\mu\text{m}$ ) to erode cell boundary to create membrane compartment	<ul style="list-style-type: none"> <li>After membrane is defined, it can be used as a compartment for subsequent organelles</li> </ul>
<b>Thickness (network)</b>			✓	Thickness ( $\mu\text{m}$ ) of individual structures of the network	<ul style="list-style-type: none"> <li>Default value is a good starting point</li> <li>Increase to identify thicker structures, decrease to identify thinner structures</li> </ul>
<b>Tubeness</b>			✓	Advanced parameter for neurite target to obtain continuous segmentation mask	<ul style="list-style-type: none"> <li>Default value is a good starting point and should be changed only when changing sensitivity and width parameters don't improve segmentation</li> <li>Higher values result in smoother segmentation masks</li> </ul>