

IN Carta Image Analysis Software

QUICK REFERENCE GUIDE

Protocol Design (Mono-Nucleated Cells)

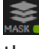
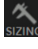
Setting Up a Protocol

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 on next page provides details on the various methods.
 - e. Click the **Measures (11)** button and select/deselect measurements for each **Target Type** as needed. See “Using the Measurements Menu” below 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 (13)**. Select **Best Focus Z** to allow the IN Carta™ software automatically determine the slice to be analyzed via image contrast analysis, for a selected channel. Click **Controls (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 time point is selected by default, and another time point selected for analysis in specimen navigation panel. Switch to the **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.

Optimizing 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 one or more of the following strategies:
 - 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.

| | 5 | 6 | 7 | 8 | 9 | 10 | 11 |
|--------------------|--|--------------|-------------|------------|--------------|-----------------|--------------|
| | Target Type | Display Name | Compartment | Wavelength | Mask Display | Segmentation | Measurements |
| 1 Applications | <input checked="" type="checkbox"/> Nuclei | Nuclei | | 5 | BLUE | Fast | Measures |
| 2 Protocols | <input checked="" type="checkbox"/> Cells | Cells | | 1 | GREEN | Robust | Measures |
| | <input checked="" type="checkbox"/> Organelles | Mitochondria | Cytoplasm | 2 | RED | Networks | Measures |
| 3 Modify | <input checked="" type="checkbox"/> Organelles 1 | Actin | Cells | 3 | FUCHSIA | Fibers | Measures |
| | <input checked="" type="checkbox"/> Organelles 2 | Centrosomes | Cells | 4 | ORANGE | Robust Puncta | Measures |
| 4 Import Protocols | 12 Run mode: Current Slice | 13 Controls | | | | 15 Run Protocol | 14 Save |

- Update **Analysis Settings** as necessary and click **Apply**. Repeat steps 2 and 3 until segmentation results are satisfactory.
- Select another well and repeat steps 2 and 3 to verify that **Analysis Settings** are appropriate across the plate.
- Click **Save (14)**.

Using the Measurements Menu


- Measures** are divided into **Count**, **Spatial**, **Intensity**, **Position**, **Coordination**, **Texture**, and **Colocalization** categories.
- Each category has some default measurements selected. Other available measures 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.

Running a Protocol

- Click **Run Protocol (15)**.
- Click and drag or click **Select All** to select wells for analysis.
Note: Press **Ctrl** during selection to select non-adjacent wells.
- Click **Run**.

Note: See Using Batch Analysis to run more than one protocol at a time.

Using Batch Analysis

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

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

| Parameter | Nuclei | Cells | Organelles | Description | Tips |
|--|--------|-------|------------|--|---|
| Area filter | ✓ | ✓ | ✓ | Post segmentation step to remove large or small objects from final data | <ul style="list-style-type: none"> Use to remove objects with an area outside the range selected (that is, small segmented background noise or large artifacts). |
| Background suppression | | | ✓ | Advanced parameter for neurite segmentation to correct for low contrast | <ul style="list-style-type: none"> Default value is a good starting point and should be changed only when changing sensitivity and width parameters don't improve segmentation. Higher values reduce noise in the segmentation mask. |
| Cell area | | ✓ | | Average cell area (μm^2) | <ul style="list-style-type: none"> Use size tools to establish appropriate value. $2 \times$ the nuclear area is a good estimate for monolayer cells. After initial segmentation use <i>Results Review</i> plot to refine. |
| Collar radius | | ✓ | | Distance (μm) to be dilated from the edges of the nuclear mask | <ul style="list-style-type: none"> Use to generate an artificial cell boundary when there is not a sufficient cell marker. |
| Connected component segmentation | | | ✓ | Advanced parameter for neurite segmentation, which triggers the use of an algorithm to segment broken neurites | <ul style="list-style-type: none"> Turn on by default and turned off only when changing sensitivity and width parameters don't improve segmentation. |
| Diameter | ✓ | | | Typical target diameter (μm) | <ul style="list-style-type: none"> Use size tools to establish appropriate value. After initial segmentation, use <i>Results Review</i> plot to refine. |
| Enhance image | | | ✓ | Advanced parameter for neurite target to improve image contrast prior to segmentation | <ul style="list-style-type: none"> Turned off by default and should be turned on only when changing sensitivity and width parameters don't improve segmentation. |
| Intensity filter | ✓ | ✓ | ✓ | A post segmentation step to remove bright or dim objects from final data | <ul style="list-style-type: none"> Use to exclude objects with a mean intensity outside the selected range. |
| Min. length | | | ✓ | Approximate length (μm) of the shortest fiber to be segmented | <ul style="list-style-type: none"> Larger values will detect longer continuous objects; smaller values will detect small fragments or break up incorrectly joint objects. This value is not a filter so the algorithm may still identify objects smaller than the specified length. |
| Minimum target area | ✓ | | | Approximate area (μm^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"> Use size tools to establish appropriate value. Default value is a good starting point. After initial segmentation, use <i>Results Review</i> plot to refine. |
| Minimum/maximum target area or target diameter | | | ✓ | Approximate area (μm^2) or diameter (μm) of the smallest and largest organelles to be segmented | <ul style="list-style-type: none"> These values adjust the steps of the algorithm to favor objects of a specified size. |
| Minimum/maximum width | | | ✓ | Approximate thickness (μm) of the smallest and largest fibers and neurites to be segmented | <ul style="list-style-type: none"> To segment individual fibers, use very small min value since actin filament width is typically close to the resolution limit. Increase Min/Max width for lower magnification images when segmenting individual fibers/neurites isn't possible. |
| Noise removal | ✓ | ✓ | ✓ | Noise removal filter to improve segmentation results for samples with few cells and/or low contrast | <ul style="list-style-type: none"> Use if background is high, illumination is uneven, or if there is debris in the field-of-view (FOV). |

| Parameter | Nuclei | Cells | Organelles | Description | Tips |
|-----------------------|--------|-------|------------|---|---|
| Ratio | | ✓ | | Controls the size of soma required for downstream neurite segmentation | <ul style="list-style-type: none"> • Similar values result in larger soma. • Decrease sensitivity first if soma is over-segmented. If soma mask still expands beyond cell borders then increase ratio value. |
| Scales | | | ✓ | Number of distinct sizes of objects to be identified | <ul style="list-style-type: none"> • Increasing the number of scales allows for detection of a larger range of organelle sizes. • When scales equal: <ol style="list-style-type: none"> 1: Only the minimum area is considered 2: Both minimum and maximum values are used 3 or more: Values between minimum and maximum are used |
| Sensitivity | ✓ | ✓ | ✓ | Determines whether pixels are assigned to objects or background based on intensity relative to local background | <ul style="list-style-type: none"> • Lower contrast images typically require higher sensitivity. • If object borders are not well defined, adjust sensitivity. • If too many small false positives are found in background, use the area filter to remove. |
| Sensitivity threshold | | | ✓ | Noise suppression filter | <ul style="list-style-type: none"> • Increase if too many organelles are identified with Sensitivity slider at lowest value. • Decrease if too few organelles are identified with Sensitivity slider at maximum value. • Increase or decrease by increments of 0.25. |
| Thickness (membrane) | | | ✓ | Distance (µm) to erode cell boundary to create membrane compartment | <ul style="list-style-type: none"> • After membrane is defined, it can be used as a compartment for subsequent organelles. |
| Thickness (network) | | | ✓ | Thickness (µm) of individual structures of the network | <ul style="list-style-type: none"> • Default value is a good starting point. • Increase to identify thicker structures, decrease to identify thinner structures. |
| Tubeness | | | ✓ | Advanced parameter for neurite target to obtain continuous segmentation mask | <ul style="list-style-type: none"> • Default value is a good starting point and should be changed only when changing sensitivity and width parameters do not improve segmentation. • Higher values result in smoother segmentation masks. |

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 | 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