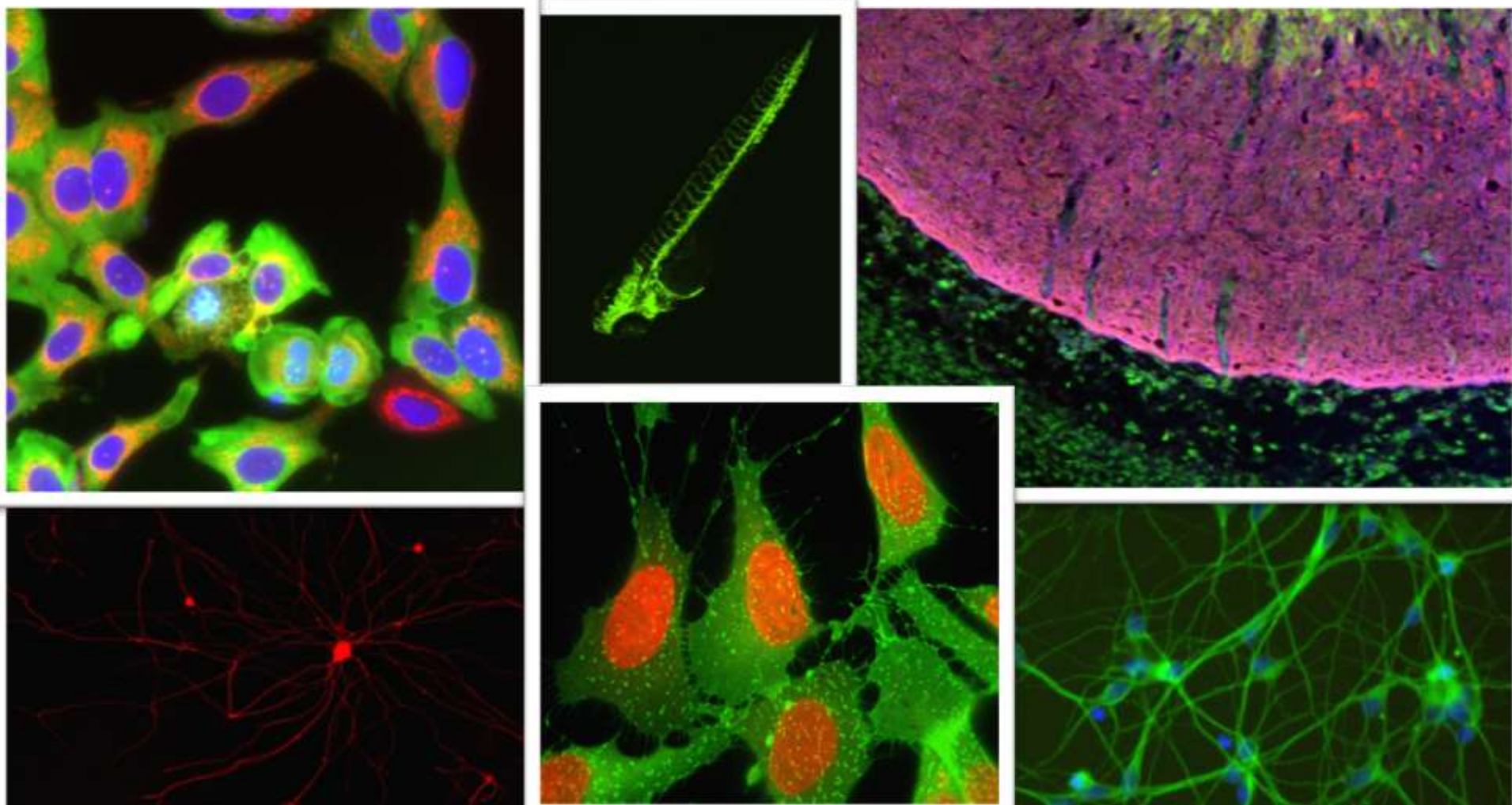




**Molecular
Devices**

Together through life sciences.

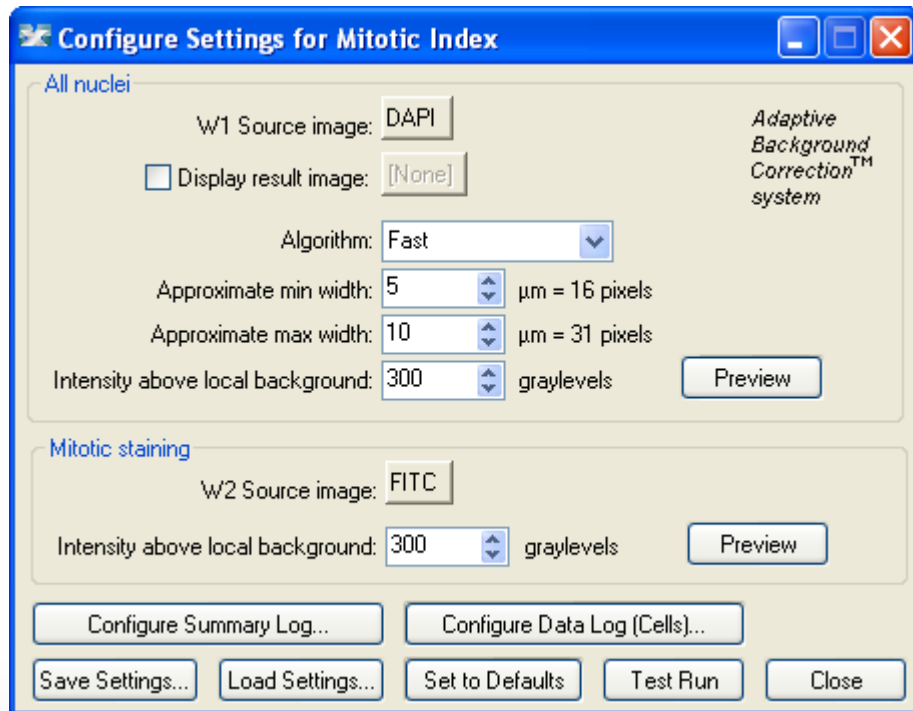
MetaXpress® Software: *Mitotic Index Module*

Together through life sciences.



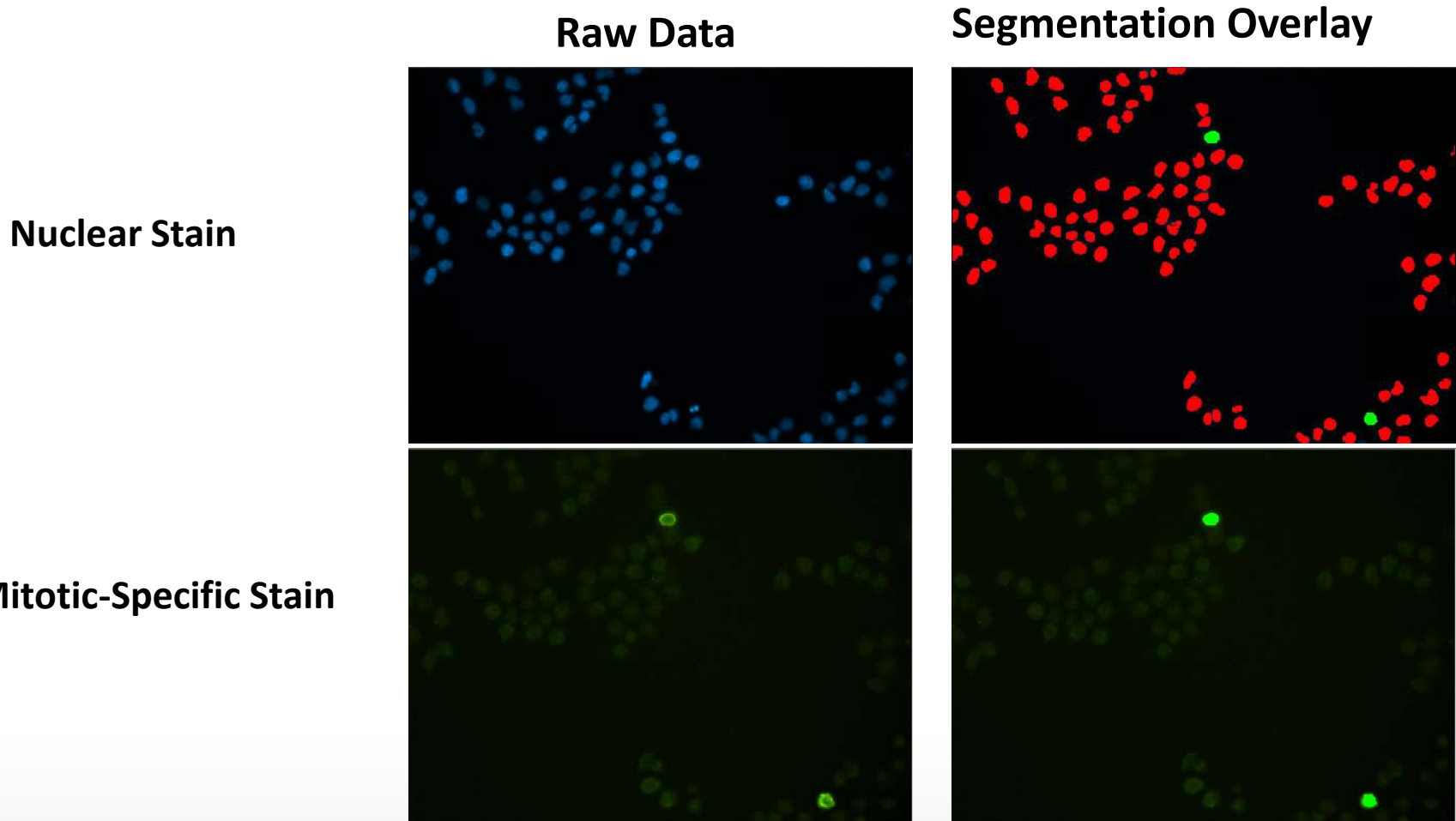
**Molecular
Devices**

Mitotic Index Module Overview



- The Mitotic Index module can be used to analyze cellular images to differentiate between Mitotic and Interphase cells in a normal cell cycle.
- This module requires a nuclear wavelength and mitotic-specific stain wavelength.
- A typical mitotic-specific stain used for this module is Histone 3 S10 phosphorylation. DNA stain labels all cells and the mitotic cells are labeled only with the mitotic-specific stain.

Mitotic Index – Example images



Interphase nuclei are shown with a red overlay and mitotic nuclei with a green overlay.

Module Settings – Nuclei classification

- Nuclei classification

Configure Settings for Mitotic Index

All nuclei

W1 Source image:

☐ Display result image:

Algorithm:

Approximate min width: $\mu\text{m} = 16$ pixels

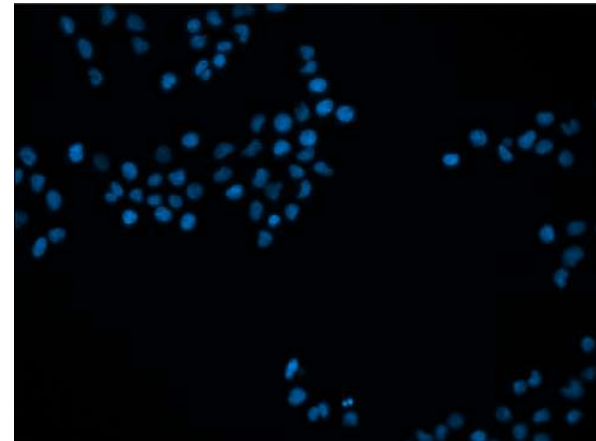
Approximate max width: $\mu\text{m} = 31$ pixels

Intensity above local background: graylevels

Mitotic staining

W2 Source image:

Intensity above local background: graylevels



2. Module Settings – DNA content

Configure Settings for Mitotic Index

All nuclei

W1 Source image: **DAPI**

☐ Display result image: [None]

Algorithm: Fast

Approximate min width: 5 μm = 16 pixels

Approximate max width: 10 μm = 31 pixels

Intensity above local background: 300 graylevels

Mitotic staining

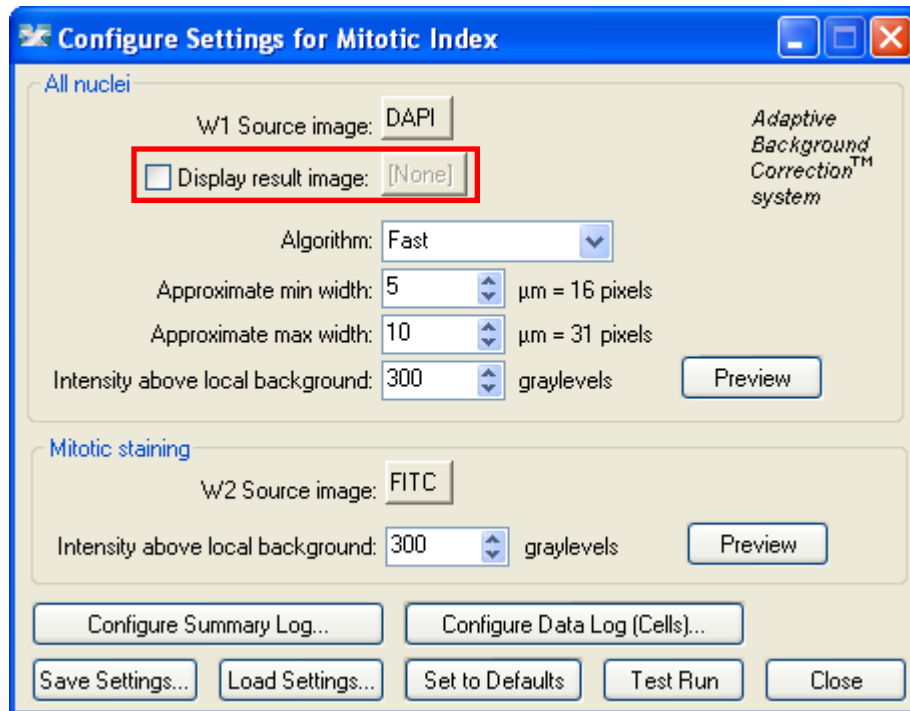
W2 Source image: FITC

Intensity above local background: 300 graylevels

Buttons: Configure Summary Log..., Configure Data Log (Cells)..., Save Settings..., Load Settings..., Set to Defaults, Test Run, Close

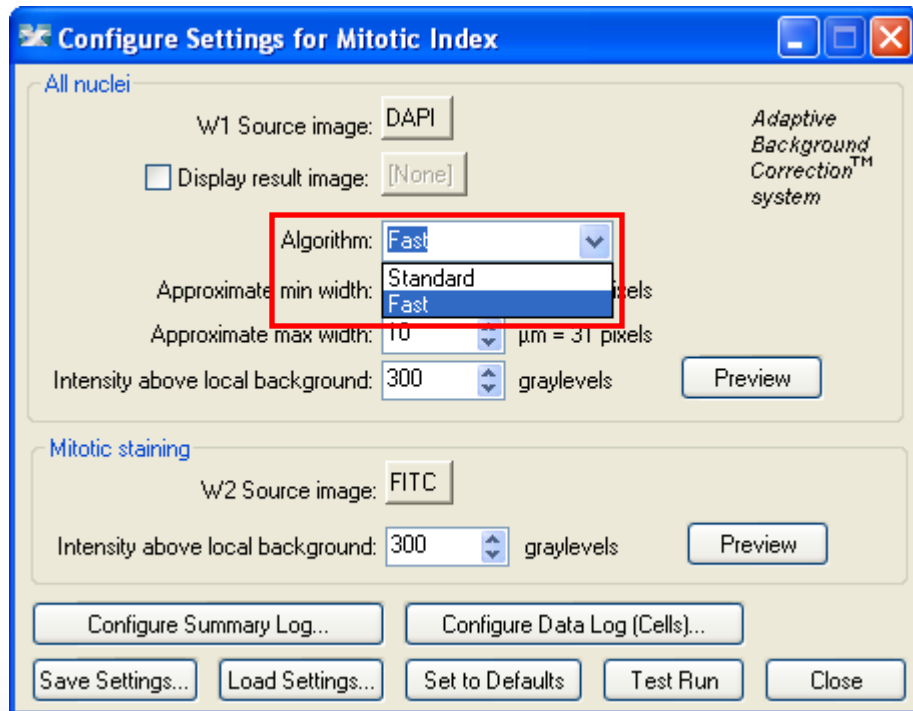
- Select the wavelength for the **DNA content** (nuclear stain)

1. Module Settings – result image



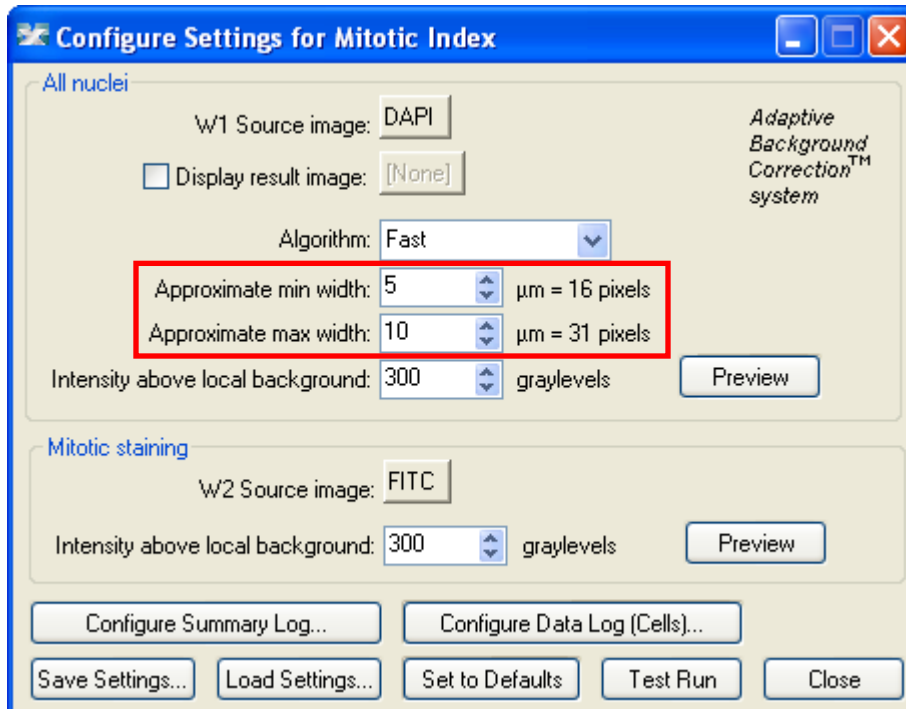
- Leave “**Display result image**” deselected (this is generally only used when journaling)

Module Settings

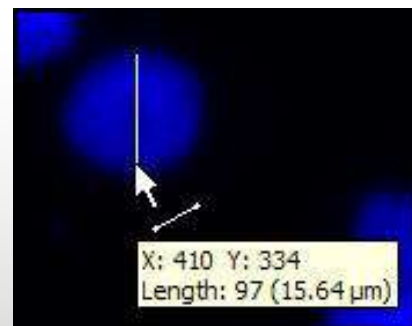


- **Algorithm**
- This option is only available in MetaXpress software version 4.0 and higher and determines how quickly the analysis is performed.
- **Fast** algorithm can perform analysis up to twice as fast as **Standard**.
- Both algorithms produce similar but not identical results.

3. Module Settings – width settings

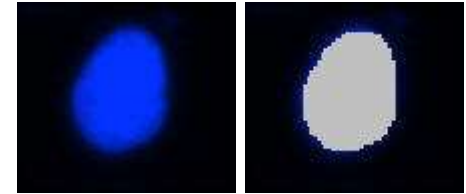


- Set the **Approximate min width** and **Approximate max width** for the range of nuclei that you want to detect
- The width is the short axis of a nucleus (in μm).
- The region tools can be used to measure widths
- Much smaller cells will be ignored
- Much larger cells will be split



3. Module Settings – width settings

Effects of varying width settings



Min width too small: splits nuclei



Min width too large: omits smaller nuclei

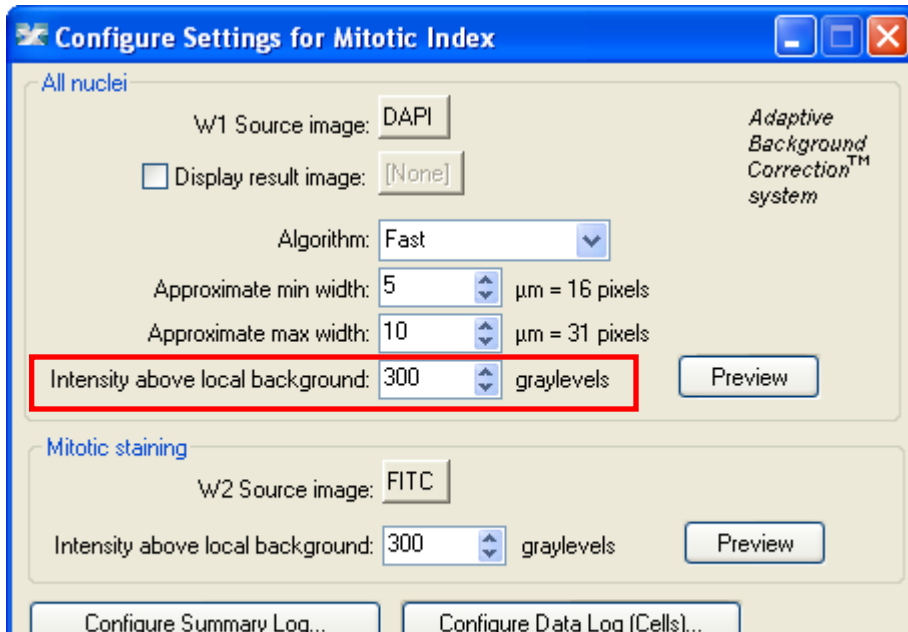


Max width too small: may shrink nuclear boundaries

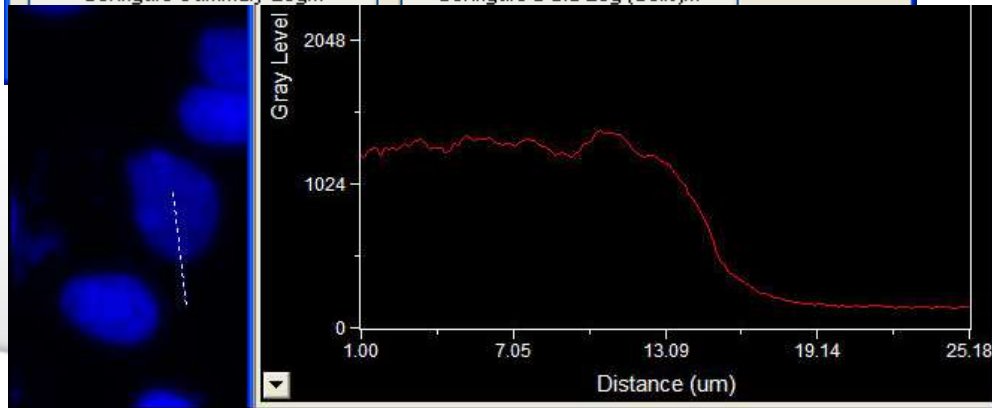


Max width too large: may slightly enlarge nuclear boundaries

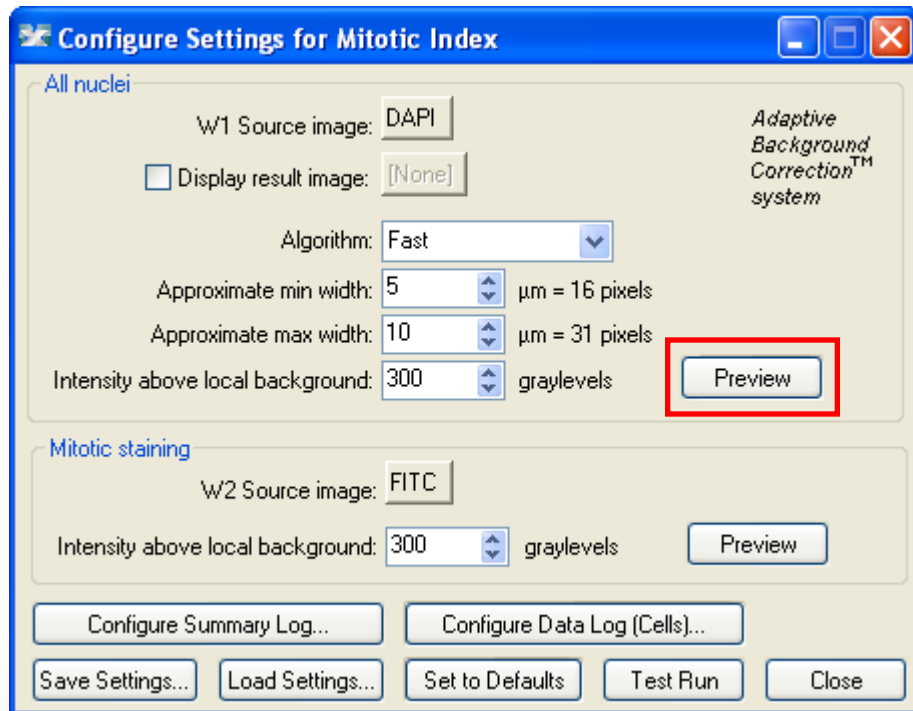
Module Settings – Intensity above background



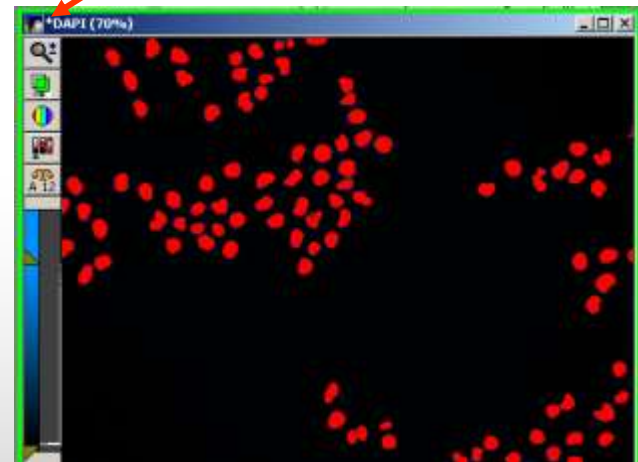
- The **intensity above local background** is used for finding the nuclei
- This value is a minimum and should be set slightly lower than the difference in intensity between a dim cell and its local background. For FAST algorithm, set this value to about half (or less) of the difference in intensity between a dim cell and local background.
- Draw a line across a cell into the background and use Measure > Linescan to determine intensity values; or simply mouse over the cell and the background and view the intensity values



4. Module Settings – Nuclei segmentation

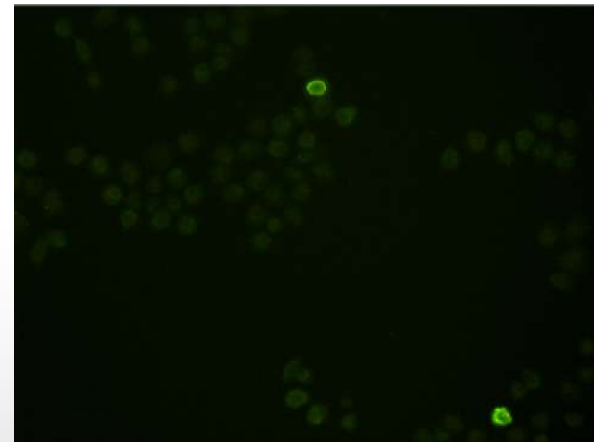
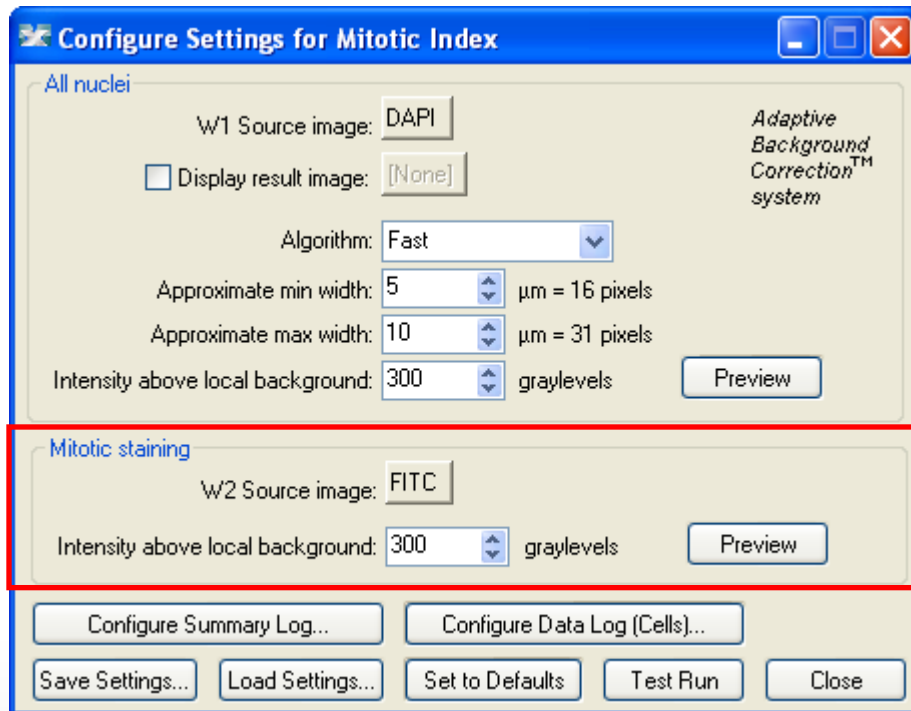


- Click on **Preview** to test the segmentation settings for the nuclei
- The Preview image will show a red overlay over the segmented nuclei. The overlay can be toggled on and off.



5. Module Settings – Mitotic Cell classification

- Mitotic classification
- The average intensity (brightness) of the mitotic-specific marker (e.g. phospho-histone H3) is used to identify mitotic cells.



6. Module Settings – mitotic classification

Configure Settings for Mitotic Index

All nuclei

W1 Source image: DAPI

☐ Display result image: [None]

Algorithm: Fast

Approximate min width: 5 $\mu\text{m} = 16$ pixels

Approximate max width: 10 $\mu\text{m} = 31$ pixels

Intensity above local background: 300 graylevels

Mitotic staining

W2 Source image: FITC

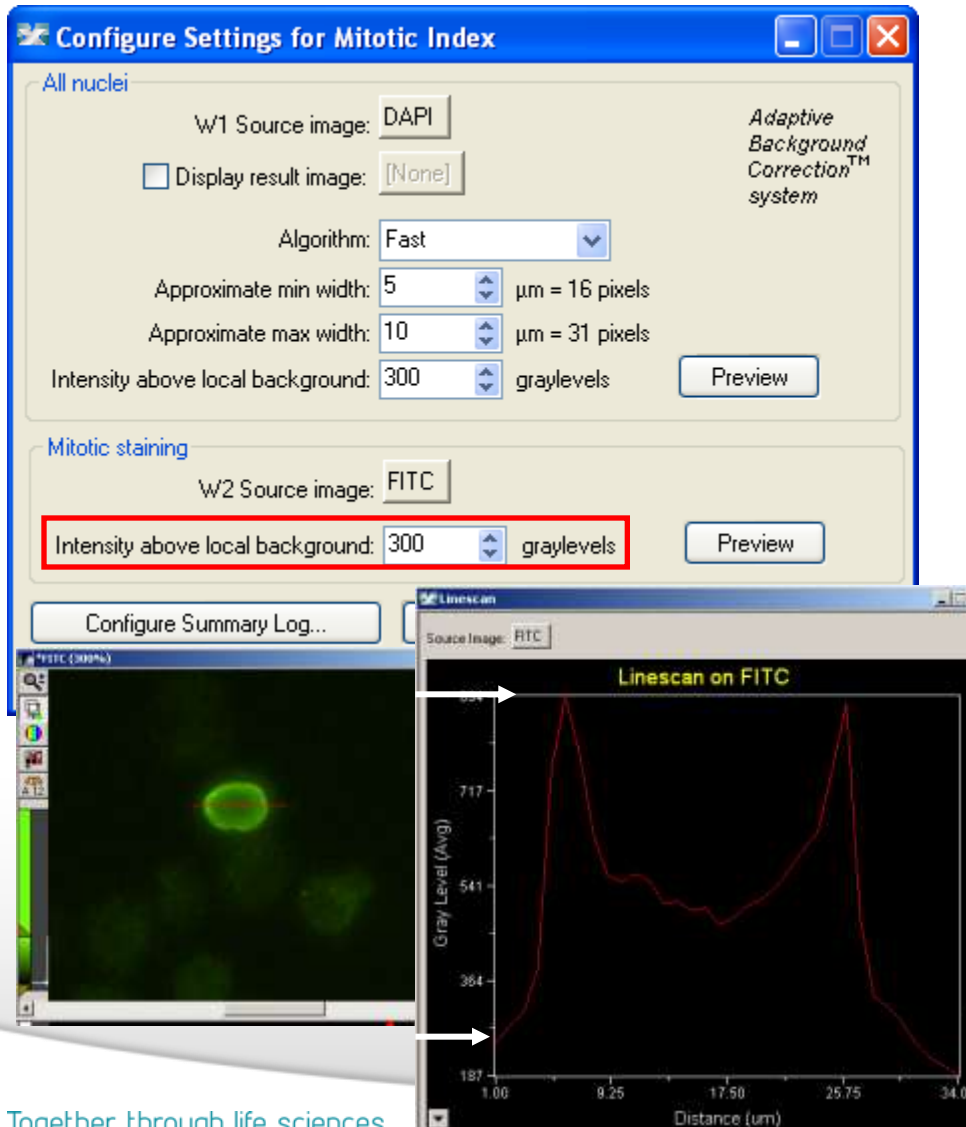
Intensity above local background: 300 graylevels

Buttons: Configure Summary Log..., Configure Data Log (Cells)...
Save Settings..., Load Settings..., Set to Defaults, Test Run, Close

- **Mitotic-specific stain**
- Select the wavelength for the mitotic marker

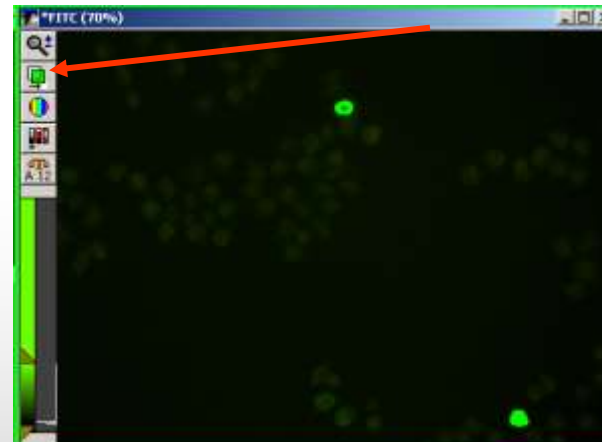
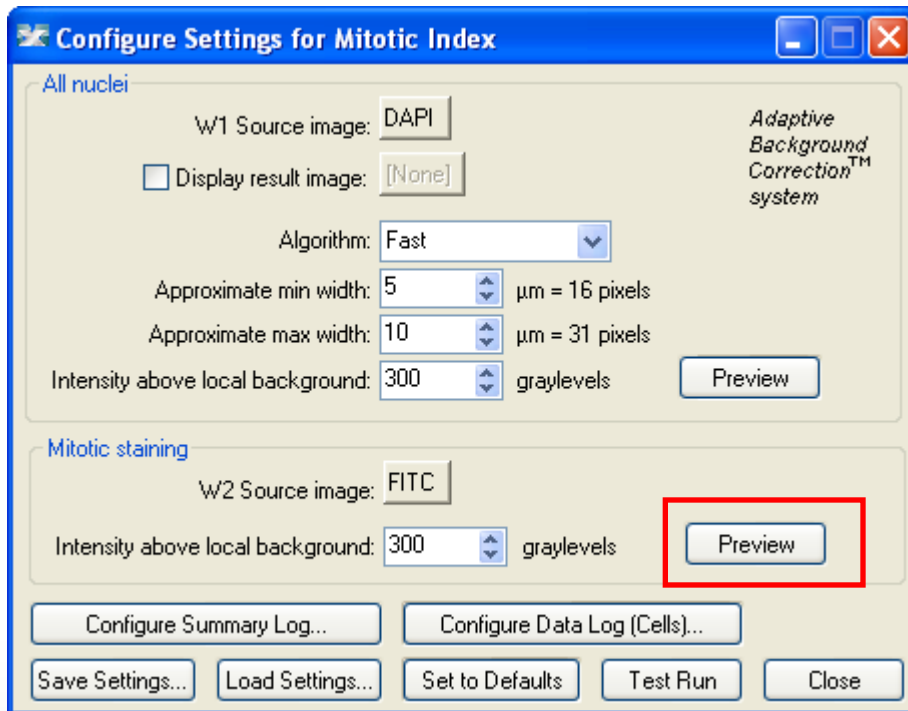
7. Module Settings – mitotic classification

- Enter a cutoff intensity value for the mitotic specific stain.
- This value is a minimum and should be set slightly lower than the difference in intensity between a dim mitotic cell and its local background. For FAST algorithm, set this value to about half (or less) of the difference in intensity between a dim cell and local background.
- Draw a line across a cell into the background and use Measure > Linescan to determine intensity values; or simply mouse over the cell and the background and view the intensity values.



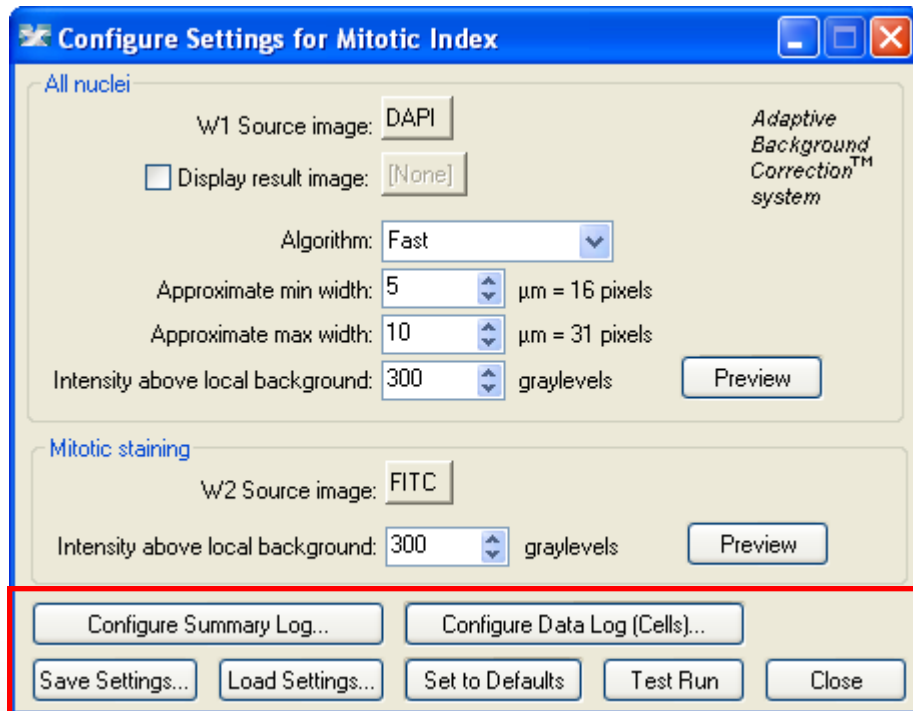
8. Module Settings – mitotic classification

- Press **Preview** to see cells with that level of staining highlighted in the image.
- The Preview image will show a green overlay on the mitotic staining image to show the segmentation derived from the image. The overlay can be toggled on and off.



9. Module Settings – final settings

- **Configure Summary Log** – select site-by-site measurements
- **Configure Data Log** – select cell-by-cell measurements
- **Save Settings** – save analysis parameters to database
- **Load Settings** – load saved analysis parameters
- **Set to Defaults** – restore default analysis parameters
- **Test Run** – test all settings together and display cell-by-cell results for this site



Summary Data (site-by-site measurements)

- ✓ Image Name
- ✓ Image Plane
- ✓ Image Date and Time
- ✓ Elapsed Time
- ✓ Stage Label
- ✓ Wavelength
- ✓ Z Position
- ✓ Total Nuclei
- ✓ Mitotic Nuclei
- ✓ Interphase Nuclei
- ✓ % Mitotic Nuclei
- ✓ % Interphase Nuclei
- ✓ All Nuclei Total Area
- ✓ All Nuclei Mean Area
- ✓ All Nuclei W1 Integrated Intensity
- ✓ All Nuclei W1 Average Intensity
- ✓ All Nuclei W2 Integrated Intensity
- ✓ All Nuclei W2 Average Intensity
- ✓ Mitotic Total Area
- ✓ Mitotic Mean Area
- ✓ Mitotic W1 Integrated Intensity
- ✓ Mitotic W1 Average Intensity
- ✓ Mitotic W2 Integrated Intensity
- ✓ Mitotic W2 Average Intensity
- ✓ Interphase Total Area
- ✓ Interphase Mean Area
- ✓ Interphase W1 Integrated Intensity
- ✓ Interphase W1 Average Intensity
- ✓ Interphase W2 Integrated Intensity
- ✓ Interphase W2 Average Intensity

- **Summary or site-by-site measurements are for the entire image and include:**

- **Total Nuclei:**

Total number cells determined from wavelength 1.

- **Mitotic Nuclei:**

Total number nuclei determined from wavelength 1 that were positive in wavelength 2.

- **Interphase Nuclei:**

Total number nuclei determined from wavelength 1 that were negative in wavelength 2.

- **% Mitotic Nuclei:**

$100 * \text{Mitotic Nuclei} / \text{Total Nuclei}$.

- **% Interphase Nuclei:**

$100 * \text{Interphase Nuclei} / \text{Total Nuclei}$.

Summary Data (site-by-site measurements)

- **Summary or site-by-site measurements are for the entire image and include:**

- ✓ Image Name
- ✓ Image Plane
- ✓ Image Date and Time
- ✓ Elapsed Time
- ✓ Stage Label
- ✓ Wavelength
- ✓ Z Position
- ✓ Total Nuclei
- ✓ Mitotic Nuclei
- ✓ Interphase Nuclei
- ✓ % Mitotic Nuclei
- ✓ % Interphase Nuclei
- ✓ All Nuclei Total Area
- ✓ All Nuclei Mean Area
- ✓ All Nuclei W1 Integrated Intensity
- ✓ All Nuclei W1 Average Intensity
- ✓ All Nuclei W2 Integrated Intensity
- ✓ All Nuclei W2 Average Intensity
- ✓ Mitotic Total Area
- ✓ Mitotic Mean Area
- ✓ Mitotic W1 Integrated Intensity
- ✓ Mitotic W1 Average Intensity
- ✓ Mitotic W2 Integrated Intensity
- ✓ Mitotic W2 Average Intensity
- ✓ Interphase Total Area
- ✓ Interphase Mean Area
- ✓ Interphase W1 Integrated Intensity
- ✓ Interphase W1 Average Intensity
- ✓ Interphase W2 Integrated Intensity
- ✓ Interphase W2 Average Intensity

- **All Nuclei Total Area:**

Total μm^2 s in wavelength 1 in all nuclei.

- **All Nuclei Mean Area:**

All Nuclei Total Area/Total Nuclei.

- **All Nuclei W1 Integrated Intensity:**

Summed grayscale values in wavelength 1 in all nuclei.

- **All Nuclei W1 Average Intensity:**

All Nuclei Integrated Intensity/All Nuclei Total (Pixel) Area.

- **All Nuclei W2 Integrated Intensity:**

Summed grayscale values in wavelength 2 in all nuclei.

- **All Nuclei W2 Average Intensity:**

All Nuclei W2 Integrated Intensity/All Nuclei Total (Pixel) Area.

Summary Data (site-by-site measurements)

- ✓ Image Name
- ✓ Image Plane
- ✓ Image Date and Time
- ✓ Elapsed Time
- ✓ Stage Label
- ✓ Wavelength
- ✓ Z Position
- ✓ Total Nuclei
- ✓ Mitotic Nuclei
- ✓ Interphase Nuclei
- ✓ % Mitotic Nuclei
- ✓ % Interphase Nuclei
- ✓ All Nuclei Total Area
- ✓ All Nuclei Mean Area
- ✓ All Nuclei W1 Integrated Intensity
- ✓ All Nuclei W1 Average Intensity
- ✓ All Nuclei W2 Integrated Intensity
- ✓ All Nuclei W2 Average Intensity
- ✓ Mitotic Total Area
- ✓ Mitotic Mean Area
- ✓ Mitotic W1 Integrated Intensity
- ✓ Mitotic W1 Average Intensity
- ✓ Mitotic W2 Integrated Intensity
- ✓ Mitotic W2 Average Intensity
- ✓ Interphase Total Area
- ✓ Interphase Mean Area
- ✓ Interphase W1 Integrated Intensity
- ✓ Interphase W1 Average Intensity
- ✓ Interphase W2 Integrated Intensity
- ✓ Interphase W2 Average Intensity

- **Mitotic Total Area:**

Total μm^2 s in wavelength 1 in mitotic nuclei.

- **Mitotic Mean Area:**

Mitotic Total Area/Mitotic Nuclei.

- **Mitotic W1 Integrated Intensity:**

Summed grayscale values in wavelength 1 in mitotic nuclei.

- **Mitotic W1 Average Intensity:**

Mitotic Nuclei W1 Integrated Intensity/Mitotic Nuclei Total (Pixel) Area.

- **Mitotic W2 Integrated Intensity:**

Summed grayscale values in wavelength 2 in mitotic nuclei.

- **Mitotic W2 Average Intensity:**

Mitotic W2 Integrated Intensity/Mitotic Total (Pixel) Area.

Summary Data (site-by-site measurements)

- ✓ Image Name
- ✓ Image Plane
- ✓ Image Date and Time
- ✓ Elapsed Time
- ✓ Stage Label
- ✓ Wavelength
- ✓ Z Position
- ✓ Total Nuclei
- ✓ Mitotic Nuclei
- ✓ Interphase Nuclei
- ✓ % Mitotic Nuclei
- ✓ % Interphase Nuclei
- ✓ All Nuclei Total Area
- ✓ All Nuclei Mean Area
- ✓ All Nuclei W1 Integrated Intensity
- ✓ All Nuclei W1 Average Intensity
- ✓ All Nuclei W2 Integrated Intensity
- ✓ All Nuclei W2 Average Intensity
- ✓ Mitotic Total Area
- ✓ Mitotic Mean Area
- ✓ Mitotic W1 Integrated Intensity
- ✓ Mitotic W1 Average Intensity
- ✓ Mitotic W2 Integrated Intensity
- ✓ Mitotic W2 Average Intensity
- ✓ Interphase Total Area
- ✓ Interphase Mean Area
- ✓ Interphase W1 Integrated Intensity
- ✓ Interphase W1 Average Intensity
- ✓ Interphase W2 Integrated Intensity
- ✓ Interphase W2 Average Intensity

- **Interphase Total Area:**

Total μm^2 s in wavelength 1 in interphase nuclei.

- **Interphase Mean Area:**

Interphase Total Area/Interphase Nuclei.

- **Interphase W1 Integrated Intensity:**

Summed grayscale values in wavelength 1 in interphase nuclei.

- **Interphase W1 Average Intensity:**

Interphase W1 Integrated Intensity/Interphase Total (Pixel) Area.

- **Interphase W2 Integrated Intensity:**

Summed grayscale values in wavelength 2 in interphase nuclei.

- **Interphase W2 Average Intensity:**

Interphase W2 Integrated Intensity/Interphase Total (Pixel) Area.

Cell Data (cell-by-cell measurements)

- ✓ Image Name
- ✓ Image Plane
- ✓ Image Date and Time
- ✓ Elapsed Time
- ✓ Stage Label
- ✓ Wavelength
- ✓ Z Position
- ✓ Cell: Assigned Label #
- ✓ Cell: Mitotic Classification
- ✓ Cell: Total Area
- ✓ Cell: W1 Integrated Intensity
- ✓ Cell: W1 Average Intensity
- ✓ Cell: W2 Integrated Intensity
- ✓ Cell: W2 Average Intensity

- **Data Log (Cell-by-Cell Measurement)** - For each nucleus, the following measurements are included in the Data Log:

- **Cell: Assigned Label #:**

Cell label number (1 through total cell number)

- **Cell: Classification:**

Interphase, Mitotic

- **Cell: Total Area:**

Total μm^2 s in the nucleus.

Cell Data (cell-by-cell measurements)

- ✓ Image Name
- ✓ Image Plane
- ✓ Image Date and Time
- ✓ Elapsed Time
- ✓ Stage Label
- ✓ Wavelength
- ✓ Z Position
- ✓ Cell: Assigned Label #
- ✓ Cell: Mitotic Classification
- ✓ Cell: Total Area
- ✓ Cell: W1 Integrated Intensity
- ✓ Cell: W1 Average Intensity
- ✓ Cell: W2 Integrated Intensity
- ✓ Cell: W2 Average Intensity

- **Data Log (Cell-by-Cell Measurement):** For each nucleus, the following measurements are included in the Data Log:
- **Cell: W1 Integrated Intensity:**
Summed grayscale values in wavelength 1 for this nucleus.
- **Cell: W1 Average Intensity:**
 $\text{W1 Integrated Intensity} / \text{Total (Pixel) Area}$
- **Cell: W2 Integrated Intensity:**
Summed grayscale values in wavelength 2 for this nucleus.
- **Cell: W2 Average Intensity:**
 $\text{W2 Integrated Intensity} / \text{Total (Pixel) Area}$



Together through life sciences.

www.moleculardevices.com