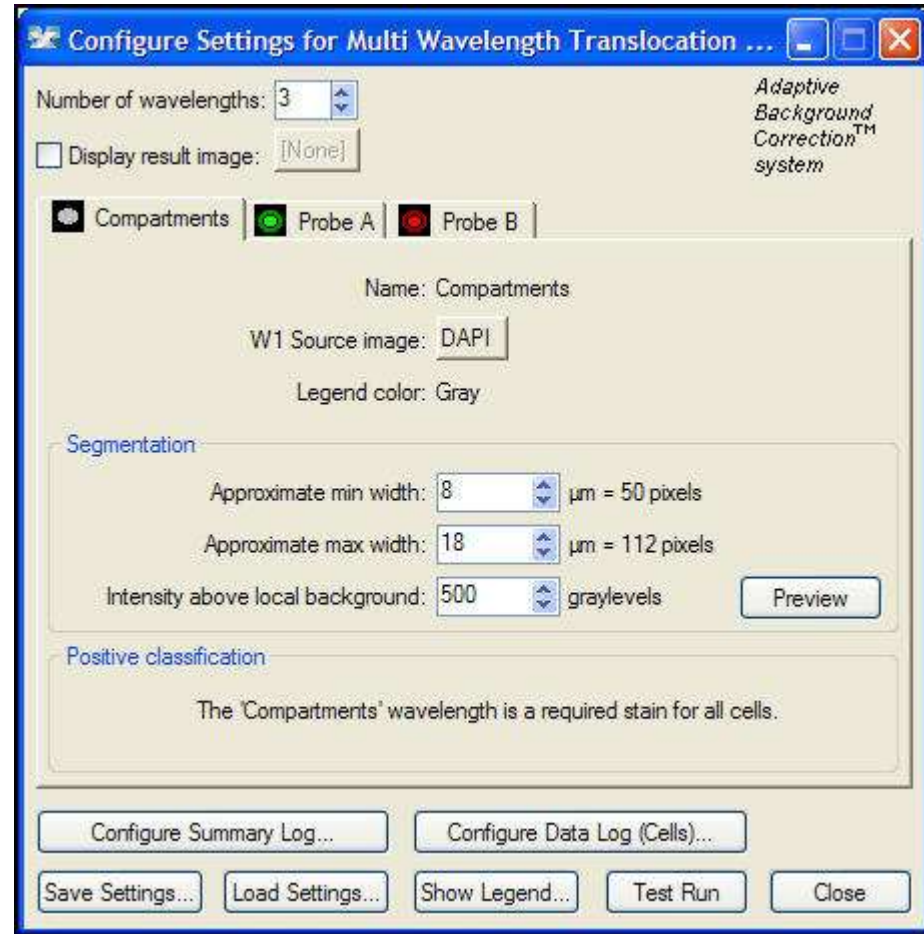


## **MetaXpress: Multi Wavelength Translocation module**

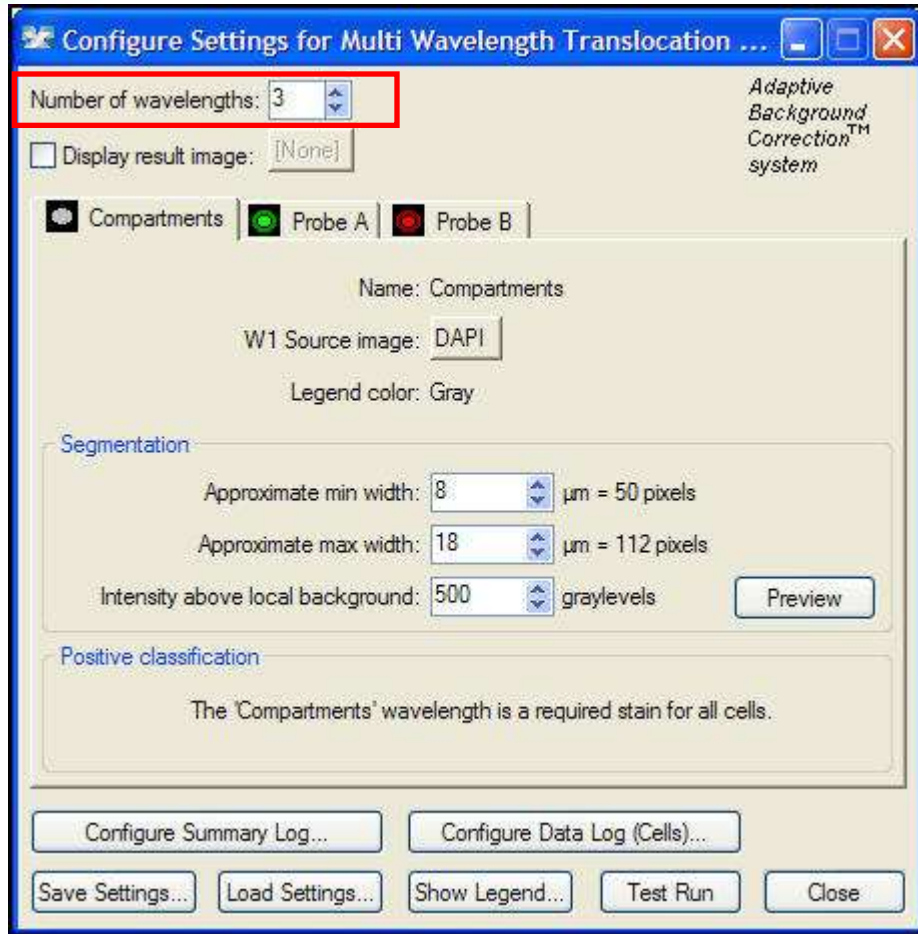
Paula Gedraitis, Ph.D.

# Multi Wavelength Translocation Module Overview

- Measures nuclear translocation in assays with a nuclear stain and 1-6 probes of interest
- Scores cells as positive or negative for each wavelength based on correlation with the nuclear stain
- Assigns profiles to cells based on their score in each wavelength

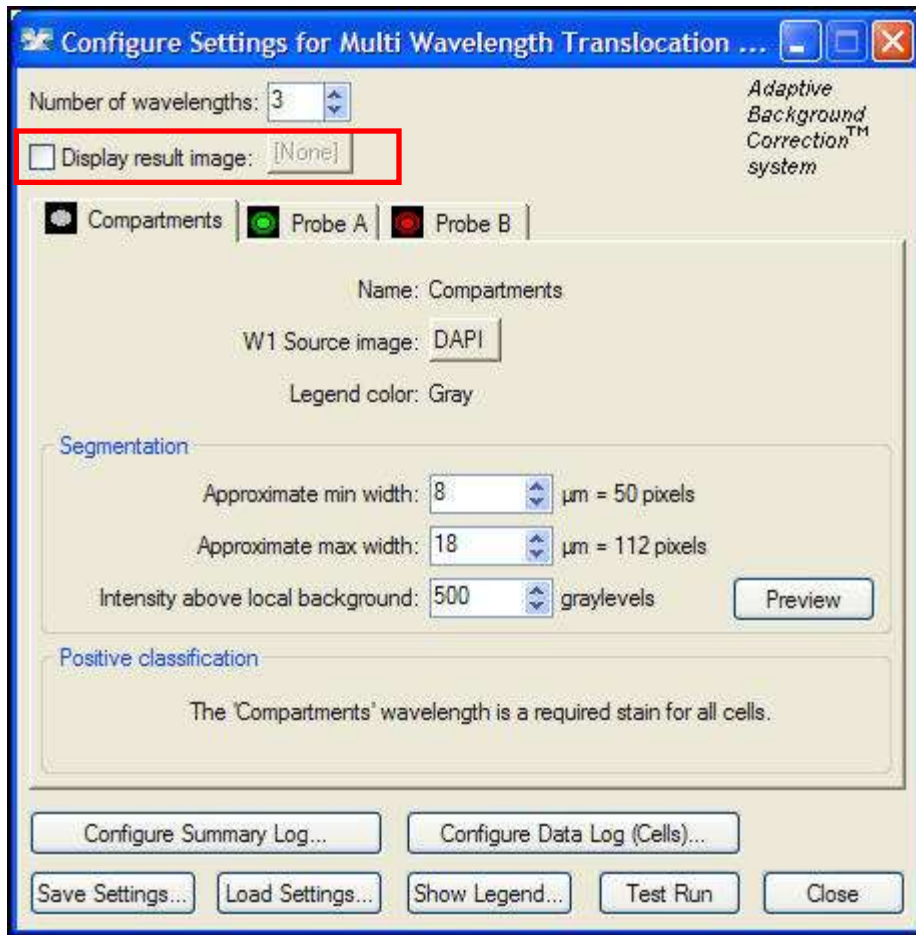


# Module Settings



- **Number of wavelengths**
- This should be set to the total number of wavelengths to be analyzed (nuclei plus probes of interest).
- Up to 7 different wavelengths can be analyzed simultaneously (wavelengths may also be repeated with different settings)

# Module Settings



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



# Module Settings: Wavelength 1

Configure Settings for Multi Wavelength Translocation ...

Number of wavelengths: 3

☐ Display result image: [None]

Adaptive Background Correction™ system

Compartments | Probe A | Probe B

Name: Compartments

W1 Source image: DAPI

Legend color: Gray

Segmentation

Approximate min width: 8  $\mu\text{m}$  = 50 pixels

Approximate max width: 18  $\mu\text{m}$  = 112 pixels

Intensity above local background: 500 graylevels

Preview

Positive classification

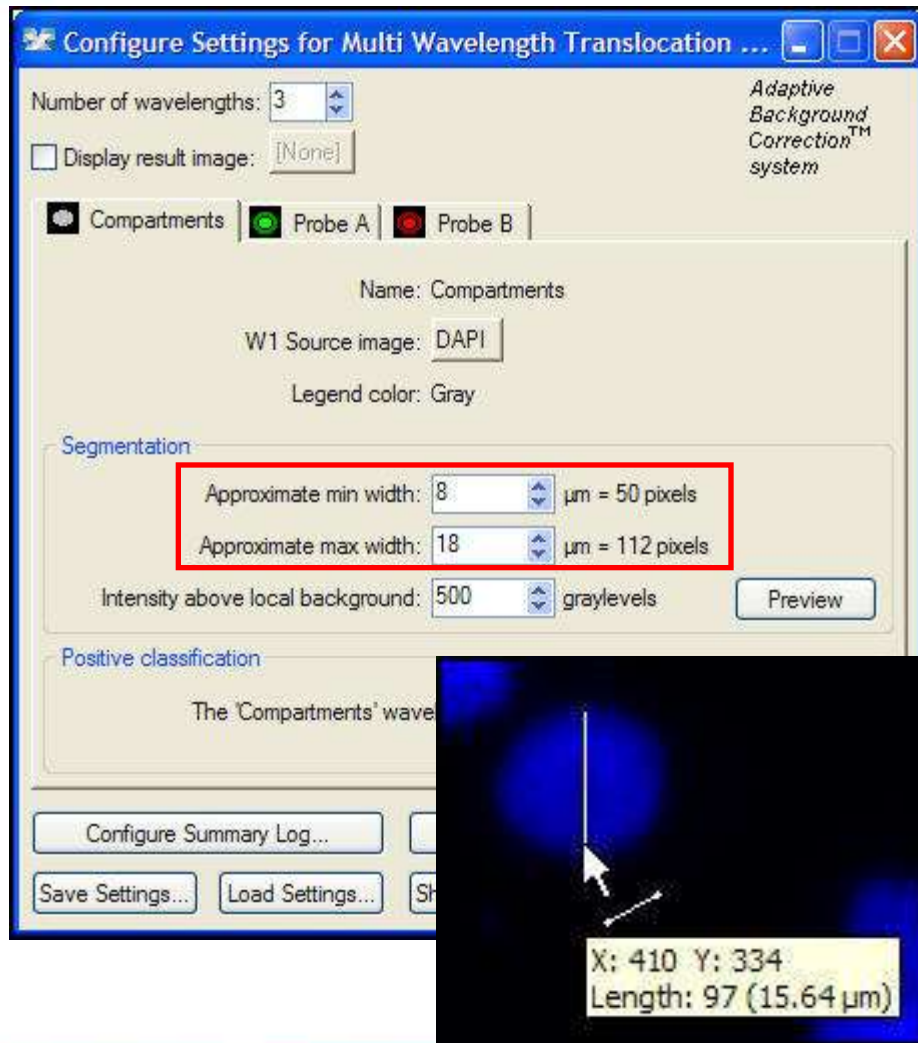
The 'Compartments' wavelength is a required stain for all cells.

Configure Summary Log... Configure Data Log (Cells)...

Save Settings... Load Settings... Show Legend... Test Run Close

- Select the wavelength for the **Compartments** (nuclei)

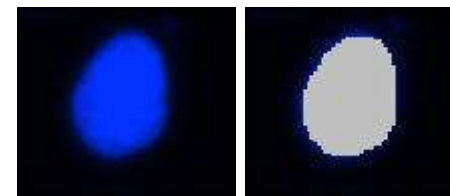
# Module Settings: Wavelength 1



- 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  $\mu\text{m}$ )
- The region tools can be used to measure widths
- Much smaller cells will be ignored
- Much larger cells will be split

# Module Settings: Wavelength 1

## Effects of varying width settings



Min width too small: splits nuclei



Min width too large: omits smaller nuclei

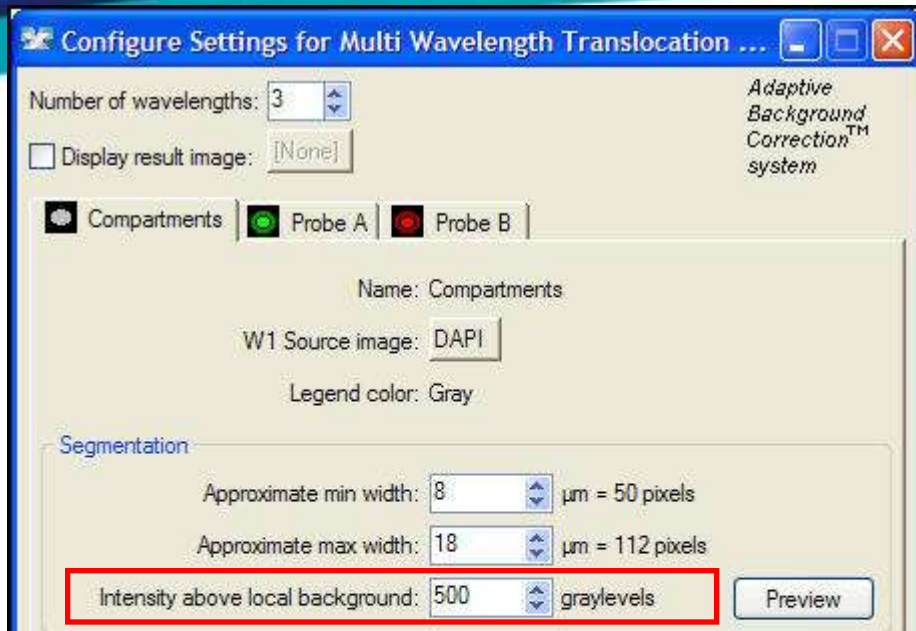
Max width too small: may shrink nuclear boundaries



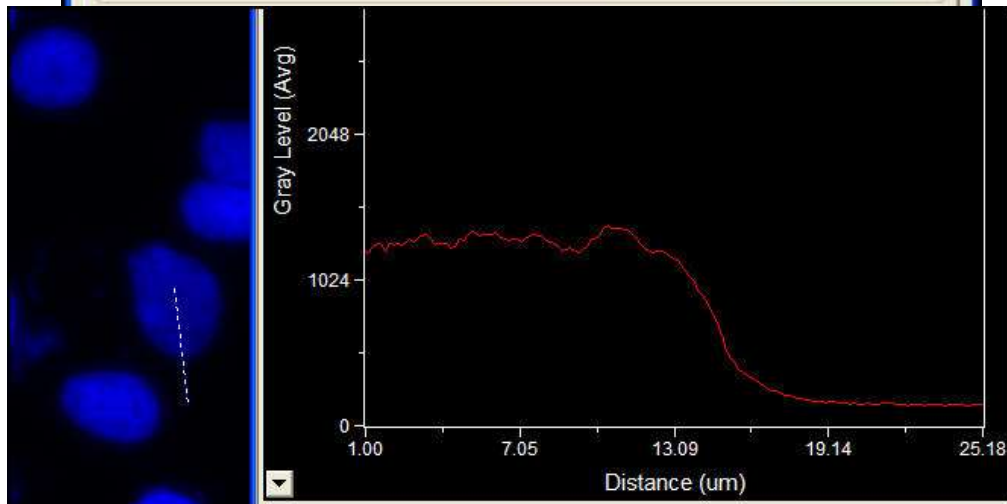
Max width too large: may slightly enlarge nuclear boundaries



# Module Settings: Wavelength 1

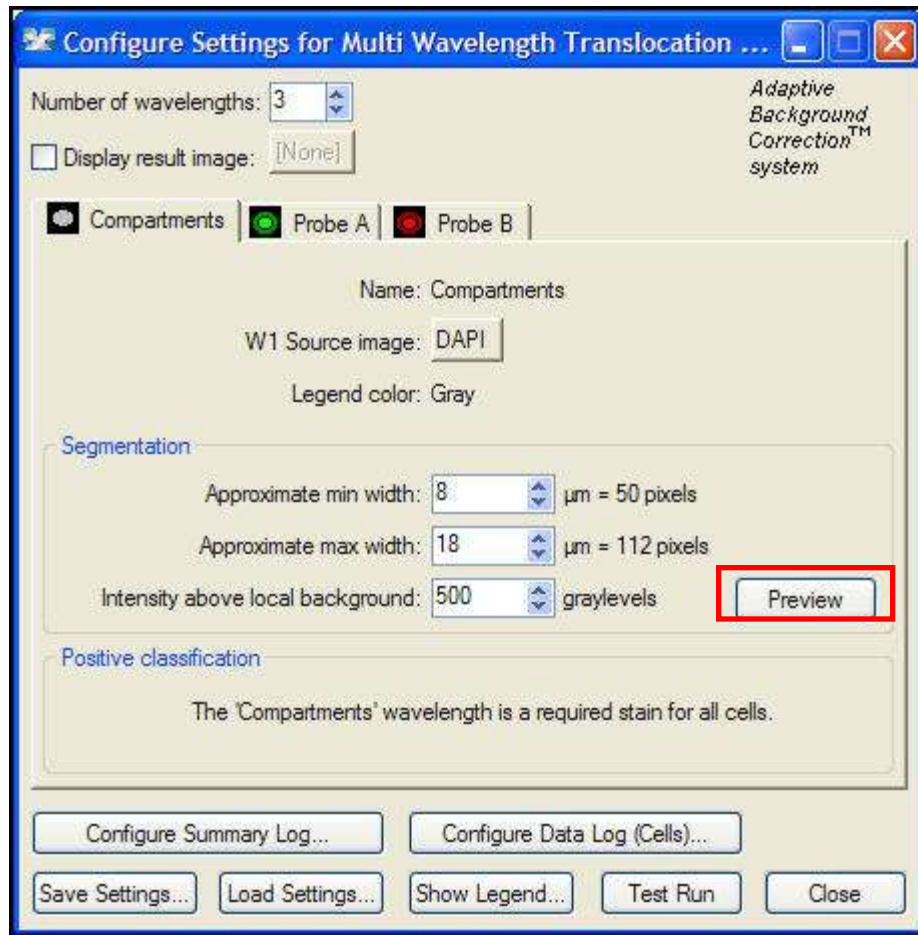


- The **intensity above local background** is used for finding the nuclei
- This value is a minimum and should be set slightly lower than the difference in intensity between a dim cell and its local background
- Draw a line across a cell into the background and use Measure > Linescan to determine intensity values; or simply mouse over the cell and the background and view the intensity values





# Module Settings: Wavelength 1



- Click on **Preview** to test settings for the current wavelength only

# Module Settings: Wavelength 2

Configure Settings for Multi Wavelength Translocation ...

Number of wavelengths: 3

☐ Display result image: [None]

Adaptive Background Correction™ system

☒ Compartments ☒ Probe A ☐ Probe B

Name: Probe A

W2 Source image: FITC

Legend color: Green

Define regions for measurement

Inner region distance in from edge: 1  $\mu\text{m}$  = 6 pixels

Outer region distance out from edge: 1  $\mu\text{m}$  = 6 pixels

Outer region width: 3  $\mu\text{m}$  = 19 pixels

Positive classification

Minimum correlation coefficient: 0.45

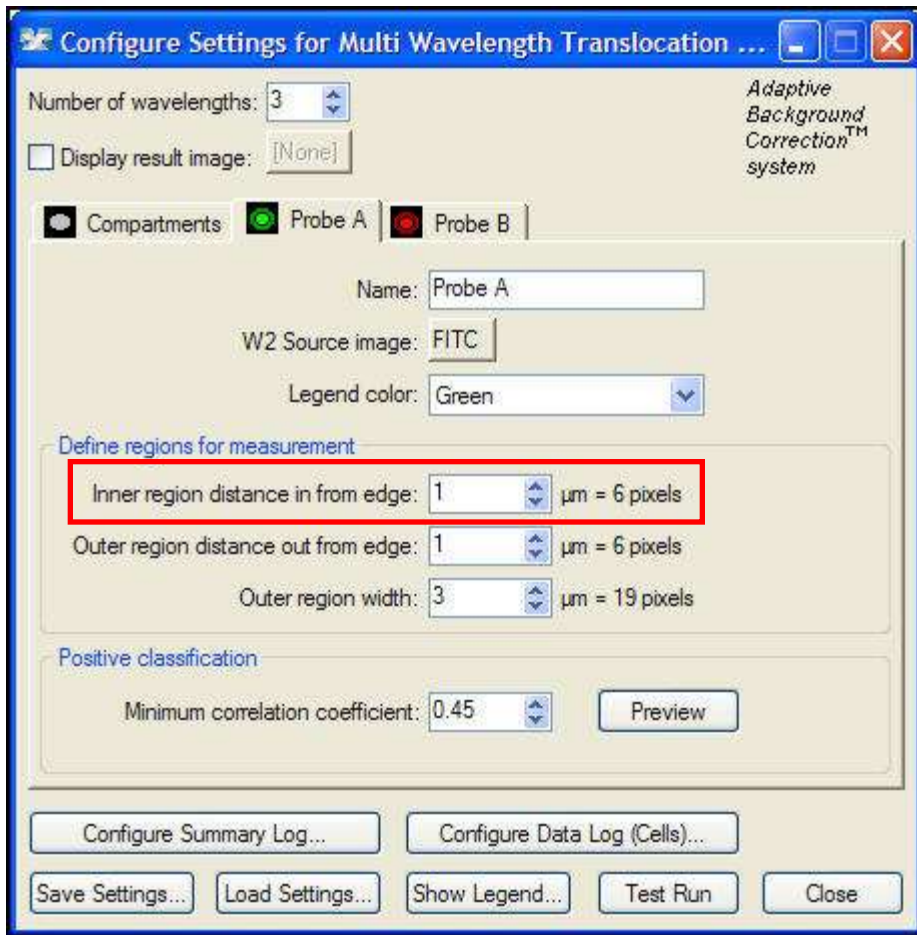
Preview

Configure Summary Log... Configure Data Log (Cells)...

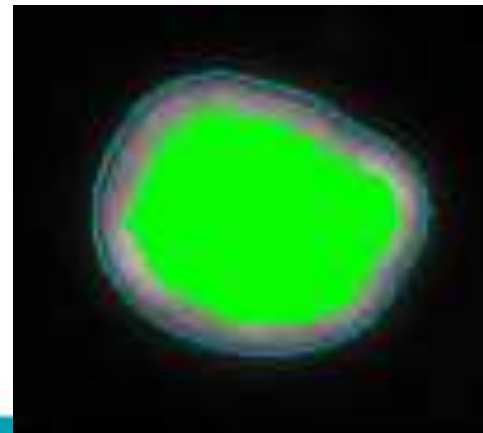
Save Settings... Load Settings... Show Legend... Test Run Close

- **Define W2:**
- Enter a name and select wavelength and legend color for your 2nd wavelength (probe A in this example)

# Module Settings: Wavelength 2



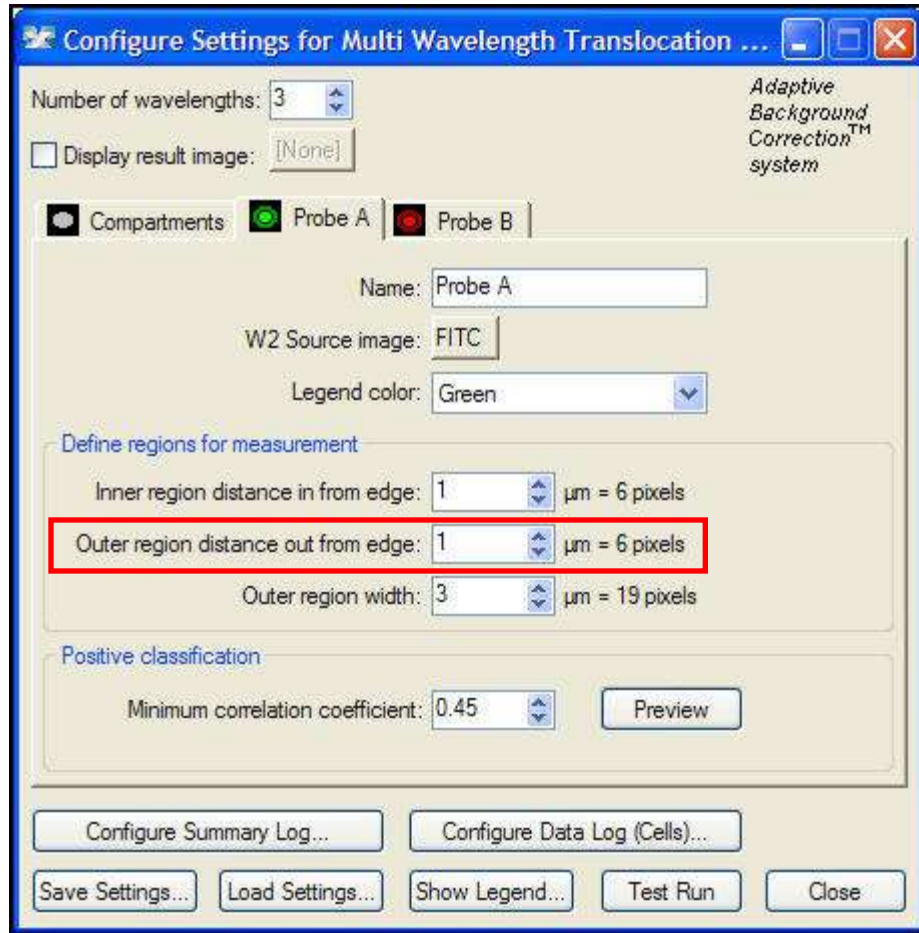
- **Define regions for measurement:**
- The "inner region" can be set to be shrunk in inner region from the detected nucleus to avoid boundary effects.
- 1  $\mu\text{m}$  is the default setting used in the Translocation module.



**2  $\mu\text{m}$  in**



# Module Settings: Wavelength 2



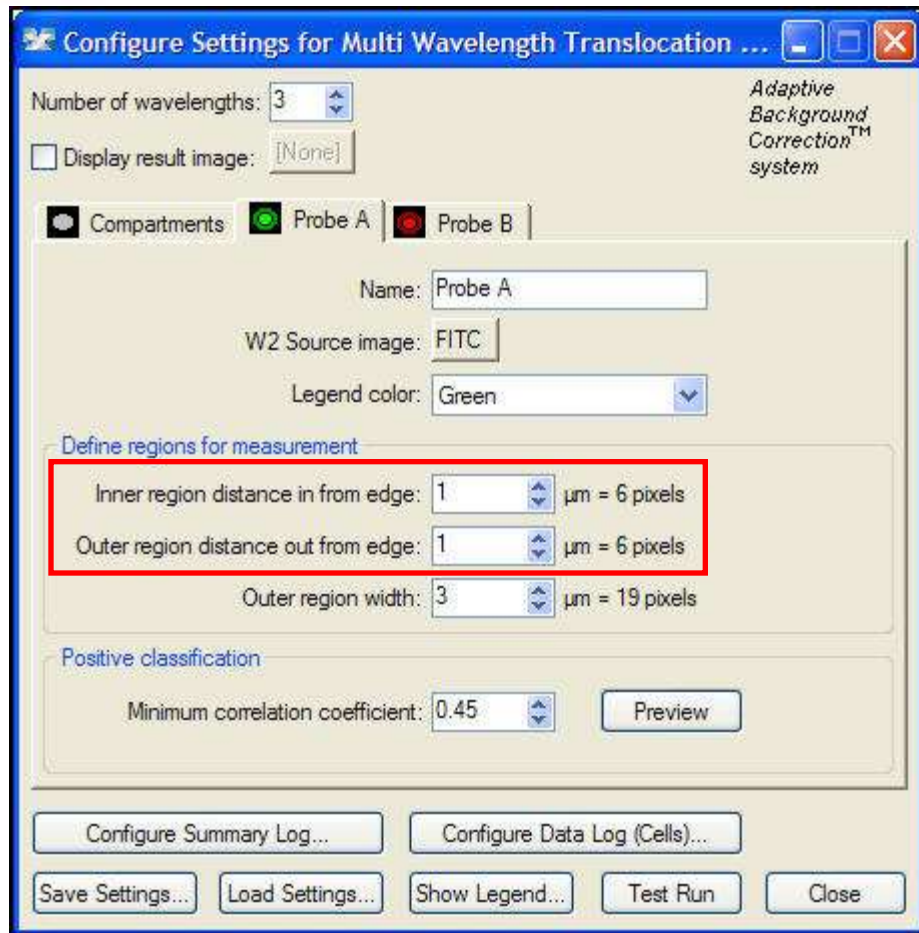
- **Define regions for measurement:**
- The “outer region” can be set to be expanded out from the detected nucleus to avoid boundary effects.
- 1  $\mu\text{m}$  is the default setting used in the Translocation module.



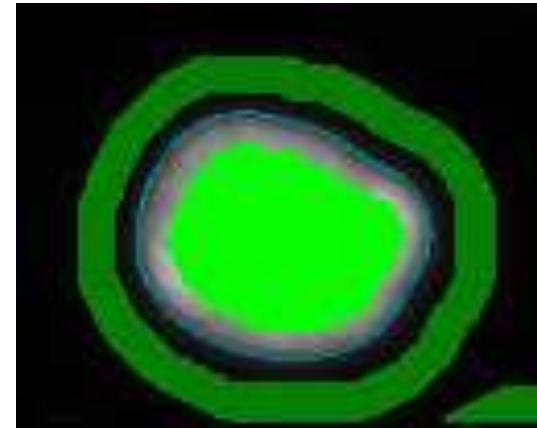
**2  $\mu\text{m}$  out**



# Module Settings: Wavelength 2



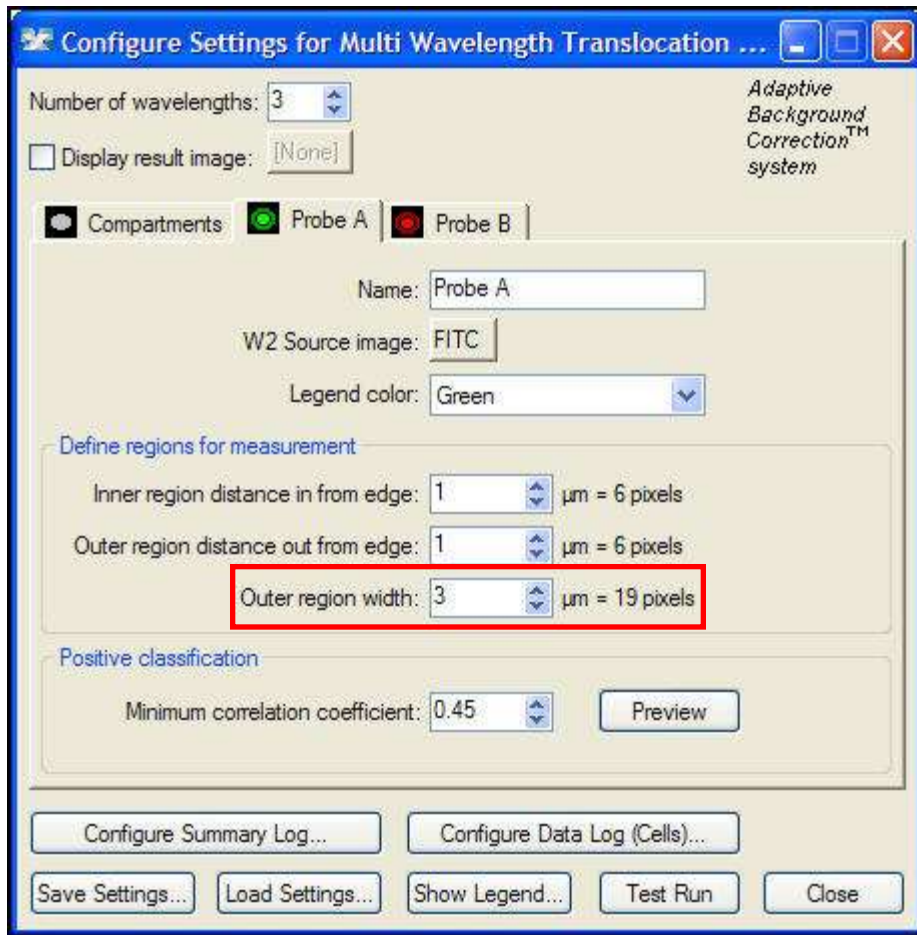
- Define regions for measurement:
- Typically the inner region and outer region distances are both utilized.



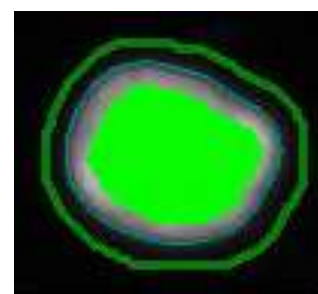
*2  $\mu\text{m}$  in and out*



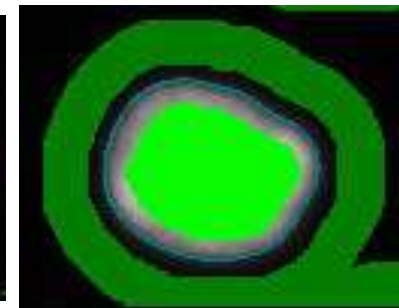
# Module Settings: Wavelength 2



- **Define regions for measurement:**
- The outer region width can be adjusted for the specific cell type (and degree of confluency).
- 3 μm is the default setting used in the Translocation module.



1 μm



5 μm

# Module Settings: Wavelength 2

Configure Settings for Multi Wavelength Translocation ...

Number of wavelengths: 3

☐ Display result image: [None]

Adaptive Background Correction™ system

☒ Compartments ☒ Probe A ☐ Probe B

Name: Probe A

W2 Source image: FITC

Legend color: Green

Define regions for measurement

Inner region distance in from edge: 1  $\mu\text{m} = 6$  pixels

Outer region distance out from edge: 1  $\mu\text{m} = 6$  pixels

Outer region width: 3  $\mu\text{m} = 19$  pixels

Positive classification

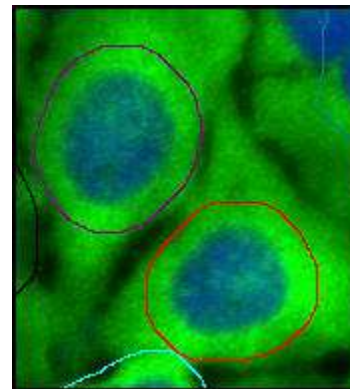
Minimum correlation coefficient: 0.45

Preview

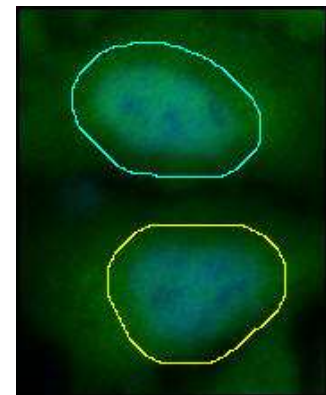
Configure Summary Log... Configure Data Log (Cells)...

Save Settings... Load Settings... Show Legend... Test Run Close

- **Positive classification:**
- Positive cells for this probe are those with nuclear staining. Negative cells for this probe are those with cytoplasmic staining.



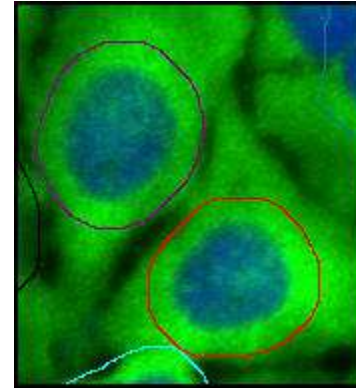
**Negative**  
(cytoplasmic)



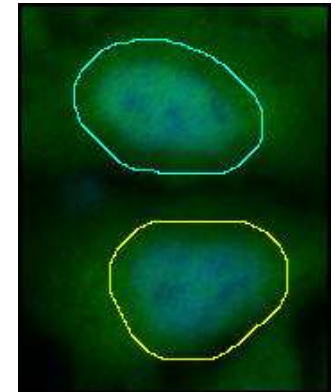
**Positive**  
(nuclear)

# Module Settings: Wavelength 2

- **Positive classification:**
- **Correlation Coefficient** - This is the Pearson's correlation coefficient of the pixel intensity of the probe of interest and the nuclear stain in the entire cell region (nucleus + gap + cytoplasm)
- This is typically the most robust method for classifying translocation
- 1.0 is perfect correlation (the two stains perfectly overlap)
- -1.0 is perfect anti-correlation (the two stains never overlap)
- 0 indicates the stains are independent



***Negative  
(cytoplasmic)***



***Positive  
(nuclear)***

# Module Settings: Additional wavelengths

- **Wavelength 3:**
- Follow the same steps to configure the settings for wavelength 3 (probe B in this example)

Configure Settings for Multi Wavelength Translocation ...

Number of wavelengths: 3

☐ Display result image: [None]

Adaptive Background Correction™ system

Compartments | **Probe A** | Probe B

Name: Probe B

W3 Source image: Texas Red

Legend color: Red

Define regions for measurement

Inner region distance in from edge: 1 µm = 6 pixels

Outer region distance out from edge: 1 µm = 6 pixels

Outer region width: 3 µm = 19 pixels

Positive classification

Minimum correlation coefficient: 0.3

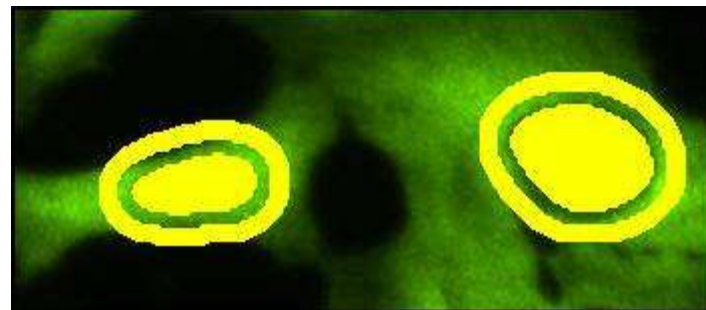
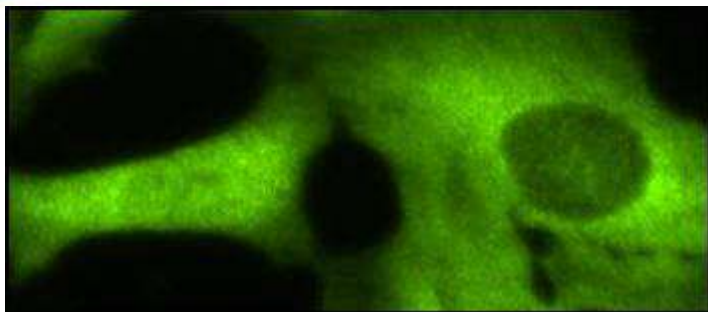
Preview

Configure Summary Log... Configure Data Log (Cells)...

Save Settings... Load Settings... Show Legend... Test Run Close



# Module Settings

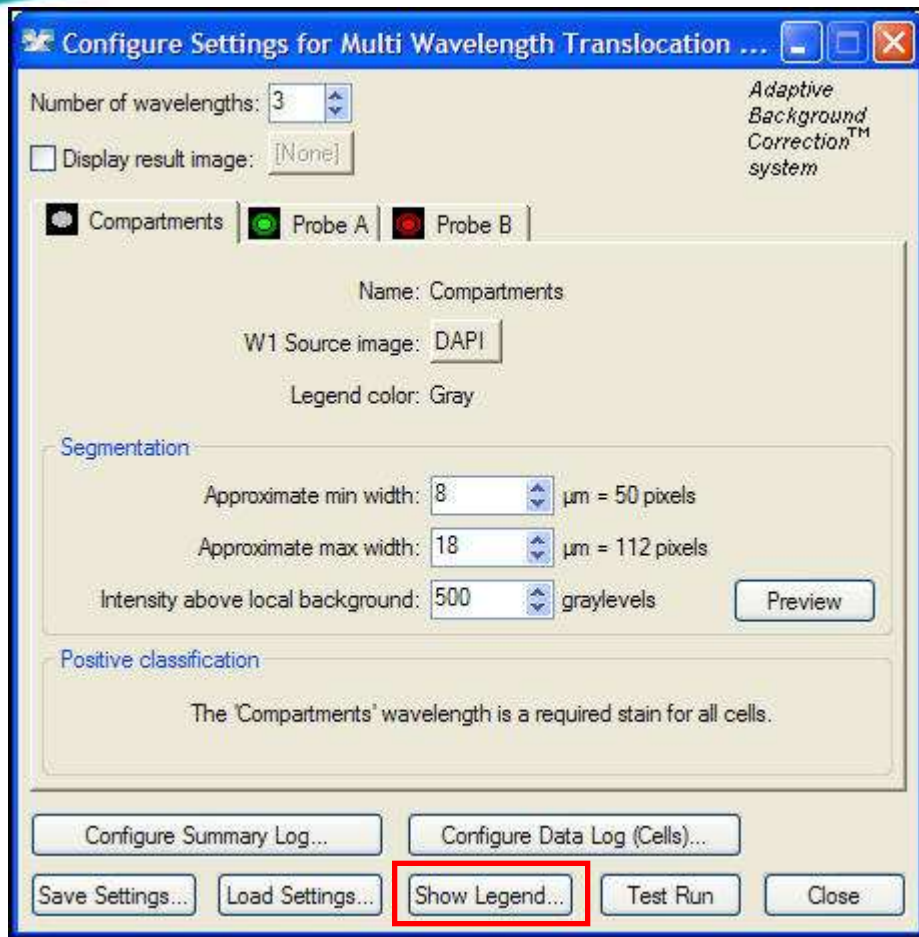


- **Setting Cell classification cutoffs:**
- Test module on positive and negative controls
- Use the interactive cellular results table to view the individual correlation or results for positive and negative cells
- An image showing both phenotypes makes it easy to compare results

Cell: Correlation Coefficient W2	Cell: Correlation Coefficient W3
-0.543586	-0.543574
-0.0849507	-0.084916
0.268204	0.268217
-0.872282	-0.872281
-0.618995	-0.618988
0.673702	0.673709
-0.635529	-0.635505
-0.852296	-0.852296
-0.262854	-0.262858
-0.277377	-0.277375
-0.859471	-0.859467
-0.103183	-0.103184
-0.409432	-0.40944
0.490158	0.49016
-0.0994213	-0.0994328
-0.760795	-0.760797
-0.508837	-0.508822
0.0343354	0.0343394
-0.780436	-0.780431
-0.532925	-0.532945
-0.681289	-0.681297
-0.692857	-0.692865
-0.717771	-0.717781



# Module Settings



- **Show Legend**
- Optionally name each profile



# Module Settings – General Settings

Configure Settings for Multi Wavelength Translocation ...

Number of wavelengths: 3

☐ Display result image: [None]

☒ Compartments ☒ Probe A ☒ Probe B

Name: Compartments

W1 Source image: DAPI

Legend color: Gray

Segmentation

Approximate min width: 8  $\mu\text{m}$  = 50 pixels

Approximate max width: 18  $\mu\text{m}$  = 112 pixels

Intensity above local background: 500 graylevels

Preview

Positive classification

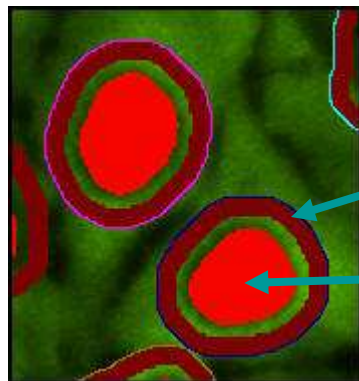
The 'Compartments' wavelength is a required stain for all cells.

Configure Summary Log... Configure Data Log (Cells)...

Save Settings... Load Settings... Show Legend... Test Run Close

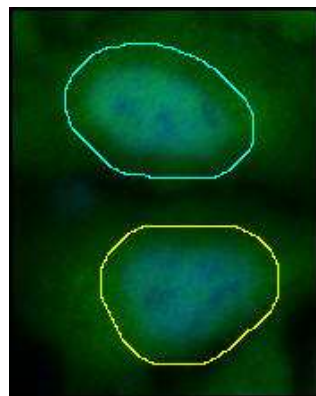
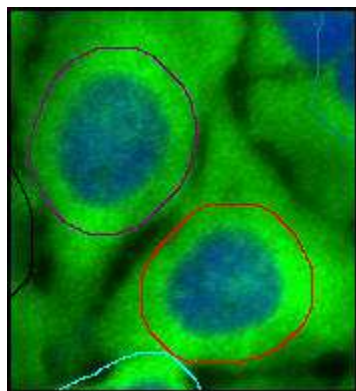
- **Configure Summary Log** – select site-by-site measurements
- **Configure Data Log** – select cell-by-cell measurements
- **Save Settings** – save analysis parameters to database
- **Load Settings** – load saved analysis parameters
- **Set to Defaults** – restore default analysis parameters
- **Test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

# Regions for Measurements



*Outer region*

*Inner region*



*Area for correlation coefficient*

- **Regions**
- **Outer region:** corresponds to cytoplasm (minus gap at boundary)
- **Inner region:** corresponds to nucleus (minus gap at boundary)
- **Marked area:** corresponds to inner + outer regions
- Area for **correlation coefficient:** inner region + gap + outer region
- Note that the regions are defined separately for each wavelength

# Summary Data (site-by-site measurements)

✓ Total Cells
✓ Mean Compartments Area
✓ Mean Integrated Compartments Intens
✓ Mean Average Compartments Intens
✓ % Positive W2
✓ % Positive W3
✓ Positive W2
✓ Positive W3
✓ Scoring Profile 1-
✓ Scoring Profile 12-
✓ Scoring Profile 1-3
✓ Scoring Profile 123
✓ Correlation Coefficient W2
✓ Correlation Coefficient W3
✓ Integrated Inner Intensity W2
✓ Integrated Inner Intensity W3
✓ Integrated Outer Intensity W2
✓ Integrated Outer Intensity W3
✓ Average Inner Intensity W2
✓ Average Inner Intensity W3
✓ Average Outer Intensity W2
✓ Average Outer Intensity W3
✓ Background Intensity W2
✓ Background Intensity W3
✓ Mean Marked Area W2
✓ Mean Marked Area W3
✓ Mean Integrated Marked Intensity W2
✓ Mean Integrated Marked Intensity W3
✓ Mean Average Marked Intensity W2
✓ Mean Average Marked Intensity W3

- **Total cells:** Total number of nuclei (cell count)
- **Mean Compartment Area:** The average nuclear area (in  $\mu\text{m}^2$ )
- **Mean Integrated Compartment Intensity:** The total pixel intensity of the nuclear stain over the nuclear area, divided by the number of cells
- **Mean Average Compartment Intensity:** The average pixel intensity of the nuclear stain over all the nuclear areas in the image



# Summary Data (site-by-site measurements)

✓ Total Cells
✓ Mean Compartments Area
✓ Mean Integrated Compartments Intens
✓ Mean Average Compartments Intens
✓ % Positive W2
✓ % Positive W3
✓ Positive W2
✓ Positive W3
✓ Scoring Profile 1--
✓ Scoring Profile 12-
✓ Scoring Profile 1-3
✓ Scoring Profile 123
✓ Correlation Coefficient W2
✓ Correlation Coefficient W3
✓ Integrated Inner Intensity W2
✓ Integrated Inner Intensity W3
✓ Integrated Outer Intensity W2
✓ Integrated Outer Intensity W3
✓ Average Inner Intensity W2
✓ Average Inner Intensity W3
✓ Average Outer Intensity W2
✓ Average Outer Intensity W3
✓ Background Intensity W2
✓ Background Intensity W3
✓ Mean Marked Area W2
✓ Mean Marked Area W3
✓ Mean Integrated Marked Intensity W2
✓ Mean Integrated Marked Intensity W3
✓ Mean Average Marked Intensity W2
✓ Mean Average Marked Intensity W3

- **Positive W2:** The number of cells positive for nuclear translocation in wavelength 2 (probe A) as defined by correlation with the nuclear stain over the cell region
- **Positive W3:** The number of cells positive for nuclear translocation in wavelength 3 (probe B) as defined by correlation with the nuclear stain over the cell region
- **% Positive W2:** The number of positive W2 cells divided by the total number of cells, times 100
- **% Positive W3:** The number of positive W3 cells divided by the total number of cells, times 100





# Summary Data (site-by-site measurements)

- ✓ Total Cells
- ✓ Mean Compartments Area
- ✓ Mean Integrated Compartments Intens
- ✓ Mean Average Compartments Intens
- ✓ % Positive W2
- ✓ % Positive W3
- ✓ Positive W2
- ✓ Positive W3
- ✓ Scoring Profile 1--
- ✓ Scoring Profile 12-
- ✓ Scoring Profile 1-3
- ✓ Scoring Profile 123
- ✓ Correlation Coefficient W2
- ✓ Correlation Coefficient W3
- ✓ Integrated Inner Intensity W2
- ✓ Integrated Inner Intensity W3
- ✓ Integrated Outer Intensity W2
- ✓ Integrated Outer Intensity W3
- ✓ Average Inner Intensity W2
- ✓ Average Inner Intensity W3
- ✓ Average Outer Intensity W2
- ✓ Average Outer Intensity W3
- ✓ Background Intensity W2
- ✓ Background Intensity W3
- ✓ Mean Marked Area W2
- ✓ Mean Marked Area W3
- ✓ Mean Integrated Marked Intensity W2
- ✓ Mean Integrated Marked Intensity W3
- ✓ Mean Average Marked Intensity W2
- ✓ Mean Average Marked Intensity W3

- **Scoring Profile 1--:** The number of cells negative for nuclear translocation in both wavelengths 2 and 3
- **Scoring Profile 12-:** The number of cells positive for nuclear translocation in wavelength 2 and negative for nuclear translocation in wavelength 3
- **Scoring Profile 1-3:** The number of cells negative for nuclear translocation in wavelength 2 and positive for nuclear translocation in wavelength 3
- **Scoring Profile 123:** The number of cells positive for nuclear translocation in both wavelengths 2 and 3

# Summary Data (site-by-site measurements)

- ✓ Total Cells
- ✓ Mean Compartments Area
- ✓ Mean Integrated Compartments Intens
- ✓ Mean Average Compartments Intens
- ✓ % Positive W2
- ✓ % Positive W3
- ✓ Positive W2
- ✓ Positive W3
- ✓ Scoring Profile 1--
- ✓ Scoring Profile 12-
- ✓ Scoring Profile 1-3
- ✓ Scoring Profile 123
- ✓ Correlation Coefficient W2
- ✓ Correlation Coefficient W3
- ✓ Integrated Inner Intensity W2
- ✓ Integrated Inner Intensity W3
- ✓ Integrated Outer Intensity W2
- ✓ Integrated Outer Intensity W3
- ✓ Average Inner Intensity W2
- ✓ Average Inner Intensity W3
- ✓ Average Outer Intensity W2
- ✓ Average Outer Intensity W3
- ✓ Background Intensity W2
- ✓ Background Intensity W3
- ✓ Mean Marked Area W2
- ✓ Mean Marked Area W3
- ✓ Mean Integrated Marked Intensity W2
- ✓ Mean Integrated Marked Intensity W3
- ✓ Mean Average Marked Intensity W2
- ✓ Mean Average Marked Intensity W3

- **Correlation Coefficient W2:**  
The Pearson's correlation coefficient between the nuclear stain and wavelength 2 over all of the pixels located in all of the cell regions (nuclei + gaps + cytoplasms) in the site
- **Correlation Coefficient W3:**  
The Pearson's correlation coefficient between the nuclear stain and wavelength 3 over all of the pixels located in all of the cell regions (nuclei + gaps + cytoplasms) in the site

# Summary Data (site-by-site measurements)

- ✓ Total Cells
- ✓ Mean Compartments Area
- ✓ Mean Integrated Compartments Intens
- ✓ Mean Average Compartments Intens
- ✓ % Positive W2
- ✓ % Positive W3
- ✓ Positive W2
- ✓ Positive W3
- ✓ Scoring Profile 1--
- ✓ Scoring Profile 12-
- ✓ Scoring Profile 1-3
- ✓ Scoring Profile 123
- ✓ Correlation Coefficient W2
- ✓ Correlation Coefficient W3
- ✓ Integrated Inner Intensity W2
- ✓ Integrated Inner Intensity W3
- ✓ Integrated Outer Intensity W2
- ✓ Integrated Outer Intensity W3
- ✓ Average Inner Intensity W2
- ✓ Average Inner Intensity W3
- ✓ Average Outer Intensity W2
- ✓ Average Outer Intensity W3
- ✓ Background Intensity W2
- ✓ Background Intensity W3
- ✓ Mean Marked Area W2
- ✓ Mean Marked Area W3
- ✓ Mean Integrated Marked Intensity W2
- ✓ Mean Integrated Marked Intensity W3
- ✓ Mean Average Marked Intensity W2
- ✓ Mean Average Marked Intensity W3

- **Integrated Inner Intensity W2:** The total pixel intensity of wavelength 2 (probe A) in all the inner W2 regions for the site after background subtraction (note this correlates with cell count)
- **Integrated Inner Intensity W3:** The total pixel intensity of wavelength 3 (probe B) in all the inner W3 regions for the site after background subtraction (note this correlates with cell count)
- **Integrated Outer Intensity W2:** The total pixel intensity of wavelength 2 (probe A) in all the outer W2 regions for the site after background subtraction (note this correlates with cell count)
- **Integrated Outer Intensity W3:** The total pixel intensity of wavelength 3 (probe B) in all the outer W3 regions for the site after background subtraction (note this correlates with cell count)



# Summary Data (site-by-site measurements)

- ✓ Total Cells
- ✓ Mean Compartments Area
- ✓ Mean Integrated Compartments Intens
- ✓ Mean Average Compartments Intens
- ✓ % Positive W2
- ✓ % Positive W3
- ✓ Positive W2
- ✓ Positive W3
- ✓ Scoring Profile 1--
- ✓ Scoring Profile 12--
- ✓ Scoring Profile 1-3
- ✓ Scoring Profile 123
- ✓ Correlation Coefficient W2
- ✓ Correlation Coefficient W3
- ✓ Integrated Inner Intensity W2
- ✓ Integrated Inner Intensity W3
- ✓ Integrated Outer Intensity W2
- ✓ Integrated Outer Intensity W3
- ✓ Average Inner Intensity W2
- ✓ Average Inner Intensity W3
- ✓ Average Outer Intensity W2
- ✓ Average Outer Intensity W3
- ✓ Background Intensity W2
- ✓ Background Intensity W3
- ✓ Mean Marked Area W2
- ✓ Mean Marked Area W3
- ✓ Mean Integrated Marked Intensity W2
- ✓ Mean Integrated Marked Intensity W3
- ✓ Mean Average Marked Intensity W2
- ✓ Mean Average Marked Intensity W3

- **Average Inner Intensity W2:** The average pixel intensity of wavelength 2 (probe A) in all the inner W2 regions for the site after background subtraction (independent of cell count)
- **Average Inner Intensity W3:** The average pixel intensity of wavelength 3 (probe B) in all the inner W3 regions for the site after background subtraction (independent of cell count)
- **Average Outer Intensity W2:** The average pixel intensity of wavelength 2 (probe A) in all the outer W2 regions for the site after background subtraction (independent of cell count)
- **Average Outer Intensity W3:** The average pixel intensity of wavelength 3 (probe B) in all the outer W3 regions for the site after background subtraction (independent of cell count)





# Summary Data (site-by-site measurements)

- ✓ Total Cells
- ✓ Mean Compartments Area
- ✓ Mean Integrated Compartments Intens
- ✓ Mean Average Compartments Intens
- ✓ % Positive W2
- ✓ % Positive W3
- ✓ Positive W2
- ✓ Positive W3
- ✓ Scoring Profile 1--
- ✓ Scoring Profile 12-
- ✓ Scoring Profile 1-3
- ✓ Scoring Profile 123
- ✓ Correlation Coefficient W2
- ✓ Correlation Coefficient W3
- ✓ Integrated Inner Intensity W2
- ✓ Integrated Inner Intensity W3
- ✓ Integrated Outer Intensity W2
- ✓ Integrated Outer Intensity W3
- ✓ Average Inner Intensity W2
- ✓ Average Inner Intensity W3
- ✓ Average Outer Intensity W2
- ✓ Average Outer Intensity W3
- ✓ Background Intensity W2
- ✓ Background Intensity W3
- ✓ Mean Marked Area W2
- ✓ Mean Marked Area W3
- ✓ Mean Integrated Marked Intensity W2
- ✓ Mean Integrated Marked Intensity W3
- ✓ Mean Average Marked Intensity W2
- ✓ Mean Average Marked Intensity W3

- **Background Intensity W2:** The average background pixel intensity of the wavelength 2 (probe A) image. This is the value that has been subtracted from other W2 intensity measurements.
- **Background Intensity W3:** The average background pixel intensity of the wavelength 3 (probe B) image. This is the value that has been subtracted from other W3 intensity measurements.
- **Mean Marked Area W2:** The total area covered by inner + outer regions as defined for wavelength 2, divided by the total cell count
- **Mean Marked Area W3:** The total area covered by inner + outer regions as defined for wavelength 3, divided by the total cell count



# Summary Data (site-by-site measurements)

- ✓ Total Cells
- ✓ Mean Compartments Area
- ✓ Mean Integrated Compartments Intens
- ✓ Mean Average Compartments Intens
- ✓ % Positive W2
- ✓ % Positive W3
- ✓ Positive W2
- ✓ Positive W3
- ✓ Scoring Profile 1--
- ✓ Scoring Profile 12-
- ✓ Scoring Profile 1-3
- ✓ Scoring Profile 123
- ✓ Correlation Coefficient W2
- ✓ Correlation Coefficient W3
- ✓ Integrated Inner Intensity W2
- ✓ Integrated Inner Intensity W3
- ✓ Integrated Outer Intensity W2
- ✓ Integrated Outer Intensity W3
- ✓ Average Inner Intensity W2
- ✓ Average Inner Intensity W3
- ✓ Average Outer Intensity W2
- ✓ Average Outer Intensity W3
- ✓ Background Intensity W2
- ✓ Background Intensity W3
- ✓ Mean Marked Area W2
- ✓ Mean Marked Area W3
- ✓ Mean Integrated Marked Intensity W2
- ✓ Mean Integrated Marked Intensity W3
- ✓ Mean Average Marked Intensity W2
- ✓ Mean Average Marked Intensity W3

- **Mean Integrated Marked Intensity W2:**  
The total pixel intensity of wavelength 2 (probe A) in the inner + outer W2 regions minus background, divided by the total cell count
- **Mean Integrated Marked Intensity W3:**  
The total pixel intensity of wavelength 3 (probe B) in the inner + outer W3 regions minus background, divided by the total cell count
- **Mean Average Marked Intensity W2:** The average pixel intensity of wavelength 2 (probe A) across all of the inner + outer W2 regions in the image minus background
- **Mean Average Marked Intensity W3:** The average pixel intensity of wavelength 3 (probe B) across all of the inner + outer W3 regions in the image minus background

# Cell Data (cell-by-cell measurements)

✓ Cell: Assigned Label #
✓ Cell: Scoring Profile
✓ Cell: Custom Profile Name
✓ Cell: Compartment Stained Area
✓ Cell: Compartment Integrated Intensity
✓ Cell: Compartment Mean Intensity
✓ Cell: Positive W2
✓ Cell: Positive W3
✓ Cell: Correlation Coefficient W2
✓ Cell: Correlation Coefficient W3
✓ Cell: Integrated Inner Intensity W2
✓ Cell: Integrated Inner Intensity W3
✓ Cell: Integrated Outer Intensity W2
✓ Cell: Integrated Outer Intensity W3
✓ Cell: Mean Inner Intensity W2
✓ Cell: Mean Inner Intensity W3
✓ Cell: Mean Outer Intensity W2
✓ Cell: Mean Outer Intensity W3
✓ Cell: Inner Area W2
✓ Cell: Inner Area W3
✓ Cell: Outer Area W2
✓ Cell: Outer Area W3
✓ Cell: Marked Area W2
✓ Cell: Marked Area W3
✓ Cell: Integrated Marked Intensity W2
✓ Cell: Integrated Marked Intensity W3
✓ Cell: Mean Marked Intensity W2
✓ Cell: Mean Marked Intensity W3

- **Cell: Assigned Label #** – Cell label number (1 through total cell number)
- **Cell: Scoring Profile** – The profile for this cell, i.e. "1—" or "12—" or "1-3" or "123"
- **Cell: Custom Profile Name** – The user-defined profile for this cell, as specified in the legend (e.g. "A positive", "Double positive")
- **Cell: Compartment Stained Area** – Total square microns of the nucleus
- **Cell: Compartment Integrated Intensity** – Total pixel intensity of the nuclear stain in the nucleus
- **Cell: Compartment Mean Intensity** – Average pixel intensity of the nuclear stain in the nucleus

# Cell Data (cell-by-cell measurements)

- ✓ Cell: Assigned Label #
- ✓ Cell: Scoring Profile
- ✓ Cell: Custom Profile Name
- ✓ Cell: Compartment Stained Area
- ✓ Cell: Compartment Integrated Intensity
- ✓ Cell: Compartment Mean Intensity
- ✓ Cell: Positive W2
- ✓ Cell: Positive W3
- ✓ Cell: Correlation Coefficient W2
- ✓ Cell: Correlation Coefficient W3
- ✓ Cell: Integrated Inner Intensity W2
- ✓ Cell: Integrated Inner Intensity W3
- ✓ Cell: Integrated Outer Intensity W2
- ✓ Cell: Integrated Outer Intensity W3
- ✓ Cell: Mean Inner Intensity W2
- ✓ Cell: Mean Inner Intensity W3
- ✓ Cell: Mean Outer Intensity W2
- ✓ Cell: Mean Outer Intensity W3
- ✓ Cell: Inner Area W2
- ✓ Cell: Inner Area W3
- ✓ Cell: Outer Area W2
- ✓ Cell: Outer Area W3
- ✓ Cell: Marked Area W2
- ✓ Cell: Marked Area W3
- ✓ Cell: Integrated Marked Intensity W2
- ✓ Cell: Integrated Marked Intensity W3
- ✓ Cell: Mean Marked Intensity W2
- ✓ Cell: Mean Marked Intensity W3

- **Cell: Positive W2** – Classification of the cell by correlation coefficient as positive for translocation (nuclear staining) in wavelength 2 (value of 1) or negative for translocation (cytoplasmic staining) in wavelength 2 (value of 0)
- **Cell: Positive W3** – Classification of the cell by correlation coefficient as positive for translocation (nuclear staining) in wavelength 3 (value of 1) or negative for translocation (cytoplasmic staining) in wavelength 3 (value of 0)
- **Cell: Correlation Coefficient W2** – The Pearson's correlation coefficient between the intensities of the nuclear stain and wavelength 2 (probe A) for all pixels in the cell region for W2 (nucleus + gap + cytoplasm). The value ranges from -1 (anti-correlated) to 1 (correlated).
- **Cell: Correlation Coefficient W3** – The Pearson's correlation coefficient between the intensities of the nuclear stain and wavelength 3 (probe B) for all pixels in the cell region for W3 (nucleus + gap + cytoplasm). The value ranges from -1 (anti-correlated) to 1 (correlated).

# Cell Data (cell-by-cell measurements)

- ✓ Cell: Assigned Label #
- ✓ Cell: Scoring Profile
- ✓ Cell: Custom Profile Name
- ✓ Cell: Compartment Stained Area
- ✓ Cell: Compartment Integrated Intensity
- ✓ Cell: Compartment Mean Intensity
- ✓ Cell: Positive W2
- ✓ Cell: Positive W3
- ✓ Cell: Correlation Coefficient W2
- ✓ Cell: Correlation Coefficient W3
- ✓ Cell: Integrated Inner Intensity W2
- ✓ Cell: Integrated Inner Intensity W3
- ✓ Cell: Integrated Outer Intensity W2
- ✓ Cell: Integrated Outer Intensity W3
- ✓ Cell: Mean Inner Intensity W2
- ✓ Cell: Mean Inner Intensity W3
- ✓ Cell: Mean Outer Intensity W2
- ✓ Cell: Mean Outer Intensity W3
- ✓ Cell: Inner Area W2
- ✓ Cell: Inner Area W3
- ✓ Cell: Outer Area W2
- ✓ Cell: Outer Area W3
- ✓ Cell: Marked Area W2
- ✓ Cell: Marked Area W3
- ✓ Cell: Integrated Marked Intensity W2
- ✓ Cell: Integrated Marked Intensity W3
- ✓ Cell: Mean Marked Intensity W2
- ✓ Cell: Mean Marked Intensity W3

- **Cell: Integrated Inner Intensity W2** – The total pixel intensity of wavelength 2 (probe A) in the inner W2 region minus background
- **Cell: Integrated Inner Intensity W3** – The total pixel intensity of wavelength 3 (probe B) in the inner W3 region minus background
- **Cell: Integrated Outer Intensity W2** – The total pixel intensity of wavelength 2 (probe A) in the outer W2 region minus background
- **Cell: Integrated Outer Intensity W3** – The total pixel intensity of wavelength 3 (probe B) in the outer W3 region minus background



# Cell Data (cell-by-cell measurements)

- ✓ Cell: Assigned Label #
- ✓ Cell: Scoring Profile
- ✓ Cell: Custom Profile Name
- ✓ Cell: Compartment Stained Area
- ✓ Cell: Compartment Integrated Intensity
- ✓ Cell: Compartment Mean Intensity
- ✓ Cell: Positive W2
- ✓ Cell: Positive W3
- ✓ Cell: Correlation Coefficient W2
- ✓ Cell: Correlation Coefficient W3
- ✓ Cell: Integrated Inner Intensity W2
- ✓ Cell: Integrated Inner Intensity W3
- ✓ Cell: Integrated Outer Intensity W2
- ✓ Cell: Integrated Outer Intensity W3
- ✓ Cell: Mean Inner Intensity W2
- ✓ Cell: Mean Inner Intensity W3
- ✓ Cell: Mean Outer Intensity W2
- ✓ Cell: Mean Outer Intensity W3
- ✓ Cell: Inner Area W2
- ✓ Cell: Inner Area W3
- ✓ Cell: Outer Area W2
- ✓ Cell: Outer Area W3
- ✓ Cell: Marked Area W2
- ✓ Cell: Marked Area W3
- ✓ Cell: Integrated Marked Intensity W2
- ✓ Cell: Integrated Marked Intensity W3
- ✓ Cell: Mean Marked Intensity W2
- ✓ Cell: Mean Marked Intensity W3

- **Cell: Mean Inner Intensity W2** – The average pixel intensity of wavelength 2 (probe A) in the inner W2 region minus background
- **Cell: Mean Inner Intensity W3** – The average pixel intensity of wavelength 3 (probe B) in the inner W3 region minus background
- **Cell: Mean Outer Intensity W2** – The average pixel intensity of wavelength 2 (probe A) in the outer W2 region minus background
- **Cell: Mean Outer Intensity W3** – The average pixel intensity of wavelength 3 (probe B) in the outer W3 region minus background

# Cell Data (cell-by-cell measurements)

- ✓ Cell: Assigned Label #
- ✓ Cell: Scoring Profile
- ✓ Cell: Custom Profile Name
- ✓ Cell: Compartment Stained Area
- ✓ Cell: Compartment Integrated Intensity
- ✓ Cell: Compartment Mean Intensity
- ✓ Cell: Positive W2
- ✓ Cell: Positive W3
- ✓ Cell: Correlation Coefficient W2
- ✓ Cell: Correlation Coefficient W3
- ✓ Cell: Integrated Inner Intensity W2
- ✓ Cell: Integrated Inner Intensity W3
- ✓ Cell: Integrated Outer Intensity W2
- ✓ Cell: Integrated Outer Intensity W3
- ✓ Cell: Mean Inner Intensity W2
- ✓ Cell: Mean Inner Intensity W3
- ✓ Cell: Mean Outer Intensity W2
- ✓ Cell: Mean Outer Intensity W3
- ✓ Cell: Inner Area W2
- ✓ Cell: Inner Area W3
- ✓ Cell: Outer Area W2
- ✓ Cell: Outer Area W3
- ✓ Cell: Marked Area W2
- ✓ Cell: Marked Area W3
- ✓ Cell: Integrated Marked Intensity W2
- ✓ Cell: Integrated Marked Intensity W3
- ✓ Cell: Mean Marked Intensity W2
- ✓ Cell: Mean Marked Intensity W3

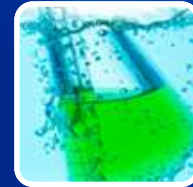
- **Cell: Inner Area W2** – Total square microns in the inner W2 region
- **Cell: Inner Area W3** – Total square microns in the inner W3 region
- **Cell: Outer Area W2** – Total square microns in the outer W2 region
- **Cell: Outer Area W3** – Total square microns in the outer W3 region
- **Cell: Marked Area W2** – The sum of the inner and outer W2 areas
- **Cell: Marked Area W3** – The sum of the inner and outer W3 areas

# Cell Data (cell-by-cell measurements)

- ✓ Cell: Assigned Label #
- ✓ Cell: Scoring Profile
- ✓ Cell: Custom Profile Name
- ✓ Cell: Compartment Stained Area
- ✓ Cell: Compartment Integrated Intensity
- ✓ Cell: Compartment Mean Intensity
- ✓ Cell: Positive W2
- ✓ Cell: Positive W3
- ✓ Cell: Correlation Coefficient W2
- ✓ Cell: Correlation Coefficient W3
- ✓ Cell: Integrated Inner Intensity W2
- ✓ Cell: Integrated Inner Intensity W3
- ✓ Cell: Integrated Outer Intensity W2
- ✓ Cell: Integrated Outer Intensity W3
- ✓ Cell: Mean Inner Intensity W2
- ✓ Cell: Mean Inner Intensity W3
- ✓ Cell: Mean Outer Intensity W2
- ✓ Cell: Mean Outer Intensity W3
- ✓ Cell: Inner Area W2
- ✓ Cell: Inner Area W3
- ✓ Cell: Outer Area W2
- ✓ Cell: Outer Area W3
- ✓ Cell: Marked Area W2
- ✓ Cell: Marked Area W3
- ✓ Cell: Integrated Marked Intensity W2
- ✓ Cell: Integrated Marked Intensity W3
- ✓ Cell: Mean Marked Intensity W2
- ✓ Cell: Mean Marked Intensity W3

- **Cell: Integrated Marked Intensity W2**  
– The total pixel intensity of wavelength 2 (probe A) in the inner + outer W2 areas (minus background)
- **Cell: Integrated Marked Intensity W3**  
– The total pixel intensity of wavelength 3 (probe B) in the inner + outer W3 areas (minus background)
- **Cell: Mean Marked Intensity W2** – The average pixel intensity of wavelength 2 (probe A) in the inner + outer W2 areas (minus background)
- **Cell: Mean Marked Intensity W3** – The average pixel intensity of wavelength 3 (probe B) in the inner + outer W3 areas (minus background)





**Thank You**