



MetaXpress® 6 Software Guide

Application Modules: Mitotic Index



Date Revised 02/03/16 Version B

Index

- Index
- [Mitotic Index Application Module Overview](#)
- Defining Application Module Settings
 - [Select an Image & Algorithm](#)
 - [Define Width of Objects](#)
 - [Define Intensity](#)
- [Module Buttons](#)
- [Configuring Summary or Data \(Cell\) Logs](#)
- Explanation of Measurement Outputs
 - [Summary Log \(Image Measurements\)](#)
 - [Data Log \(Cell Measurements\)](#)
- [Support Resources](#)



Mitotic Index Application 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 (a.k.a. phospho-histone H3). The DNA stain labels all cells and the mitotic cells are labeled only with the mitotic-specific stain.



Module Settings: Selecting an Image & Algorithm

All nuclei (W1 Source image)

- Select the image for all nuclei.

Mitotic staining (W2 Source image)

- Select the image for the mitotic marker.

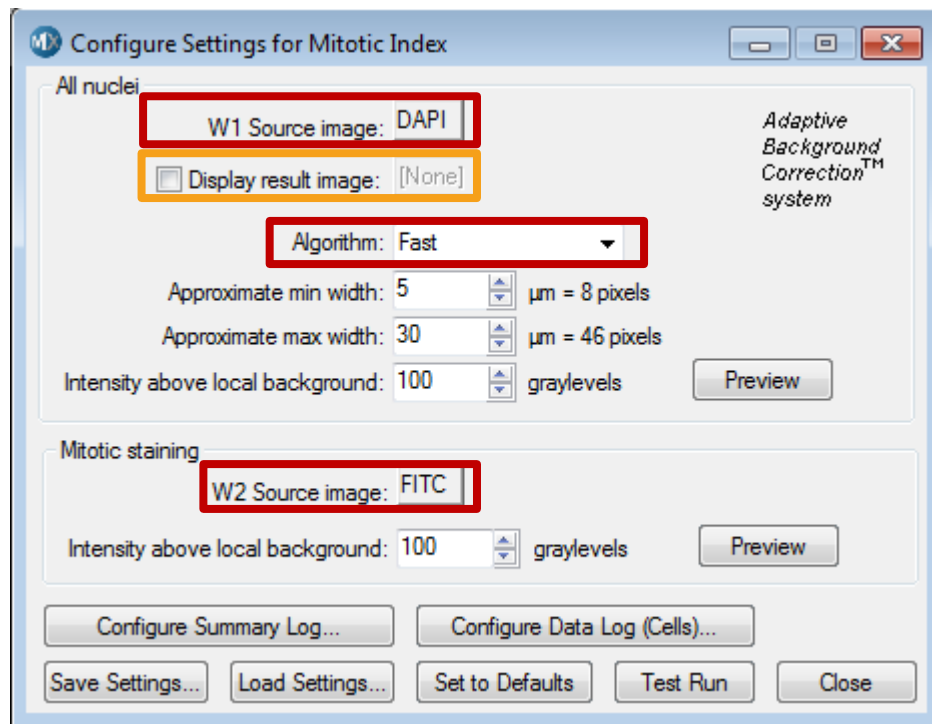
**NOTE* Do not choose images with “HTS” in the name*

Display result image:

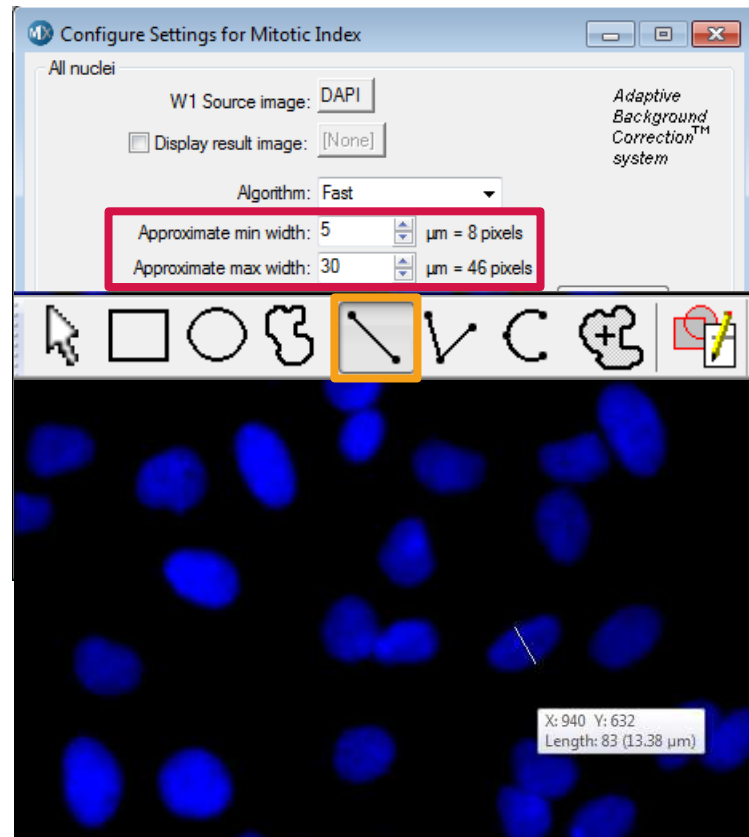
- Leave **Display result image** deselected (this is generally only used when creating a journal).

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**.



Module Settings: Defining the Size of Objects

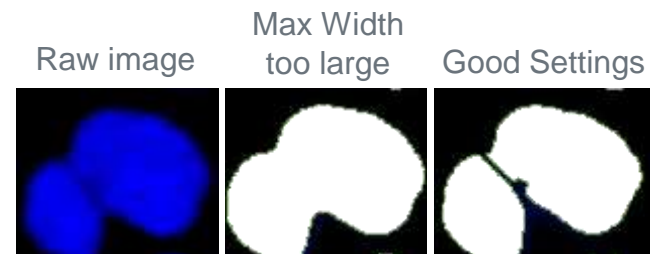


Approximate min and max widths

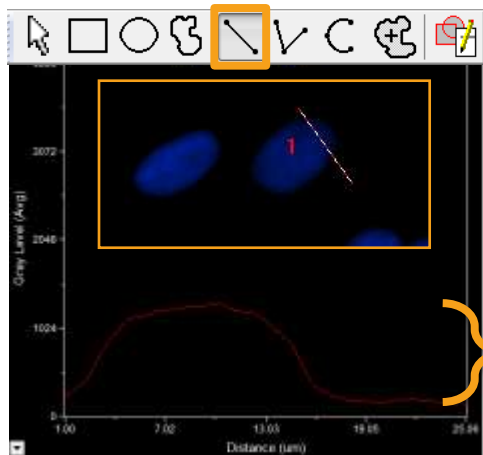
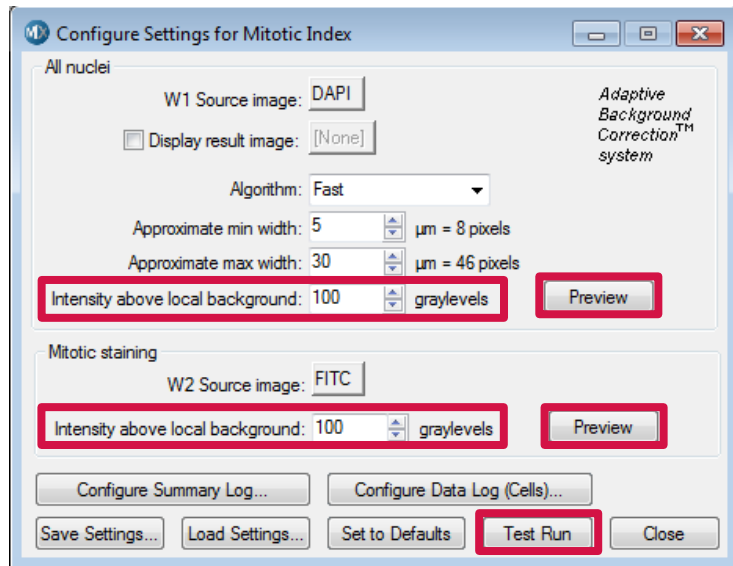
- Select the single line region from **Regions Tools**.
- Click and drag across a representative small and large object; a tooltip will show the length of the line. (single click to start a line, single click to terminate).

**NOTE* Do not click the image again. This will cause the tooltip to disappear. If the tooltip disappears, repeat the drawing procedure.*

- The width is the short axis of a object (in μm).
- Much smaller or much larger cells will be ignored.
- Click **Apply** and adjust values as necessary
- Repeat for all desired objects (all nuclei, mitotic staining)



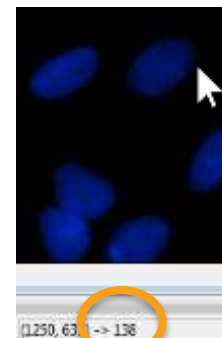
Module Settings: Defining the Intensity



Use the difference in intensity values to set **Intensity above local background**

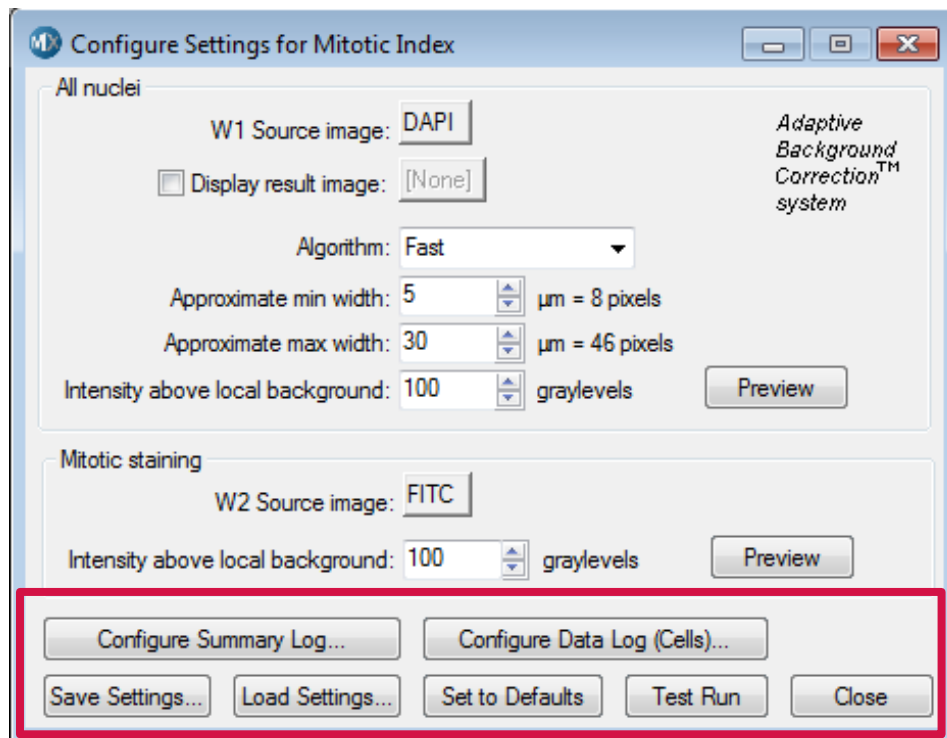
Intensity above local background

- Draw a line over the dimmest object of interest that covers both the object and background using the Single line tool
- In the main menu, select **Measure > Intensities > Linescan**. This will display a histogram of intensity values; or simply mouse over the dimmest object of interest and the background and view the intensity values.
- For **Fast algorithms**, set **Intensity above local background** to half (or less) of the difference in intensity between an object and background.
 - Example: $1180 - 138 = 1042 / 2 = 521$ (or less)
- For **Standard algorithms**, set this value slightly lower than the difference
 - Example: $1180 - 138 = 1042 - 100$. Start setting to 942
- Click on **Preview** to test individual wavelength or **Test Run** to test all settings based on the W1 and W2 markers. Adjust values as necessary by changing one parameter at a time.



Mouse over the dimmest object and outside to measure intensity values. Use the difference to set **Intensity above local background**

Module Buttons



Configure Summary Log: Select image-by-image measurements.

Configure Data Log: Select cell-by-cell measurements.

Save Settings: Save application module settings.

Load Settings: Load saved application module settings.

Set to Defaults: Restore default application module settings.

Test Run: Test all settings together and display cell-by-cell results for the displayed image.



Configuring Summary or Data (Cell) Logs

Double click on a measurement to select or deselect it for logging into the database.

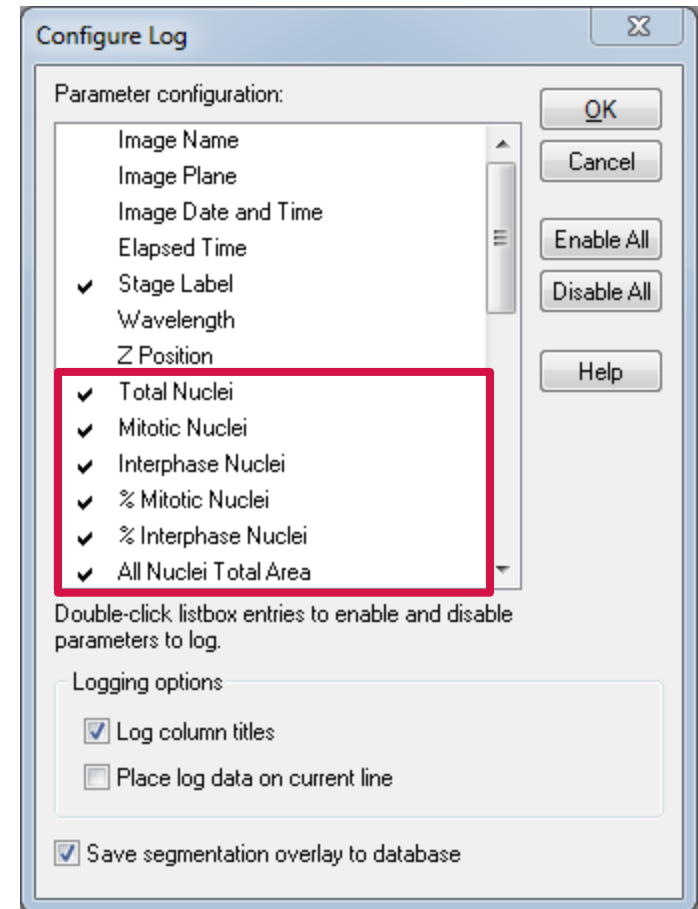
- ✓ Indicates a parameter that will be logged into the database (or Excel/text file log if open)

Log column titles: Does not affect database logging. If you have an Excel or text file log open, this records the parameter name as the column header for easy review. We recommend **enabling** this option.

Place log data on current line: Does not affect database logging. If you have an Excel or text file log open, this records the data into the last row used, to the right of the previous data. We recommend **disabling** this option to ensure that new data is recorded into a new row.

Save segmentation overlay to database: A mask (or binary image) is temporarily created for every raw image analyzed on a plate (see Transfluor overview for an example of a mask). This option allows you to save these masks to the database. We recommend **enabling this option for assay development and disabling it for screening.**

- Pro: Allows you to quickly review your segmentation results after analysis has been run across the entire plate
- Con: These masks take up a significant amount of space in a database, which may be limited in size. Saving the masks may also slow down analysis.



Configure Summary Log (Image 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

- **Total Nuclei:** Total number of cells determined from wavelength 1.
- **Mitotic Nuclei:** Total number of nuclei determined from wavelength 1 that were positive in wavelength 2.
- **Interphase Nuclei:** Total number of nuclei determined from wavelength 1 that were negative in wavelength 2.
- **% Mitotic Nuclei:** $100 * \# \text{ Mitotic Nuclei} / \# \text{ Total Nuclei}$.
- **% Interphase Nuclei:** $100 * \# \text{ Interphase Nuclei} / \# \text{ Total Nuclei}$.

Configure Summary Log (Image 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

- **All Nuclei Total Area:** Total μm^2 in wavelength 1 of All Nuclei.
- **All Nuclei Mean Area:** All Nuclei Total Area/#All 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.

Configure Summary Log (Image 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 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.

Configure Summary Log (Image 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 in wavelength 1 in all Interphase nuclei.
- **Interphase Mean Area:** Interphase Total Area/Total # 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.

Configure Data Log (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

- **Cell: Assigned Label #:** Cell label number (1 through total cell number).
- **Cell: Classification:** Labels cell as Interphase or Mitotic.
- **Cell: Total Area:** Total μm^2 in the nucleus of a given cell.
- **Cell: W1 Integrated Intensity:** Summed grayscale values in wavelength 1 for this nucleus.
- **Cell: W1 Average Intensity:** W1 Integrated Intensity/Total (Pixel) Area for this nucleus.
- **Cell: W2 Integrated Intensity:** Summed grayscale values in wavelength 2 for this nucleus.
- **Cell: W2 Average Intensity:** W2 Integrated Intensity/Total (Pixel) Area for this nucleus.

Support Resources

- F1 / HELP within MetaXpress® Software
- Support and Knowledge Base: <http://mdc.custhelp.com/>
- User Forum: <http://metamorph.moleculardevices.com/forum/>
- Request Support: <http://mdc.custhelp.com/app/ask>
- Technical Support can also be reached by telephone:
 - 1 (800) 635-5577
 - Select options for Tech Support → Cellular Imaging Products → ImageXpress Instruments





ADVANCING PROTEIN AND CELL BIOLOGY