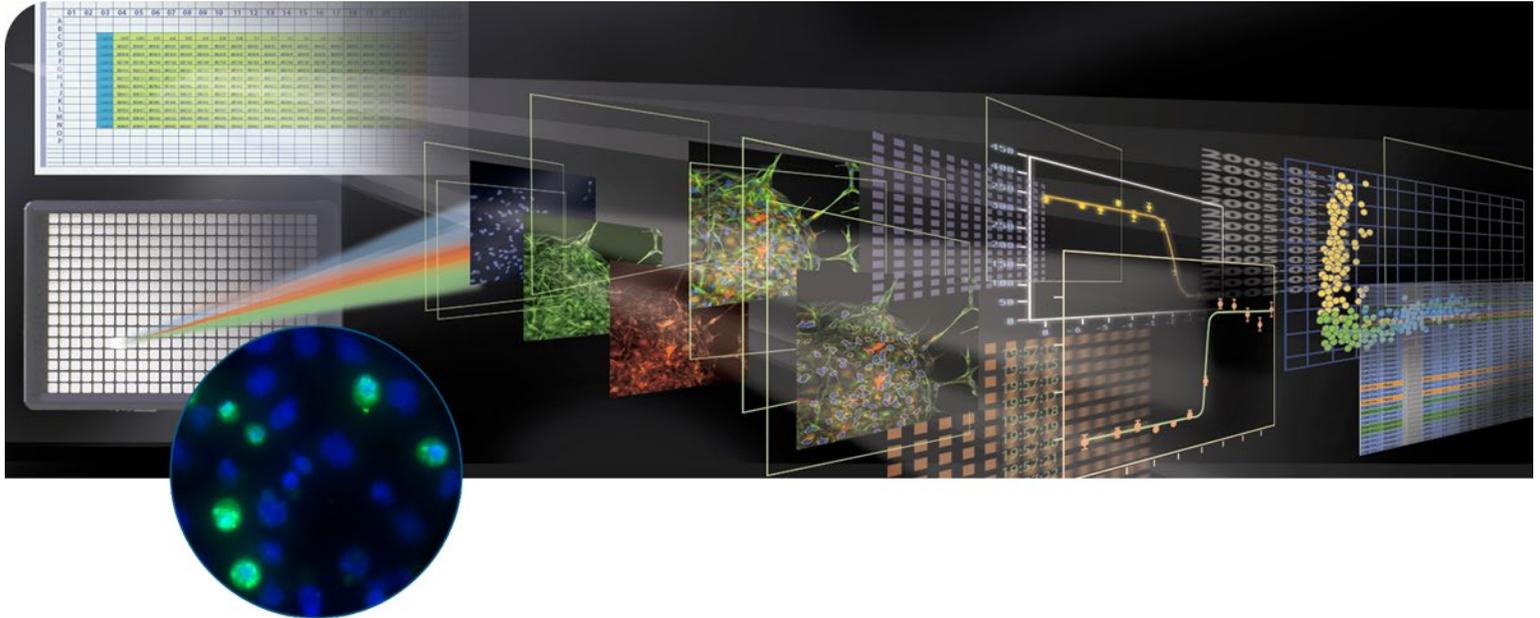


MetaXpress Software Mitotic Index Application Module

ANALYSIS SOFTWARE DROP-IN FOR METAXPRESS SOFTWARE



→ QUANTITATION OF MITOTIC AND INTERPHASE CELLS

→ MULTI-PARAMETER ANALYSIS OF TWO PROBES

→ ADAPTIVE BACKGROUND CORRECTION FOR IMPROVED SEGMENTATION

→ CELL-BY-CELL MEASUREMENTS AND DATA LOGGING

Cell-based assays for cell cycle status have become critical for oncology drug discovery programs. For example, potential anti-cancer therapeutics rely on arresting mitosis in cancerous cells to prevent uncontrolled proliferation.

In some forms of cancer, progression through mitosis can be stopped by merely disrupting the mitotic spindle. In others, the feedback is missing and cells proceed through mitosis without proper chromosome segregation and accumulate mutations. Cell-based assays provide an efficient method for modeling the effects of drugs on the cell cycle and quantitating the accumulation of cells in mitosis.

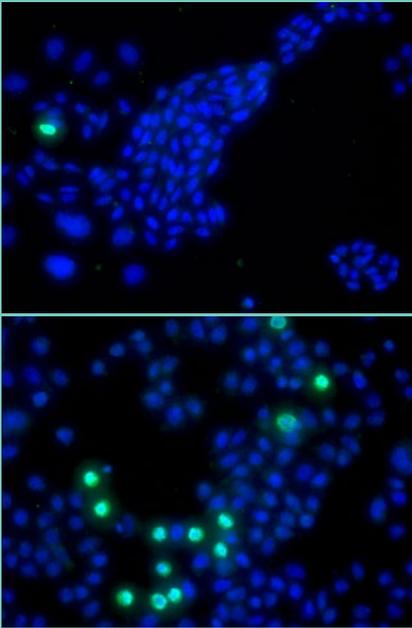
The Mitotic Index Application Module for MetaXpress® Software from Molecular Devices is designed for the quantitative discrimination of mitotic and interphase cells.

Cells are labeled with a DNA stain and a fluorescent mitosis-specific marker, such as immunofluorescence staining for Histone 3 S10 phosphorylation. Two different wavelengths are acquired and the images are analyzed with the Mitotic Index module.

A simple interface minimizes setup efforts and settings can be saved for future use. Segmentation parameters are set for each wavelength detection and the assay is run for a site, selected wells, the entire plate or multiple plates.

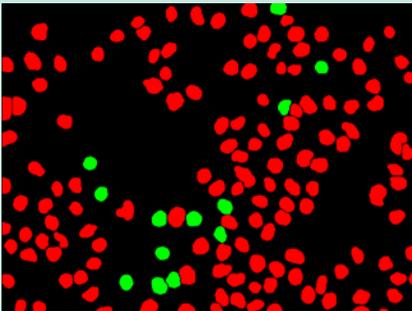
The module utilizes Adaptive Background Correction (ABC) which adapts the detection algorithm to the local intensity ranges between and within cells to provide the most robust segmentation available in an image-based screening system. ABC enables probe detection even with highly variable background fluorescence within a single image.

Multiple Probe Acquisition



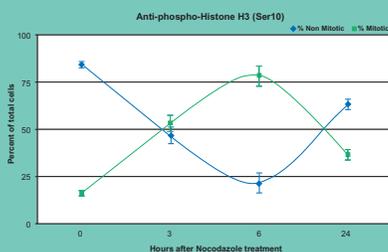
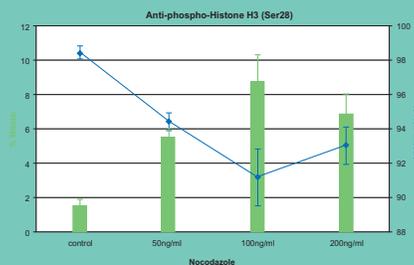
CHO-K1 cells treated with Nocadazole for 18 hours before staining with anti-phospho-Histone H3 (Ser28) and acquired with Molecular Devices' Discovery-1™ System. Top: control, bottom: 50 ng/mL Nocodazole.

Image Analysis



The Mitotic Index module identifies mitotic cells (green) and interphase cells (red).

Export Data for Further Analysis



EASY CONFIGURATION FOR ANALYSIS

1. Select the nuclear stain image
2. Specify the cell size range
3. Set the intensity above local background
4. Select the mitotic nuclear stain image
5. Set the intensity above local background
6. Optionally choose the reporting parameters

INTERACTIVE DATA DISPLAY

Once the analysis is run, the Cellular Results table allows you to interactively view individual cells' data. Clicking a cell in the image highlights the data for the selected cell in the table.

Cell Mitotic Classification	Cell Total Area	Cell W1 Integrated Intensity	Cell W1 Average Intensity	Cell W2 Integrated Intensity	Cell W2 Average Intensity	
43	Interphase	168.643	89009	211.423	70810	168.195
44	Interphase	169.444	89187	210.844	77830	183.995
45	Mitotic	171.046	114743	268.719	206074	482.609
46	Mitotic	171.046	148726	348.304	268296	628.328
47	Interphase	171.046	105017	245.941	73514	172.164
48	Interphase	171.847	93174	217.189	78084	182.014
49	Interphase	172.649	124008	287.722	85750	198.956
50	Interphase	173.049	105020	243.102	70969	164.28
51	Interphase	173.45	108178	249.834	69106	159.598
52	Mitotic	176.254	141599	321.816	209759	476.725
53	Mitotic	177.456	141173	318.675	325994	734.975
54	Interphase	179.458	111718	249.371	69390	154.799
55	Interphase	179.859	109715	244.354	79252	176.508
56	Interphase	182.663	91144	199.877	71049	155.809
57	Interphase	182.663	104651	229.498	80520	176.579
58	Interphase	184.265	95559	207.737	70712	153.722
59	Interphase	185.066	97737	211.552	71664	155.117

CUSTOMIZATION THROUGH MACROS

MetaXpress Software is seamlessly integrated with the power and flexibility of MetaMorph® Software and its sophisticated and powerful macros that record and perform a series of tasks without the need for a programming language.

VALIDATED DATA

Development of application modules includes research and testing with a library of in-house and third-party data sets.

MULTI-PARAMETER ANALYSIS

The application module generates a number of field or cell-by-cell parameters, including:

- Count of mitotic and non-mitotic nuclei
- Percentage of mitotic and non-mitotic nuclei
- Total and mean nuclei area of mitotic and non-mitotic cells
- Wavelength-specific integrated and average intensities of mitotic and non-mitotic nuclei

ORDERING INFORMATION

Mitotic Index Application Module for MetaXpress Software

Part Number: 9500-0036

FOR MORE INFORMATION

For more details, please visit

www.moleculardevices.com/mx

SALES OFFICES

- USA & Canada +1-800-635-5577
- Brazil +55-11-3616-6607
- China (Beijing) +86-10-6410-8669
- China (Shanghai) +86-21-6887-8820
- Germany 00800 665 32860
- Japan (Osaka) +81-6-7174-8831
- Japan (Tokyo) +81-3-6360-5260
- South Korea +82-2-3471-9531
- United Kingdom +44-118-944-8000

Check our web site for a current listing of our worldwide distributors.

www.moleculardevices.com

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

The trademarks used herein are the property of Molecular Devices, LLC or their respective owners.

Specifications subject to change without notice.