

ImageXpress Nano Automated Imaging System

BENEFITS

- Integration into MDCStore
 Data Management tool for easy access to needed files
- Bidirectional image viewer facilitates data quality control by connecting images to numerical data
- Multi-parametric analysis and visualization
- Dose response analysis and curve fitting with over 30 different predefined functions

Acquire

Widefield screening with automated microscopy allows researchers to scale up their experiments to multi-slide or microplate-level evaluations. The ImageXpress® Nano Automated Imaging System features a long life, solid state, light engine, and optics to reliably deliver the right assay sensitivity.

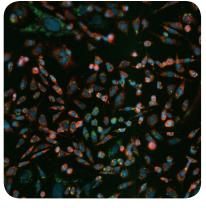
Analyze

Expedite image analysis and data processing with application modules that allow you to quickly and easily analyze your data. Our modular toolbox set up hundreds of routinely used HCS assays whether you're a novice or an expert. MetaXpress® software application modules give you an interactive environment to quantify 2D and 3D measurements including size, intensities, and distance.

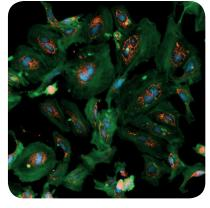
Intelligent segmentation

The MetaXpress software application modules are equipped with a proprietary adaptive background correction algorithm, which adjusts image segmentation to the local intensity ranges and shape features within and between cells. This capability gives you the most robust segmentation available in an image-based screening system and improves detection on both phase contrast and fluorescent labeled images.

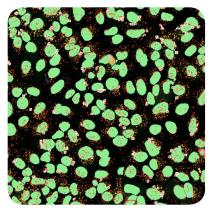




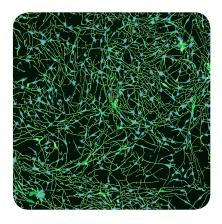
Autophagy



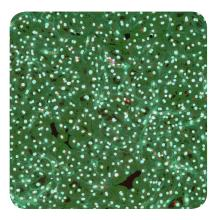
3 color toxicity



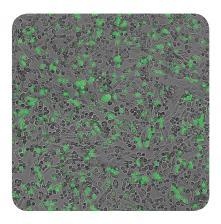
Intracellular vesicle mask



Neurite tracing



Nuclei and cytoplasm mask



Transfection efficiency with overlay

Software features

- Integration into MDCStore[™] Data Management tool for easy access to needed files
- Bidirectional image viewer that facilitates data quality control by connecting images to numerical data
- Multi-parametric analysis and visualization
- Dose response analysis and curve fitting with over 30 different predefined functions

System features

- · Widefield imaging
- Large field of view (A single 4x image captures whole 384 well)
- Solid state light engine with expected lifetime of 20,000 hours
- Acquire from a variety of multi-well plate formats and large format samples (e.g. slides)
- 4 position software-controlled objective changer
- 5 position software-controlled filter cube changer
- High speed laser autofocus technology
- Fully automated X, Y, and Z stages with resolution better than 25 nm

Options

- Transmitted light for label free brightfield imaging
- Cell incubation (temperature, humidity, CO₂) with environmental control feature
- Desktop computer with pre-installed software
- Objectives: 2x to 60x magnification
- · Select from a variety of filter cubes

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.



ImageXpress® Nano Automated
Imaging System

