

NEW ImageXpress Micro 4 High-Content Imaging System

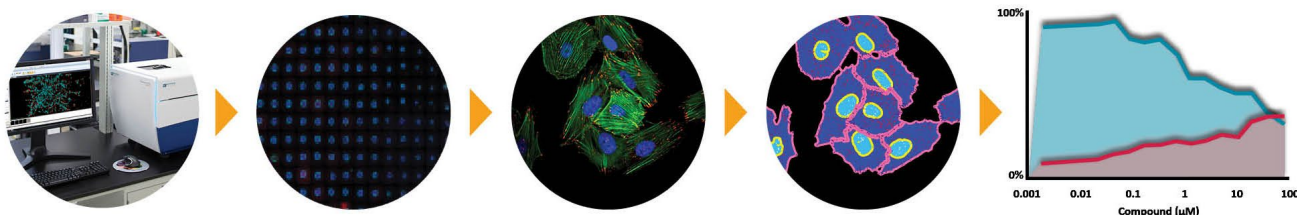
Position your research—with breakthrough, intelligent imaging solutions

BENEFITS

- Capture excellent images today with the capability to upgrade to confocal in the field if needed
- Tackle a wide range of assays with objectives ranging from 1x to 100x, and a configurable 5-position filter cube changer
- Acquire statistically relevant data faster with large field of view and >3 log dynamic range
- Expand your research capabilities with transmitted light, liquid handling, and environmental control options

The ImageXpress® Micro 4 High-Content Imaging System represents the culmination of four generations of imaging expertise. The latest agile design allows you to boost your research with this faster-than-ever system, while providing the option to upgrade to confocal in the future to align with your research needs.

Combined with MetaXpress® High-Content Image Acquisition and Analysis Software, the ImageXpress Micro 4 system provides you with an intelligent, multi-dimensional, high-throughput screening solution to help you discover your next landmark scientific breakthrough.



Combined with MetaXpress software, the ImageXpress Micro 4 system is a complete solution that enables you to interpret your images, understand your data, and explore new ideas.

Applications

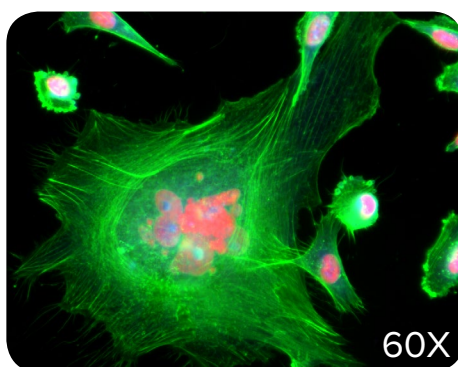
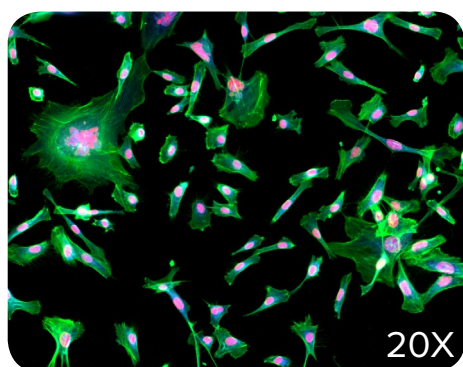
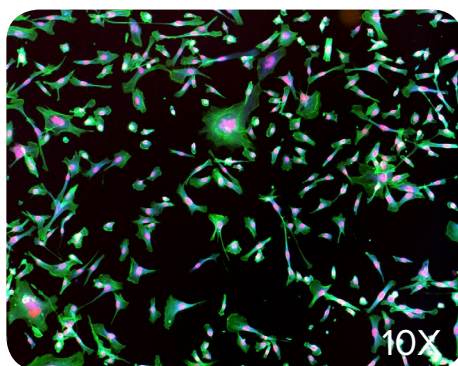
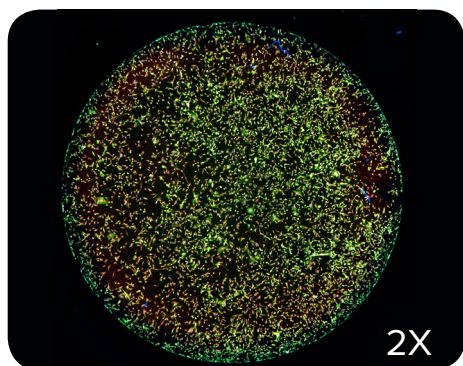
Capture images from Apoptosis to Zebrafish with the ImageXpress Micro 4 system configured to suit your specific biological needs.

We offer unlimited configurability with:

- User exchangeable filter cubes enabling multiplexed assays
- The widest range of objective lenses to capture a variety of cellular and intracellular events

- Environmental control for live cell monitoring
- Transmitted light for stain-free imaging
- Fluidics options for kinetic cellular response studies

Built on over 30 years of cell-based imaging experience, the ImageXpress Micro 4 system is engineered for performance to reliably accelerate the pace of your research.



Automatically acquire images 1-100x objective magnification in slides and multi-well plates up to 1536-well format. Acquire >210,000 wells per day at 4x magnification with the system's unique multi-well crop feature. A 4x objective captures 11 mm² of a 384-well plate to be imaged in one field-of-view while increasing magnifications allows the resolution of fine detail for analysis of subcellular events.

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.



Key features

- Widefield imaging with digital confocal imaging
- Field upgradeable to AgileOptix™ confocal technology
- Large field of view captures 1.97 mm² at 10X
- > 3 log dynamic range
- Acquire from a variety of multi-well plate formats up to 1536 including unusual formats developed for 3D assays with round bottom or trans-well plates
- 4 position objective changer and 5 position filter cube changer
- High speed laser and image autofocus
- Simplified Z stack acquisition and analysis
- Widest selection of objectives
 - » 1x to 100x magnification
 - » Air objectives: up to 0.95 NA
 - » Oil objectives: up to 1.4 NA
- Fully automated X-Y and Z stages with resolution better than 100 nm

Learn more about the ImageXpress Micro 4 High-Content Imaging System. Visit www.moleculardevices.com/ixm4 or contact your local sales representative.

Additional enhancements

Available options include field upgradeable Confocal imaging, Environmental Control (temperature, humidity and CO₂ control for multi-day live-cell experiments), Transmitted Light True Phase contrast option and Fluidics (on board disposable pipetting).