



# Molecular Devices Imaging

Elegant workflows, crystalline images, compelling data

# Faster discoveries – More relevant data – Less time to image



## SpectraMax® i3x Multi-Mode Microplate Reader with MiniMax™ imaging cytometer

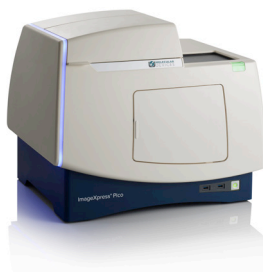
Multi-mode microplate reader  
+ basic imaging

### Features

- SoftMax Pro Software for capture and analysis of plate reader and imaging data
- Customer upgradeable to include reagent injection and western blot detection capabilities

### Key specifications

- 4X microscope objective
- 2 color fluorescence + Brightfield



## ImageXpress® Pico Automated Cell Imaging System

Affordable, compact system with  
intelligent image acquisition and  
analysis capabilities

### Features

- Compact system that can be easily installed on lab benches
- Easy-to-use CellReporterXpress software for image acquisition and analysis
- 20+ pre-built analysis protocols
- Access data from a browser—anytime, anywhere

### Key specifications

- 4X – 63X objective imaging
- White light/brightfield, colorimetric, and fluorescence imaging
- Optional CO<sub>2</sub>, humidity
- Standard temperature control
- Optional dual injectors



## ImageXpress® Nano Automated Imaging System with CellReporterXpress software

The freedom to do more with a  
widefield fluorescence and label-free  
automated imaging to help easily  
automate your cellular or tissue  
based assays.

### Features

- CellReporterXpress software is very easy to use for all lab members
- 20+ pre-built analysis routines
- Access data from a browser – anytime, anywhere
- Tablet friendly

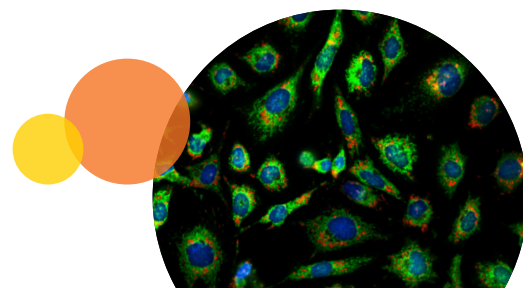
### Key specifications

- 2X–60X microscope objectives available
- UV-VIS fluorescence
- Optional CO<sub>2</sub>, humidity, temperature control
- Optional Brightfield imaging

## See how Mimetas, the organ-on-a-chip company, uses the ImageXpress Micro Confocal system

Mimetas uses their OrganoPlate® to develop disease, toxicology, and transport models for research, development and drug screening. For development and application of tissue models, they use a range of image-based readouts, such as phase-contrast time lapse, fluorescence microscopy, confocal microscopy, in combination with automated 3D image analysis. The team uses the **ImageXpress Micro Confocal High-Content Imaging System** on a daily basis, serving as their standard high-throughput imager for most in-chip assays.

See this story and more at [www.moleculardevices.com/customer-stories](http://www.moleculardevices.com/customer-stories).





## ImageXpress® Micro 4 Widefield High-Content Imaging System

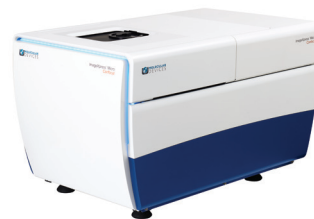
An agile solution that provides stunning widefield and label-free images at high throughput with software that transforms those images into intriguing results.

### Features

- Sophisticated and flexible software that transforms your incredible images into intriguing results
- High-content imaging workhorse that can be configured to fit your assay needs

### Key specifications

- 1X–100X air or oil microscope objectives
- Multi-threaded parallel processing for faster analysis speeds
- Optional on-board liquid handling
- Optional environmental control for CO<sub>2</sub>, humidity and temperature control
- Brightfield and phase imaging options
- Optically-encoded voice coil driven stages with encoders that measure position with better than 25 nm resolution
- Solid state light engine with expected lifetime of >20,000 hours



## ImageXpress® Micro Confocal High-Content Imaging System

Expand the boundaries of your research with the ultimate combination of speed, sensitivity and flexibility in a turnkey solution.

### Features

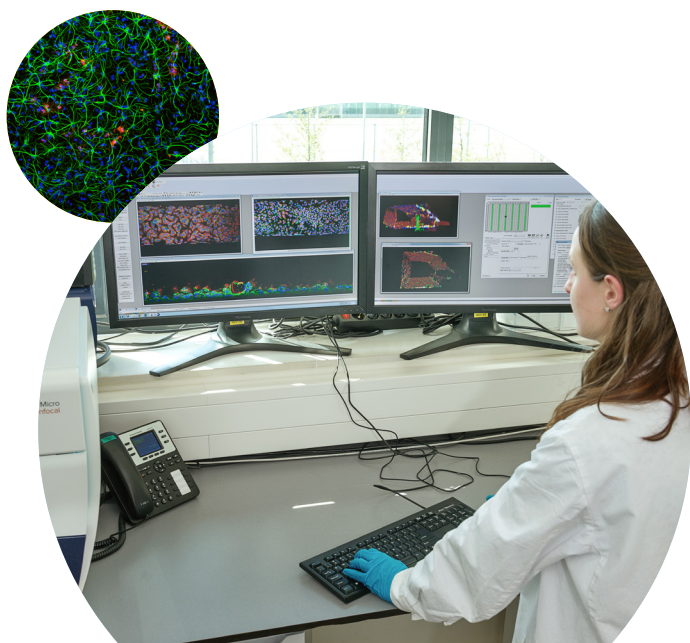
- Sophisticated and flexible software that transforms your incredible images into intriguing results
- High-content imaging workhorse that can be configured to fit your assay needs

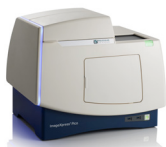
### Key specifications

- 1X–100X objective imaging
- Multiple disc geometries for AgileOptix™ Spinning Disc Technology provides the ability to trade resolution and speed with a selection of disc geometries
- Multi-threaded parallel processing for faster analysis speeds
- Optional on-board liquid handling
- Optional environmental control for CO<sub>2</sub>, humidity and temperature control
- Brightfield and phase imaging options
- Optically-encoded voice coil driven stages with encoders that measure position with better than 25 nm resolution
- Solid state light engine with expected lifetime of >20,000 hours

### Explore modifications with AWES

The Molecular Devices Advanced Workflow Engineering Solution (AWES) team has successfully tailored the ImageXpress Micro system for some customers to include a variety of light engines to address ultra violet (UV) to near infrared (NIR) applications, environmental control with gas mixers for CO<sub>2</sub> and Hypoxia, fluidics with simultaneous imaging and dispensing as well as integration of other lab components such as incubators, liquid handlers, robotics for a fully automated work cell. The AWES team is available to explore these modifications with you. Price, time to deliver, and specifications will vary based on mutually agreed technical requirements. Solution requirements may cause adjustment to standard performance. Purchase Terms available at [www.moleculardevices.com/custom-products-purchase-terms](http://www.moleculardevices.com/custom-products-purchase-terms).





System Specifications	SpectraMax® i3x Multi-Mode Microplate Reader with MiniMax™ imaging cytometer	ImageXpress Pico Automated Cell Imaging System	ImageXpress Nano Automated Imaging System <sup>#</sup>	ImageXpress Micro 4 High-Content Imaging System	ImageXpress Micro Confocal High-Content Imaging System
Tablet and touchscreen friendly		•	•		
Measures cell confluency	•	•	•	•	•
Microplate types	6 to 1536 wells for plate reader, 96 and 384 wells for imaging	Up to 384 wells	Up to 384 wells	Up to 1536 wells	Up to 1536 wells
Microscope slides	•	•	•	•	•
Label-free imaging	•	•	•	•	•
Number of fluorescent channels	2	4	5	5	5
Objective range	4X	4–63X	2–60X	1–100X	1–100X
Timelapse imaging		•	Version 2 of software	•	•
Z stacking				•	•
Optional environmental control for CO <sub>2</sub> , humidity and temperature control		•	•	•	•
Optional automated 3D analysis with volumetric output				•	•
Optional on-board liquid handling				•	•
AgileOptix Spinning Disc Technology				Upgradeable	•
Optional dual injectors	•	•			

<sup>#</sup>Features available with CellReporterXpress software

Contact your Molecular Devices Technical Sales Specialist and get the perfectly right imager for your lab.

#### Contact Us

Phone: +1.800.635.5577  
 Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
 Email: [info@moldev.com](mailto:info@moldev.com)

Check our website for a current listing of worldwide distributors.

#### Regional Offices

USA and Canada  
 +1.800.635.5577  
 United Kingdom  
 +44.118.944.8000  
 Europe\*  
 00800.665.32860

China (Beijing)  
 +86.10.6410.8669  
 China (Shanghai)  
 +86.21.3372.1088  
 Hong Kong  
 +852.2248.6000

Japan (Osaka)  
 +81.6.7174.8331  
 Japan (Tokyo)  
 +81.3.6362.5260  
 South Korea  
 +82.2.3471.9531

\*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland