



## ImageXpress Micro Confocal High-Content Imaging System

The confocal solution for your complex biology

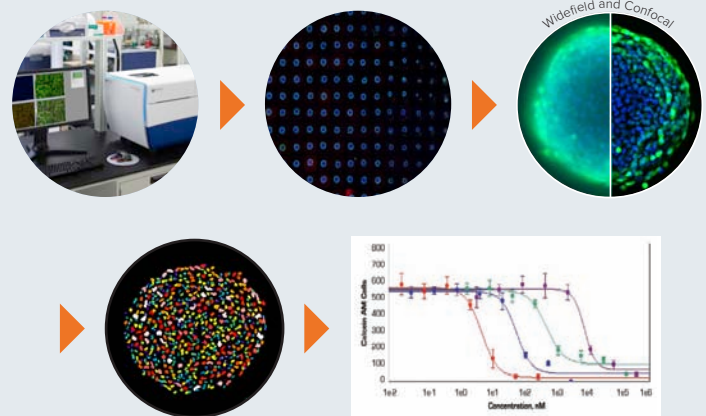
### Higher quality images, faster throughput, and more powerful analysis

The ImageXpress® Micro Confocal High-Content Imaging System provides improved quantification for live or fixed cell assays. This versatile imaging system features a unique confocal technology which allows you to explore more physiologically relevant, complex 3D models including spheroids, tissues, and whole organisms and to generate publication quality images at high throughput for samples in slides or one to 1536-well microplates.

For researchers looking to expand their laboratory's capabilities, the ImageXpress Micro Confocal system leverages large field-of-view optics to map macrostructures with minimal tiling. In addition, querying of large cell populations is accelerated, speeding up the characterization of highly heterogeneous samples or identification of rare subpopulations.

### Key features

- Acquire statistically relevant data quickly with an advanced scientific CMOS detector, enabling >3 log dynamic range
- Improve visualization and quantitation with 3D assay models
- Achieve excellent image quality without sacrificing throughput using our unique optical path technology
- Expand your research capabilities with water immersion objectives, transmitted light, phase contrast optics, on-board liquid handling, and environmental control options



Combined with MetaXpress® High-Content Image Acquisition and Analysis Software, the ImageXpress Micro Confocal system is a complete solution that enables you to interpret your images, understand your data, and explore new ideas—in both widefield and confocal modes.

## Specifications

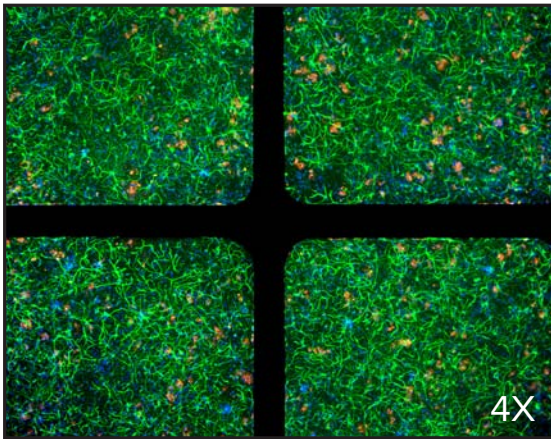
### System

- High-speed laser autofocus with integrated image autofocus option
- Linear encoded voice coil driven X, Y, and Z stages with better than 25 nm resolution
- 4-position automated objective changer\*
- 5-position software selectable dichroic filter wheel\*
- 8-position software selectable emission filter wheel\*
- Sample compatibility: slides and one to 1536-well microplates, round or flat bottom, low to high profile, and Transwell® plates

### AgileOptix optical path

- AgileOptix™ technology enables the ImageXpress Micro Confocal system to deliver the sensitivity and throughput needed for demanding applications by combining a powerful solid-state light engine, high-quantum efficiency 16-bit, >4 megapixel scientific CMOS sensor, and selectable unique confocal geometries
- >3 log dynamic range is available in both widefield and confocal modes
- Large field of view (1.96 mm<sup>2</sup> at 10X) imaging maximizes collection of publication quality images and statistically relevant data

\*User changeable



Acquire greater than 200,000 wells per day at 4X magnification with the system's unique multi-well crop feature.

### 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	China (Beijing)	+86.10.6410.8669	Japan (Osaka)	+81.6.7174.8331
United Kingdom	+44.118.944.8000	China (Shanghai)	+86.21.3372.1088	Japan (Tokyo)	+81.3.6362.5260
Europe*	00800.665.32860	Hong Kong	+852.3971.3530	South Korea	+82.2.3471.9531

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

Option	Feature
Water Immersion Objectives	<ul style="list-style-type: none"> <li>• 20X, 40X, and 60X (up to 1.2 NA)</li> <li>• Increase signal up to 4 times for brighter intensity at lower exposure times</li> <li>• Increase in penetration depths dependent on sample</li> <li>• Improve z-resolution and decrease optical aberrations</li> <li>• Auto water replenishment enables screening or imaging across a plate</li> </ul>
Environmental Control	<ul style="list-style-type: none"> <li>• Multi-day, live cell time-lapse imaging</li> <li>• Provides appropriate atmospheric conditions (e.g. 5% or 10% CO<sub>2</sub>)</li> <li>• Mimics physiological environment (30–40 °C ± 0.5 °C)</li> <li>• Controls humidity and minimizes evaporation (0.5 µL/well/hour for 96- or 384-well formats)</li> </ul>
Phase Contrast	<ul style="list-style-type: none"> <li>• High contrast imaging where unstained cells are easily viewed or separated from background (4X–60X)</li> <li>• Ideal for non-fluorescent histochemically stained samples</li> <li>• Nikon 100W Pillar Diascopic Illuminator with TE-C ELWD Condenser</li> <li>• 0.3 NA with 65 mm WD and PhL, Ph1, and Ph2 selectable phase rings</li> <li>• Fluorophore-independent morphology visualization with fluorescent imaging overlay</li> </ul>
On-board Fluidics	<ul style="list-style-type: none"> <li>• Single-channel pipettor</li> <li>• Dispense volumes from 3 µL to 200 µL (±1 µL; ±5%)</li> <li>• Compatible with 96- or 384-well format FLIPR System pipette tips</li> <li>• Holds two plates for compound addition or media exchange</li> <li>• Optional plate heating</li> <li>• Environmental control</li> </ul>

Note: all options, filters, and objectives are available at point of sale or as after market upgrades. Configurations shown herein do not encompass all configurations available. Contact your sales and support team today to identify the system configuration most suitable for your applications.

## Implement a solution that works for you

Molecular Devices can successfully tailor the ImageXpress Micro Confocal High-Content Imaging System to include customized software and hardware including the features described herein, as well as integration of other lab components such as incubators, liquid handlers, and robotics for a fully automated workflow. With over 30 years of experience in the life science industry, you can count on us to deliver quality products and provide worldwide support.

Sale is subject to our Custom Product Purchase Terms available at [www.moleculardevices.com/custom-products-purchase-terms](http://www.moleculardevices.com/custom-products-purchase-terms).