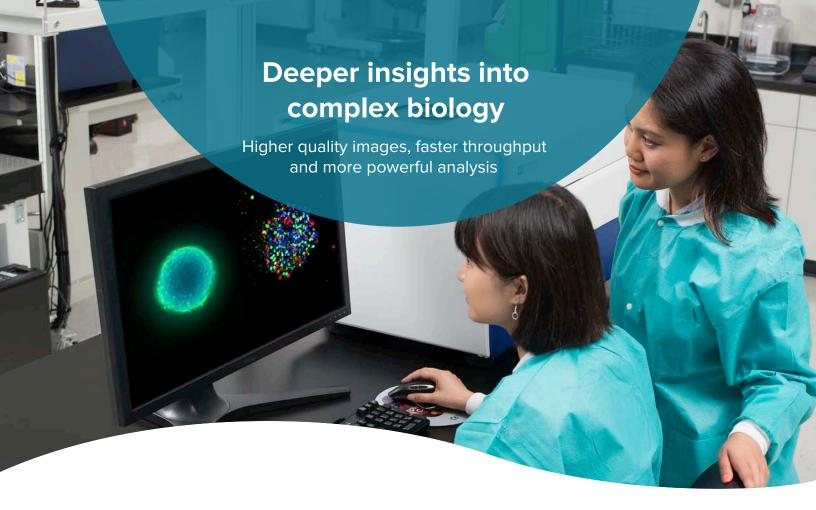


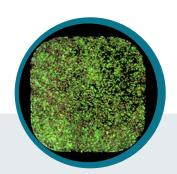
# **ImageXpress Micro Confocal**

High-Content Imaging System

The confocal solution for your complex biology



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 three dimensional 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.



### **Key capabilities**

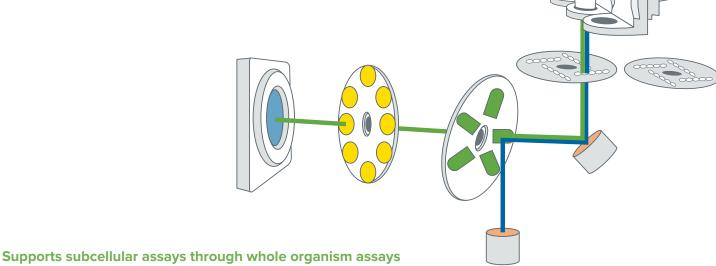
- Select a confocal geometry optimal for your assay and throughput needs
- Achieve excellent image quality without sacrificing throughput via our unique optical path technology
- Acquire statistically relevant data quickly with an advanced scientific CMOS detector, enabling >3 log dynamic range
- · Large field of view enables whole-well imaging
- Expand your research capabilities with water immersion objectives, transmitted light, phase contrast optics, on-board liquid handling, and environmental control options

## Confocal technology at the speed of widefield imaging

- Capture an entire well of a 384-well plate with a single image at 4X magnification
- Capture four wells of a 1536-well plate in a single image at 4X magnification
- Throughput of >160K wells/day confocal, >200K wells/day widefield

### AgileOptix technology at the heart of the ImageXpress Micro Confocal system

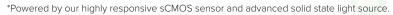
Software-selectable configurations right for your research

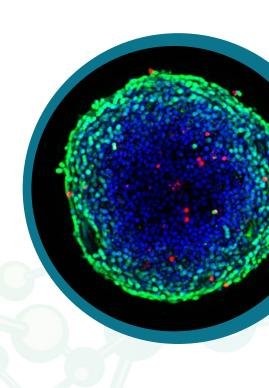


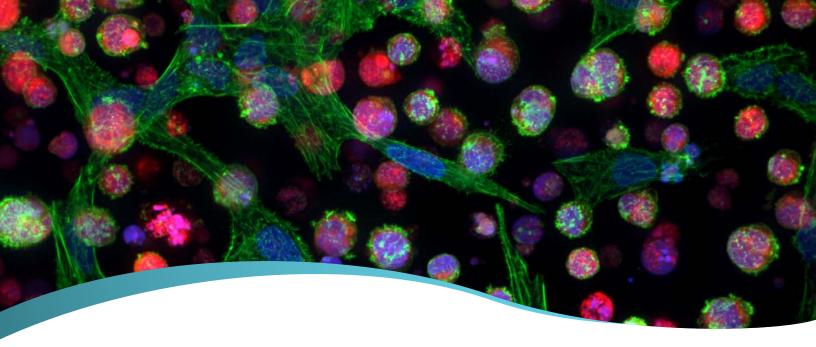
- Widest selection (> 25) of objectives
- Oil objectives with up to 1.4 NA
- Air objectives with 0.05 to 0.95 NA
- Water objectives with up to 1.2 NA

## Select a spinning disk confocal geometry matched to your assay requirements.

Spinning disk geometry	60 μm pinhole (single disk)	60 μm dual disk with 50 μm slit	60 μm dual disk with 42 μm pinhole
High-sensitivity detection	•	•	•
Fast acquisition	•	•	•
>3 log dynamic range*	•	•	•
Widefield mode for flat biology	•	•	•
Most confocal applications	•	•	•
Highest resolution imaging			•
High throughput applications		•	



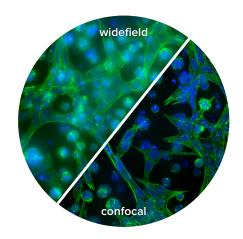




### **Expand your research into a new dimension**

More relevant results with 3D assay models into complex biology

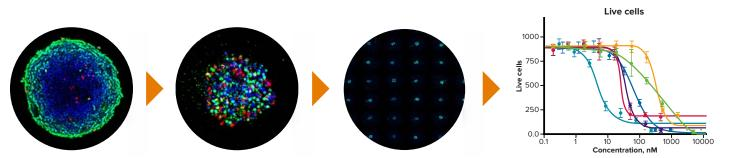
Complex, 3D cellular models yield more predictive, physiologically relevant results versus monocultures or other 2D cellular models. Explore the complexities of these models faster and gain better results even in samples grown in a thick extracellular matrices using the ImageXpress Micro Confocal system. It offers flexible options to meet your specific research needs and ensure the best images for your assay. Get more insight into 3D cells with options like water immersion objectives to improve signal up to 4 times at depth and decreasing optical aberrations for better data quality or monitor cell health kinetics over time using environmental control.



# **Cell stained for nuclei and actin growing in a 3D gel.** Projection image of seven planes acquired with a 40X Plan Apo objective.

### Clearer images and improved quantitative screening for:

- · Spheroids
- Zebrafish and C. elegans
- · Thick tissue samples
- · Homogenous no-wash assays
- Organoids



**Spheroids in round bottom plates.** Dose–dependent effects of selected compounds. Image montage of the 384 plate. Entire spheroid is captured with one image, example segmentation of spheroid and 4-parametric curve fits for the number of live cells in spheroids.

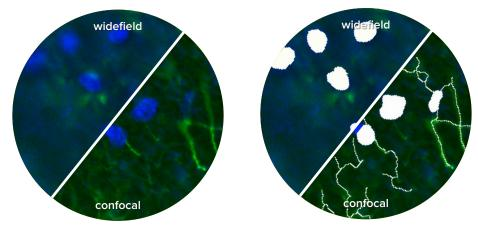
# Improve visualization and quantitation with 3D assay models

### Confocal capability improves image clarity and data quality

Compared with widefield imaging, images of thick samples captured with confocal have reduced background and improved sharpness, resulting in improved image segmentation. Further improve signal and sharpness by using water immersion objectives to capture brighter images at lower exposure times.

### **Enhanced imaging of tissues** and **3D matrices**

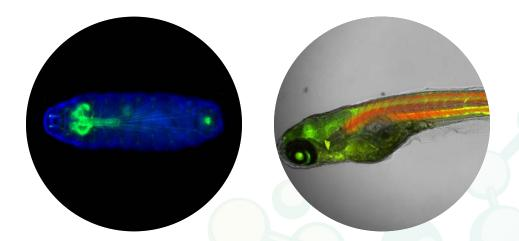
- Select specific cells of interest in 3D matrices such as neurons and stem cells
- Reject high-background fluorescence in thick tissue samples
- Acquire Z-stacks easily with 3D reconstruction capability



Rat brain section, stained for nuclei and neural outgrowths. Images taken with a 20X Plan Fluor ELWD objective, confocal with 60 µm pinhole. Left Without segmentation and right with image segmentation with Neurite Outgrowth module.

## Clearer imaging of whole organisms

- Large field of view enables imaging of an entire well of a 384-well plate with a single image
- Organisms remain in focus for the duration of the experiment
- Perform sophisticated analysis of images and create time-lapse videos



# A complete solution for screening your most complex biological questions

Delivers seamless workflows from image acquisition to data analysis

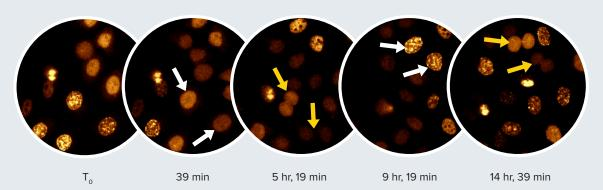
Enjoy the benefits of a streamlined high-content screening (HCS) workflow in a fully integrated environment with our complete imaging solution for your most complex biological questions.

#### **Acquire**

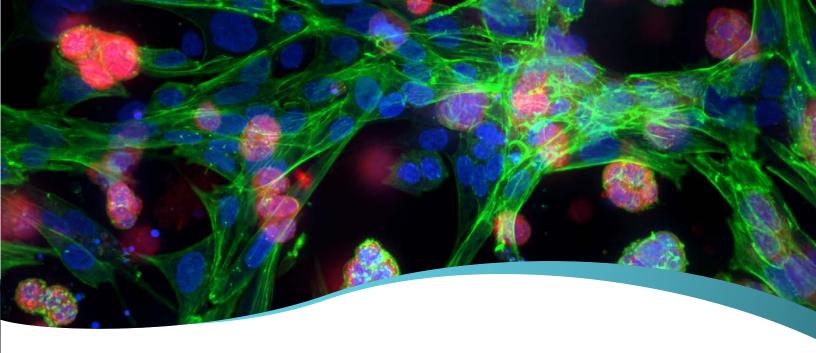
MetaXpress® High-Content Image Acquisition and Analysis Software powers our ImageXpress Micro Confocal system, giving you precise control over image acquisition and analysis, all within a unified interface.

- Acquisition wizard for the entire workflow guides protocol setup for ease of use in multi dimensional imaging
- Laser-based and software configurable image-based auto-focusing system ensures robust focus across a range of sample types
- Acquisition of live cell images enables monitoring of cell growth, death, differentiation, and migration; viral or bacterial invasion, cancer metastasis, chemotaxis, drug toxicity, or translocation





**Live cell acquisition and analysis.** HeLa cells expressing Cell Cycle Chromobody undergoing normal cell division while being imaged in confocal mode. In G1, the cells have a homogeneous fluorescence signal. During S phase, signal accumulates in the nucleus with formation of foci. In G2, the morphology returns to homogeneous and the cell divides. White arrows indicates cell before cell division and yellow arrows indicates daughter cells after division event.



### **Analyze**

Avoid delays in image analysis and data processing using our MetaXpress software with application modules that allow you to quickly and easily analyze your data.

- · Plug-and-play application modules can be adapted to hundreds of image-based analysis workflows
- · Custom module editor empowers you to further tailor your image analysis routines for a perfect fit
- Adaptive Background Correction™ adjusts image segmentation to the local intensity ranges and features within and between cells for better quantitation

Software

- 2D projection algorithms include best focus, maximum and minimum, and sum projection for easy interpretation of 3D image data
- 3D volumetric analysis evaluates volume, XYZ location, distance to neighboring objects, diameter, depth, various intensity measurements, texture, or number of objects
- Save as cell-by-cell and/or image-by-image data

AcuityXpress™ Informatics Software, data visualization, mining, and hit selection are ready to use upon system installation.

#### **Store**

Regardless of the acquisition system used, images taken can be stored in the secure MDCStore™ Data Management Solution.

- Accessible for sophisticated analysis by the MetaXpress® software
- Data migration portal for integration with third-party imaging systems or analysis tools to external host databases or third-party applications



### **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

#### 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 confocal geometries
- · Large field of view (1.96 mm<sup>2</sup> at 10X) imaging maximizes collection of publication quality images and statistically relevant data
- >3 log dynamic range is available in both widefield and confocal modes
- Confocal can be purchased in one of the following three configurations:
  - Single-disk configuration with 60 μm confocal pinhole and widefield modes
  - High-throughput dual disk configuration with 60 μm confocal pinhole, unique and exclusive 50 µm slit confocal and widefield modes
  - High-resolution dual disk configuration with 60 μm and 42 μm confocal pinholes and widefield modes
- · Laser illumination or custom light sources available

Option	Feature		
Water Immersion Objectives	<ul> <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	• Multi-day, live cell time-lapse imaging   • Provides appropriate atmospheric conditions (e.g. 5% or 10% $CO_2$ )   • Mimics physiological environment (30–40 °C ± 0.5 °C)   • Controls humidity and minimizes evaporation (0.5 $\mu$ L/well/hour for 96- or 384-well formats)		
Phase Contrast	High contrast imaging where unstained cells are easily viewed or separated from background (4X–60X)  Ideal for non-fluorescent histochemically stained samples  Nikon 100W Pillar Diascopic Illuminator with TE-C ELWD Condenser  O.3 NA with 65 mm WD and PhL, Ph1, and Ph2 selectable phase rings  Fluorophore-independent morphology visualization with fluorescent imaging overlay		
On-board Fluidics	<ul> <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.

#### **Contact Us** Regional Offices

Phone: +1.800.635.5577 Web: www.moleculardevices.com Email: info@moldev.com Check our website for a current listing of worldwide distributors.

USA and Canada +1.800.635.5577 United Kingdom +44.118.944.8000 00800.665.32860 China (Beijing) +86.10.6410.8669 China (Shanghai) +86.21.3372.1088

+852.3971.3530

Japan (Osaka) +81.6.7174.8331 Japan (Tokyo) +81.3.6362.5260

+82.2.3471.9531 South Korea

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

Hong Kong

Europe\*

<sup>\*</sup>User changeable