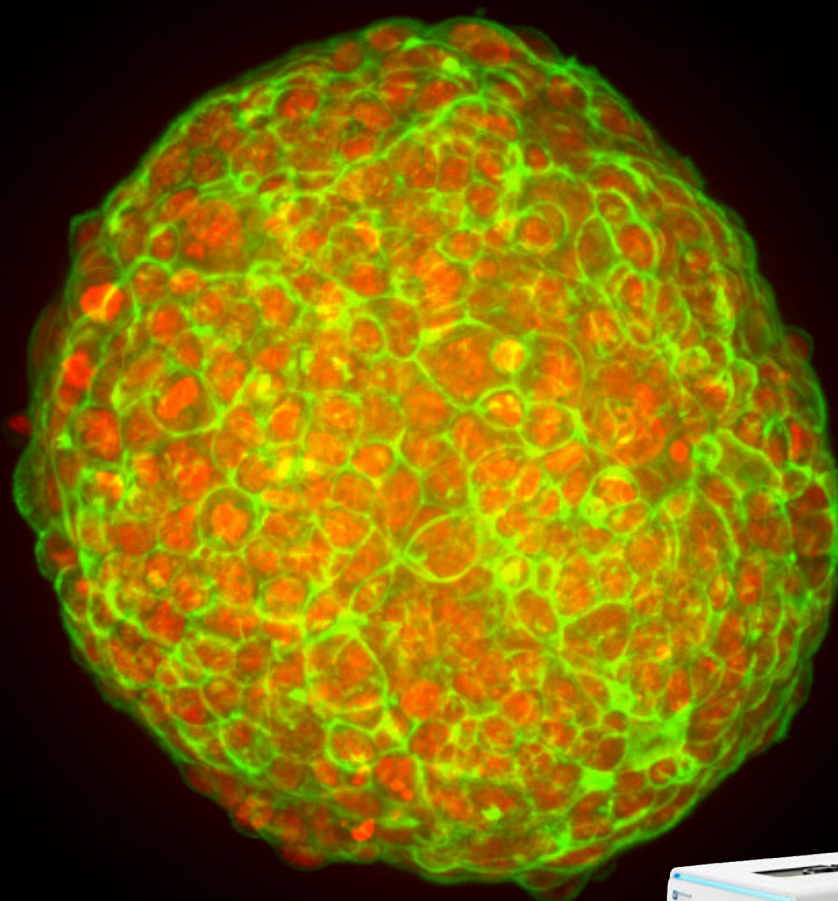


Capture more data at greater depths for 3D and thick tissue samples

Scalable, high-performance, high-content
screening solutions with the ImageXpress®
Micro Confocal system



NEW

High-performance Customizations

- Water immersion objectives
- 5- and 8-channel high-intensity lasers
- Deep tissue penetrating, confocal disk module
- Turnkey high-throughput, long-term kinetics
- Real-time dose response pipettor

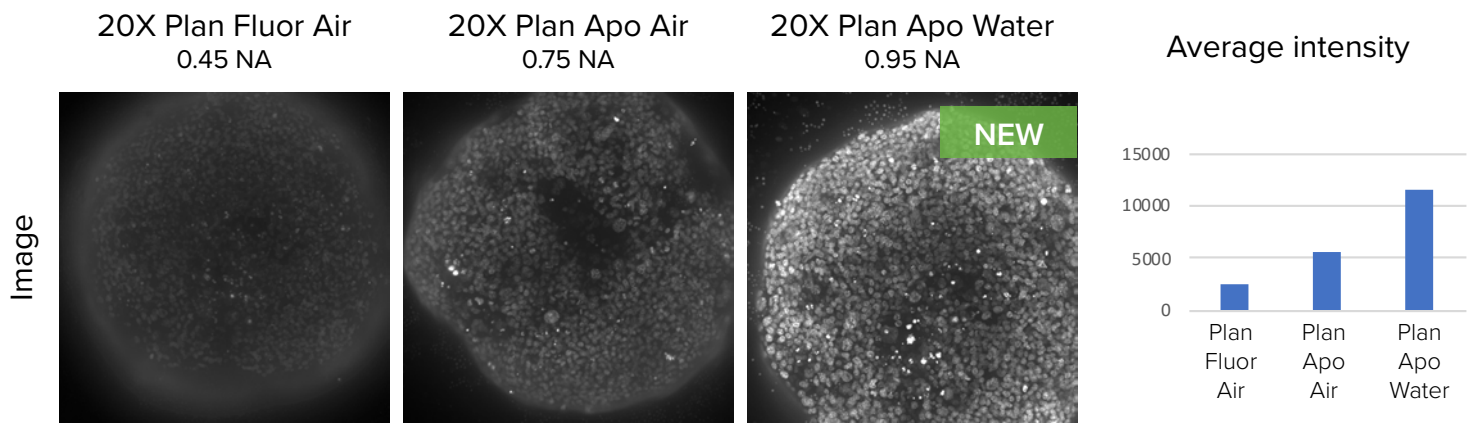




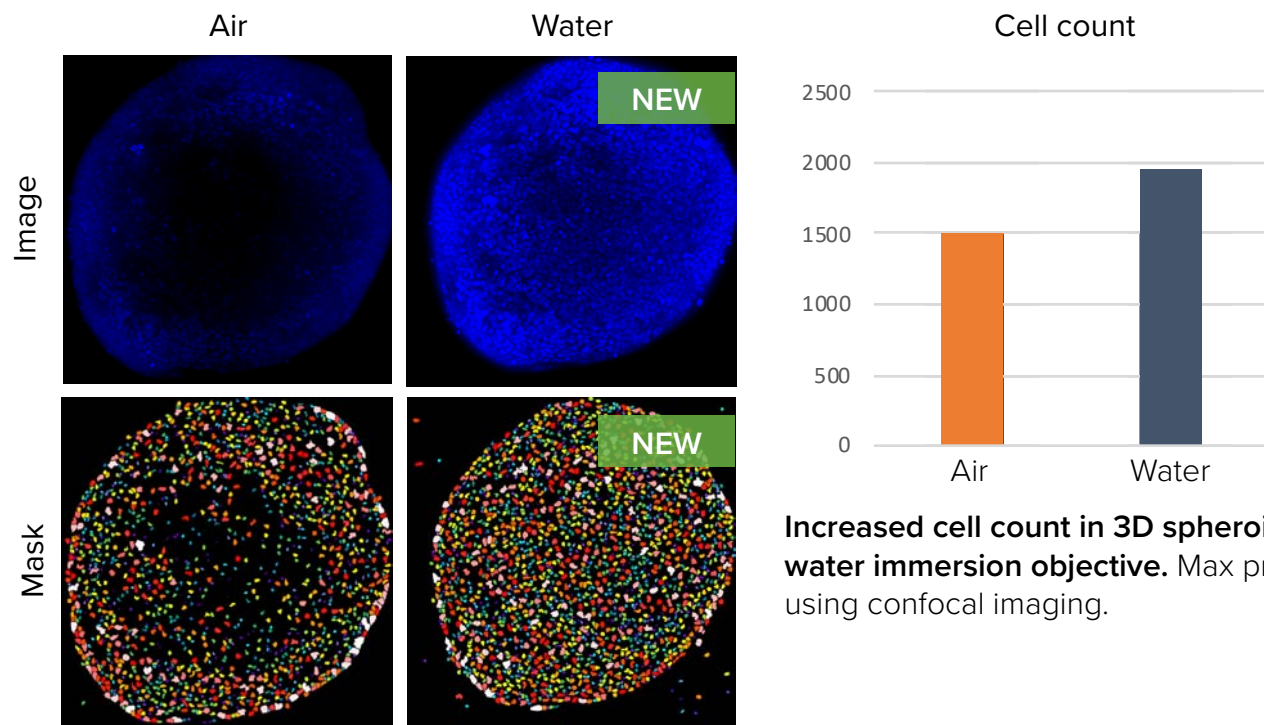
NEW | Water immersion objectives

Water immersion objectives can offer signal increases of up to 4X⁺ which can help scientists see deeper into 3D and thick tissue samples at lower exposure times.

- Increase penetration depths into samples
- Improve z-resolution and spherical aberrations
- Gain sharper, crisper images
- Access more data than ever before



Increased average intensity in 3D assay using water immersion objective. Nuclei of spheroids at 50ms exposure, maximum projection using confocal imaging.



Increased cell count in 3D spheroid using water immersion objective. Max projection using confocal imaging.

*Data and images were acquired during development using customer samples. Results may vary. Highlighted features' price, time to deliver, and specifications will vary based on mutually agreed technical requirements. Solution requirements may cause adjustment to standard performance.

NEW | High-intensity lasers

High output laser excitation can reduce exposure times by up to 75%.[†] The 5-channel laser light source has an output of 400–1,000mw/channel. The 8-channel laser light source (including nIR) with 1,000mw/channel is ideal for customers with increased multiplexing requirements.

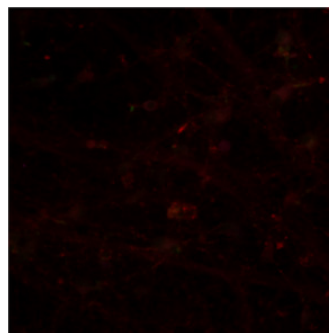
- Obtain sharper images with higher signal-to-noise
- Generate up to a 2X[†] boost in scan speed attributed to significantly reduced exposure times
- Run FRET experiments using lasers for CFP and YFP

NEW | Deep tissue penetrating, confocal disk module

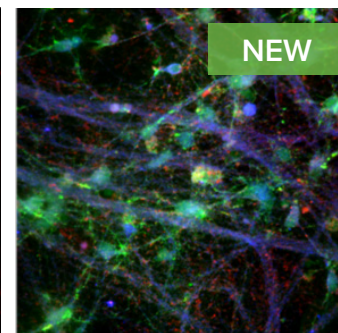
An upgrade to the deep tissue penetrating, confocal disk module, combined with a laser light source, can improve resolution when imaging thick tissue samples[†].

- Improve suppression of out-of-focus light
- Reduce haze (pinhole crosstalk)
- Penetrate deeper into thick tissue samples for sharper images

Standard

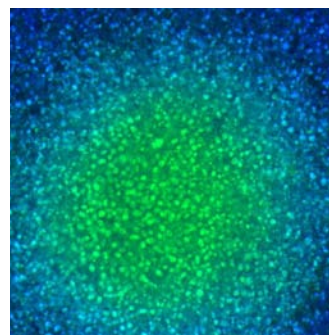


Laser

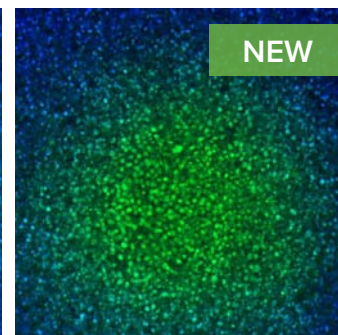


Images taken at the same exposure.

Standard spinning disk



Deep tissue penetrating, confocal disk module

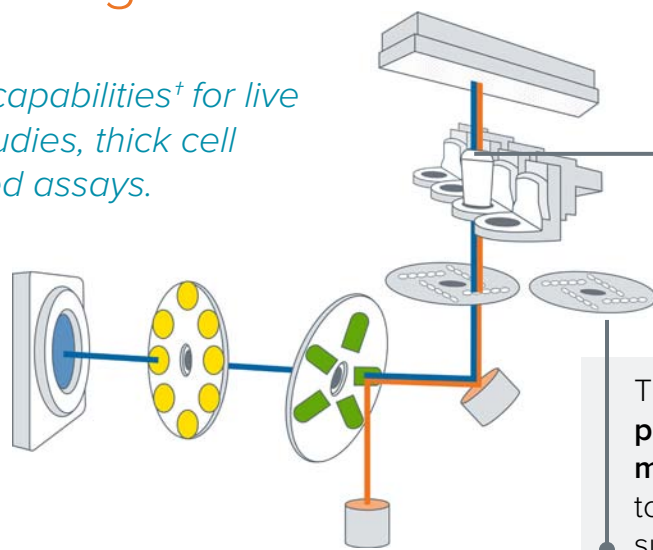


Images taken at the same exposure.

Maximize flexibility in application workflow and experimental design

Get optimized imaging capabilities[†] for live cell imaging, hypoxia studies, thick cell samples, and multiplexed assays.

Expand experiment capabilities with **5- or 8-channel high-intensity lasers.**



20X, 40X, and 60X water immersion objectives improve the geometric accuracy during acquisition and reduce light refraction for brighter intensity at lower exposure times.

The **deep tissue penetrating, confocal disk module** reduces crosstalk to improve out-of-focus light suppression and penetrates deeper into tissue.

[†]Data and images were acquired during development using customer samples. Results may vary. Highlighted features' price, time to deliver, and specifications will vary based on mutually agreed technical requirements. Solution requirements may cause adjustment to standard performance.



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 work-cell. 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.

Additional high-performance customizations



Scale-up automation

Increase throughput, eliminate human errors, maintain sterility, and achieve consistent sample handling. Modular automation design—components can be added in modules and are upgradeable.



Customized optics and light sources

Improve image quality, system sensitivity, confocality, and ability to image at depth using water immersion objectives. Improve speed and image quality with multiple laser upgrade options.



Real-time dose response

Automatic pipettor enables compound addition while simultaneously live streaming at >100 frames per second.



Turnkey high-throughput long-term kinetics

Schedule and image multiple plates over long periods of time while keeping consistent temperature, O₂ (Hypoxia), CO₂, and humidity conditions. Expand live cell walk-away capacity to 200+ plates.

**Add media exchange for a turnkey high throughput organoid screening solution.

Contact Us

Phone: [+1.800.635.5577](tel:+18006355577)
Web: www.moleculardevices.com
Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada
[+1.800.635.5577](tel:+18006355577)
United Kingdom
[+44.118.944.8000](tel:+441189448000)
Europe*
[00800.665.32860](tel:0080066532860)

China (Beijing)
[+86.10.6410.8669](tel:+861064108669)
China (Shanghai)
[+86.21.3372.1088](tel:+862133721088)
Hong Kong
[+852.2248.6000](tel:+85222486000)

Japan (Osaka)
[+81.6.7174.8331](tel:+81671748331)
Japan (Tokyo)
[+81.3.6362.5260](tel:+81363625260)
South Korea
[+82.2.3471.9531](tel:+82234719531)

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