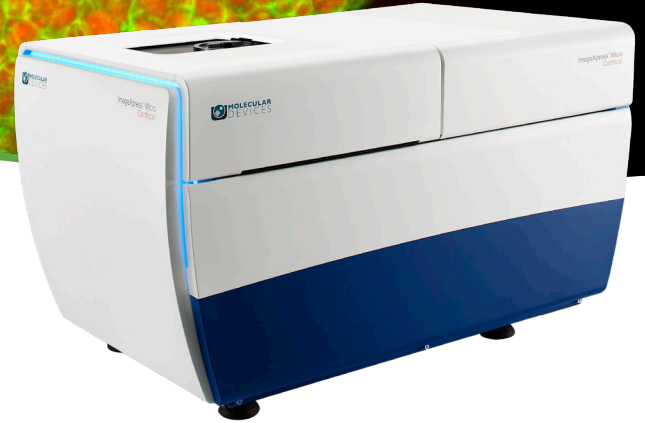
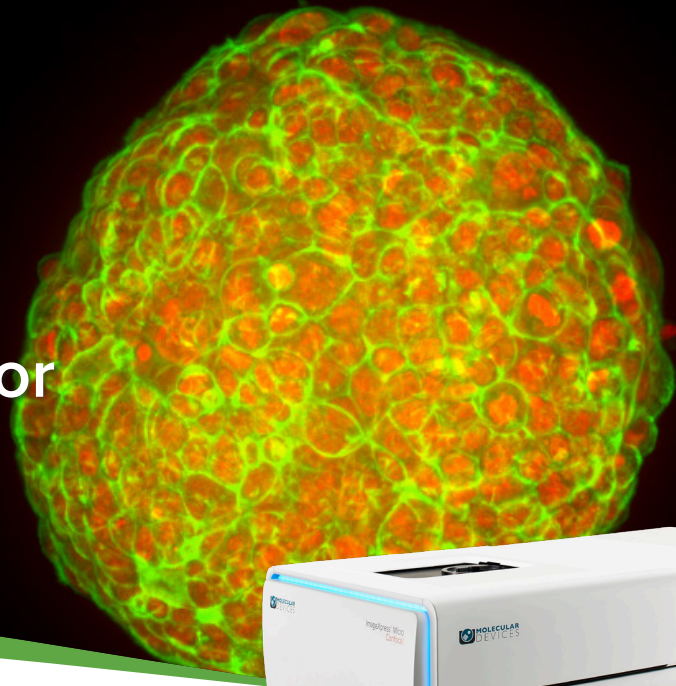


# Capture more data at greater depths for large 3D samples

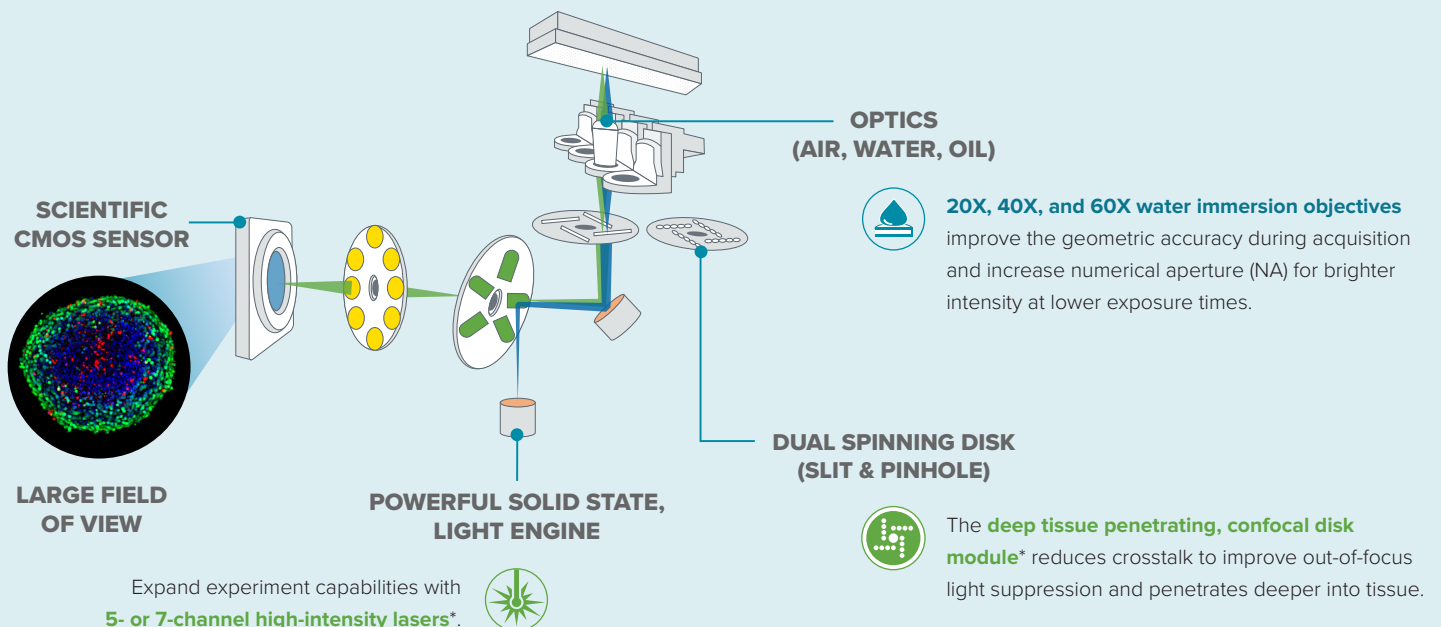


## Scalable, high-performance, high-content screening solutions with the ImageXpress® Micro Confocal system

- Water Immersion objectives
- 5- and 7-channel high-intensity lasers
- Deep tissue penetrating, confocal disk module
- Turnkey high-throughput, long-term kinetics
- Real-time dose response pipettor

## Maximize flexibility in application workflow and experimental design

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

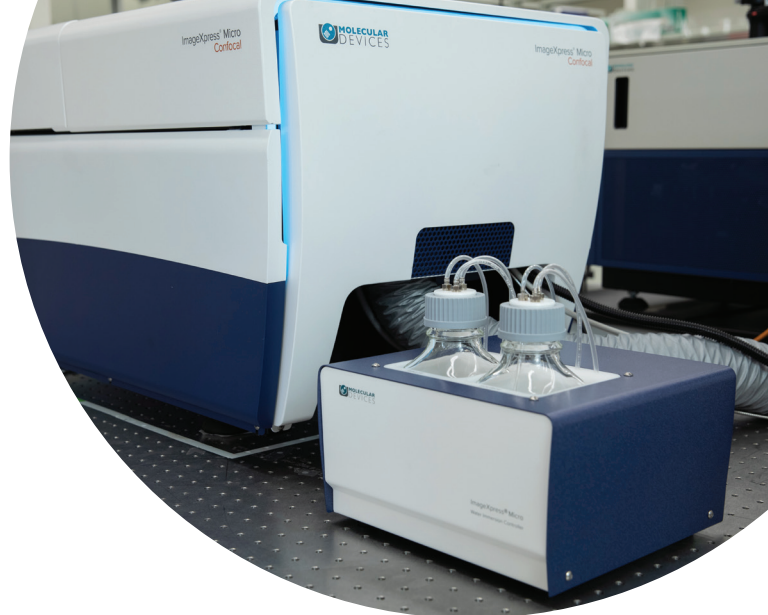




## Water immersion objectives

Water immersion objectives can offer signal increases of up to 4X<sup>+</sup> 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

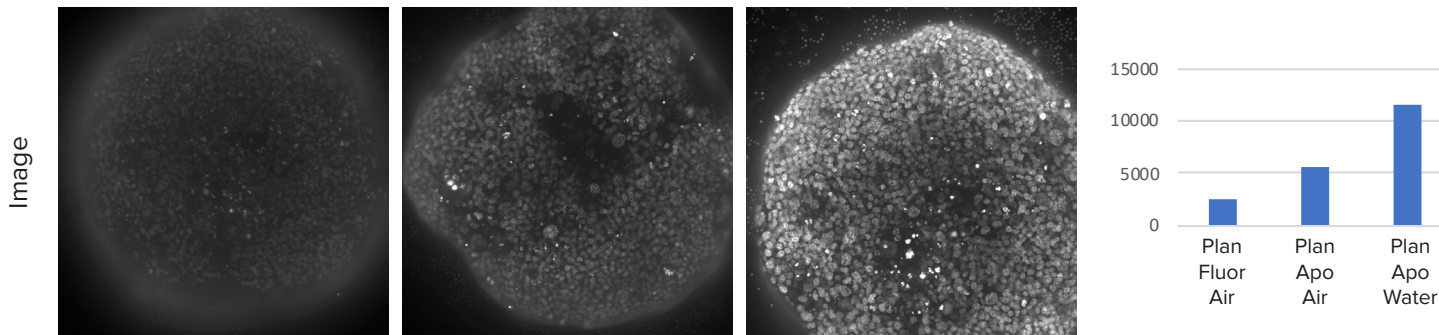


20X Plan Fluor Air  
0.45 NA

20X Plan Apo Air  
0.75 NA

20X Plan Apo Water  
0.95 NA

Average intensity

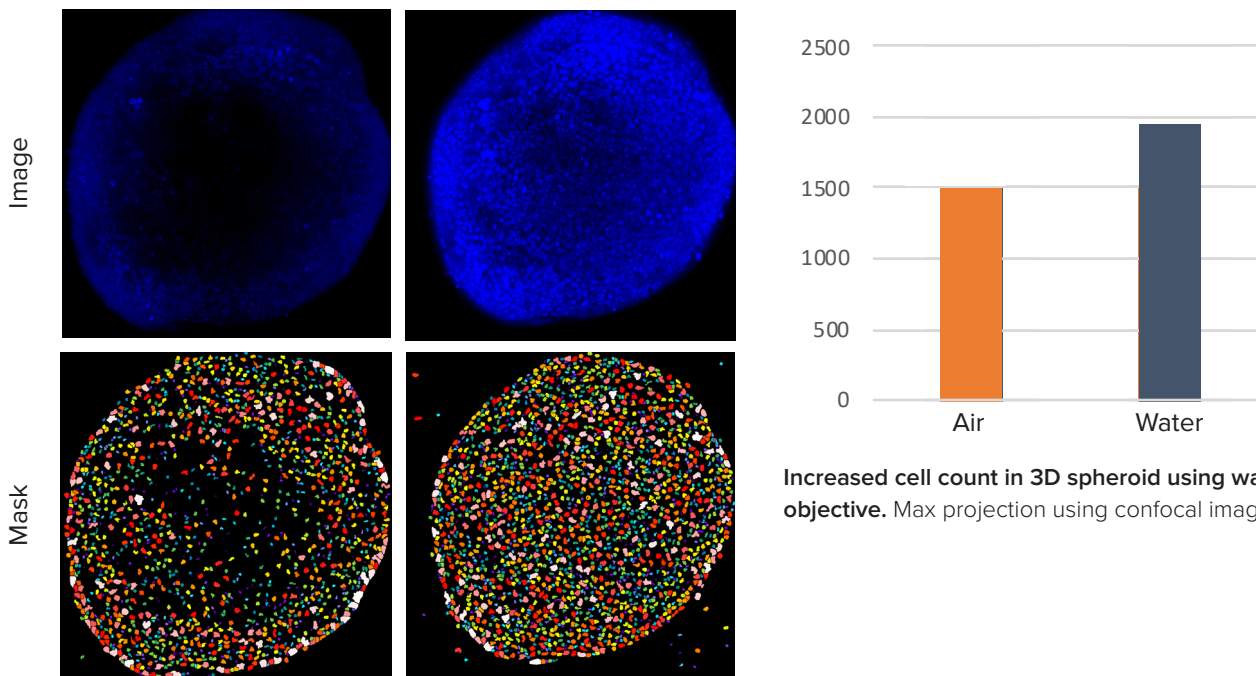


Increased average intensity in 3D assay using water immersion objective. Nuclei of spheroids at 50ms exposure, maximum projection using confocal imaging.

Air

Water

Cell count



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.

## \*High-performance customizations

Molecular Devices can successfully tailor the ImageXpress Micro Confocal High-Content Imaging System to include customized software and hardware including the features described below, 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](http://www.moleculardevices.com/custom-products-purchase-terms).

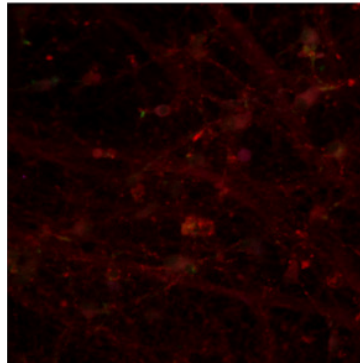


### High-intensity lasers

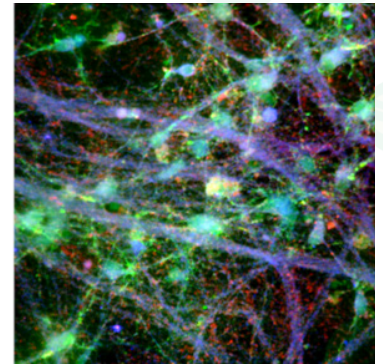
High output laser excitation can reduce exposure times by up to 75%.<sup>†</sup> The laser light source is available as either a 5- or 7-channel light source with outputs of 400–1,000 mW/channel. The 7-channel laser light source includes near IR and is ideal for customers with increased multiplexing requirements.

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

Standard



Laser



Images taken at the same exposure.

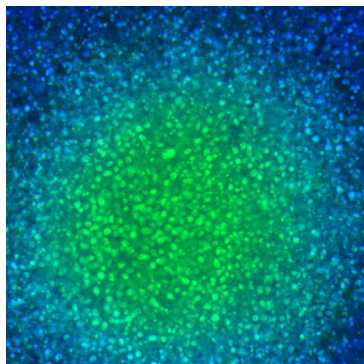


### Deep tissue penetrating, confocal disk module

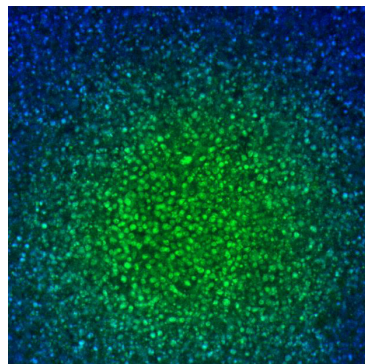
An upgrade to the deep tissue penetrating, confocal disk module, combined with a laser light source, can improve resolution when imaging thick tissue samples<sup>†</sup>.

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

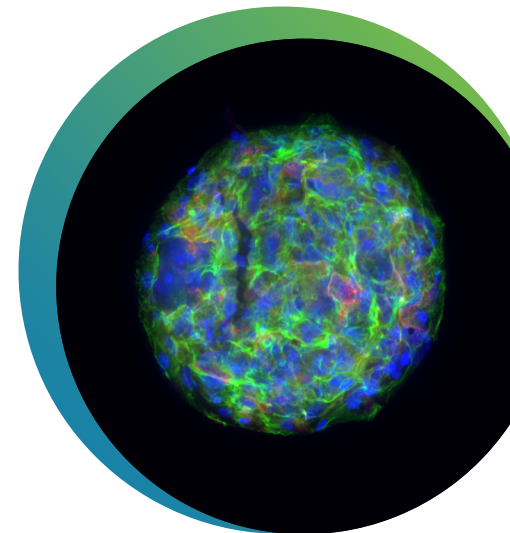
Standard spinning disk



Deep tissue penetrating, confocal disk module

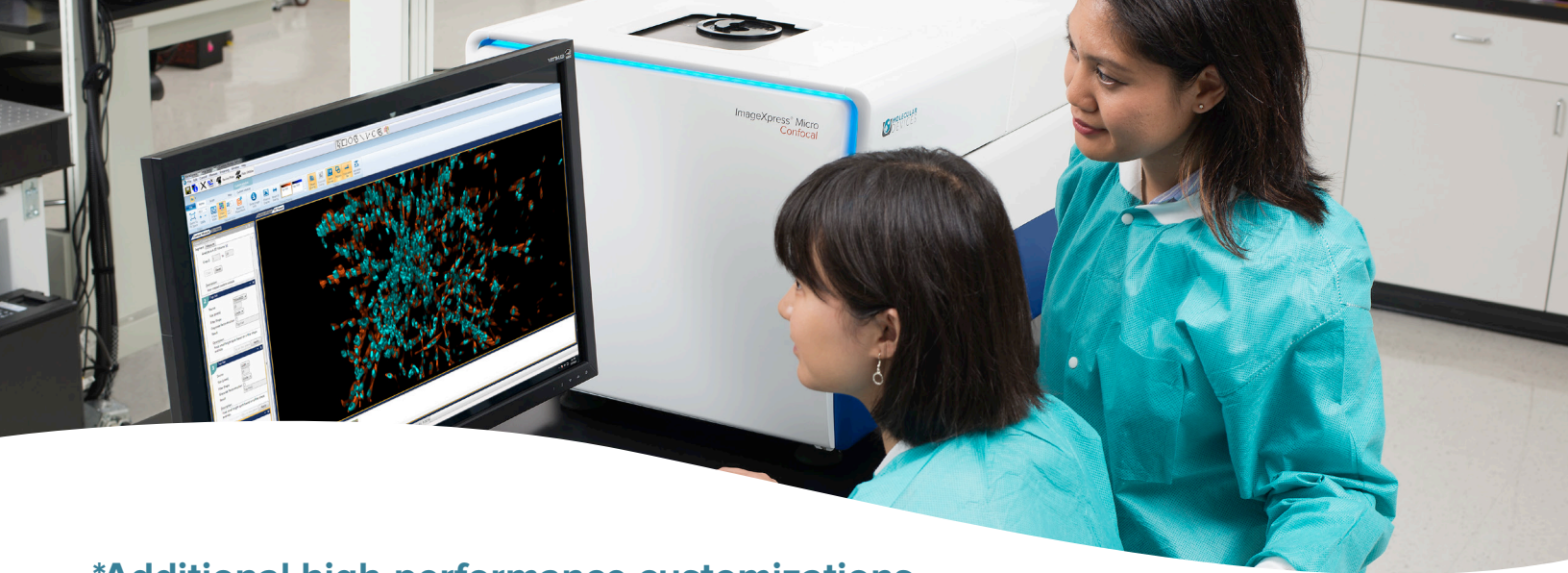


Images taken at the same exposure.



<sup>†</sup>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.





## \*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.



### Real-time dose response

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



### Customized light sources

Improve speed and image quality with multiple laser upgrade options.



### Turnkey high-throughput long-term kinetics

Schedule and image multiple plates over long periods of time while keeping consistent temperature, O<sub>2</sub> (Hypoxia), CO<sub>2</sub>, and humidity conditions. Expand live cell walk-away capacity to 200+ plates.

## 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 above, 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](http://www.moleculardevices.com/custom-products-purchase-terms).

### 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	+81.3.6362.9109
United Kingdom	+44.118.944.8000	China (Shanghai)	+86.21.3372.1088	South Korea	+82.2.3471.9531
Europe*	00800.665.32860	Hong Kong	+852.3971.3530		

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