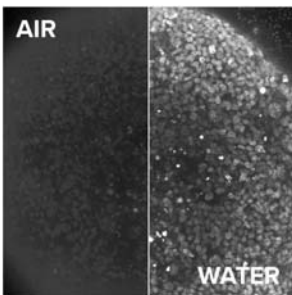


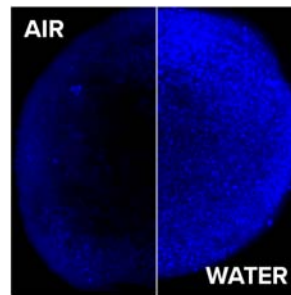
## Water Immersion Objectives for the ImageXpress Micro Confocal system

Gain the sensitivity to capture more phenotypic data at greater depths. Our high-content imagers and expert team are here to help your lab simplify imaging and analysis of complex, 3D assay workflows and fast-track your discoveries.



### Gain up to 4 times more signal at depth

Improve z-resolution and decrease optical aberrations for more accurate reconstruction of 3D samples.

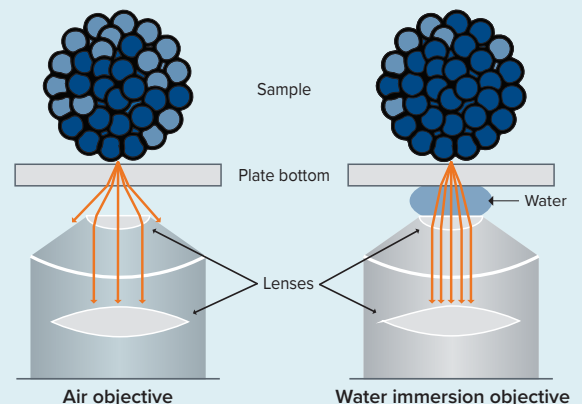


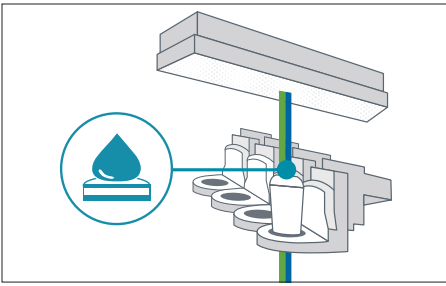
### Gain sharper, crisper images

Better light collection results in images with brighter intensity at lower exposure times, contributing to better quality of data.

## Benefits

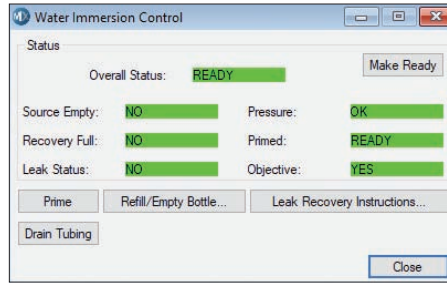
- Increase in penetration depths dependent on sample
- Improve z-resolution and decrease optical aberrations
- Gain sharper, crisper images
- Acquire 3D samples at high throughput
- Shorten image acquisition time





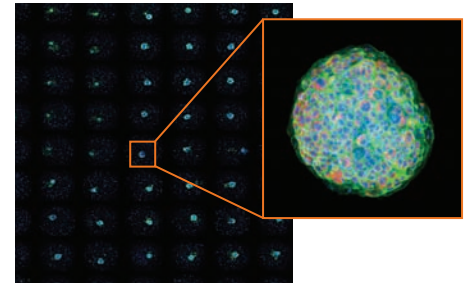
### Streamline set up to get started more quickly

Water immersion objectives are field upgradeable and can be implemented with minimal change to current workflows.



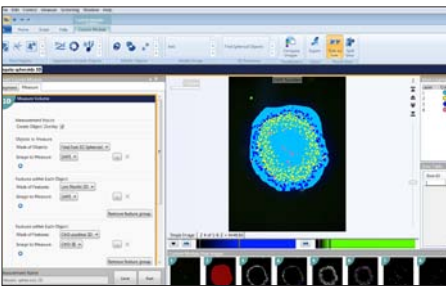
### Run your samples with confidence

MetaXpress software includes integrated, user-intuitive sensors and alerts to indicate the system health of the water controller system to give you peace of mind when running your samples.



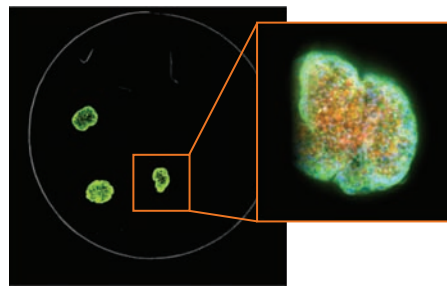
### Acquire 3D samples at high throughput

The ImageXpress Micro Confocal system with water immersion objectives automatically replenishes the water during acquisition, so the entire microplate sample can be acquired at high throughput.



### Simplify viewing and quantitation of 3D structures

Analyze spheroids, microtissues, cells in a 3D matrix, and small organisms that are acquired as a stack of z-planes. Utilize the MetaXpress 3D Analysis Module to evaluate volume, XYZ location, distance to neighboring objects, diameter, depth, various intensity measurements, texture, or number of objects.



### Shorten image acquisition time

QuickID targeted acquisition images at low magnification to find objects of interest or rare events, then, automatically acquires high-magnification images with the required wavelengths, z-stack, or time points. This analysis process shortens acquisition time and reduces imaging storage requirements by acquiring selected targets at high magnification.



Specifications	
Available water objectives	20X, 40X, and 60X (up to 1.2 NA)
Signal (Brightness/intensity)	Increase up to 4 times
Auto-water replenishment	Yes
Automation-friendly	Yes
Additional options	Transmitted Light Tower, Environmental Control

Available high-performance customizations <sup>†</sup>	Laser Light Source, Deep Tissue Penetrating Confocal Spinning Disk Module
Field upgradeable	Yes
Compatible systems	ImageXpress <sup>®</sup> Micro Confocal High-Content Imaging System

<sup>†</sup>For terms and conditions, please visit [moleculardevices.com/ixmc#options](http://moleculardevices.com/ixmc#options)

#### 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 +852.3971.3530

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