



# MetaXpress 6 Software

Transform your high-content imaging workflow

## BENEFITS

- **Integrated Z-stack acquisition and analysis**
- **Faster-than-acquisition 2D Z analysis enabled by MetaXpress PowerCore software**
- **3D analysis module simplifies 3D data quantification and throughput**
- **New peak analysis for cardiac beating assays and curve fitting**

## Simplify workflow

Reimagine your current workflow with MetaXpress® High-Content Image Acquisition and Analysis Software. The latest layout simplifies plate acquisition setup by consolidating key functions in one place. Always visible and always accessible plate layout views make acquiring your data easy and quick including the capability of appending time points to already acquired plates for asynchronous kinetic experiments.

## Intuitively set up 3D acquisition

Easy-to-use graphical interface enables you to rapidly set up 3D imaging. Interactive tools help you determine optimal settings for Z acquisition. Save images from each and every plane or use best focus, maximum, minimum, or sum software algorithms to generate and save only a single 2D projection image.

## Customize analysis with ease

Extend your assay capabilities as your needs grow and change using Custom Module Editor. Quickly render and record 3D visualization of raw data and segmentation as well as quantify your 3D structures with measurements such as volume, intensity, distance and more.

## Strong foundation, great new features

Key features	Versions 3 & 4	Version 5	Versions 6 & 6.5
<b>MetaXpress® Acquisition and Offline Analysis Software</b>			
Time-series acquisition and analysis	With Journals	Integrated with Acquisition and Analysis Wizards	Integrated with Acquisition and Analysis Wizards
Z-series acquisition	With Journals	With Journals	Integrated with Acquisition Wizard
Z-series analysis	With Journals	With Journals	<ul style="list-style-type: none"> <li>• Integrated into workflow</li> <li>• Compatible with MetaXpress PowerCore software</li> </ul>
Customized analysis	With Journals	External Custom Module Editor	Integrated Custom Module Editor
3D analysis	N/A	N/A	Available within Custom Module Editor
Transmitted light segmentation	With Journals	With Journals	Simple step cards available with Custom Module Editor (6.5)
Asynchronous kinetics	N/A	N/A	Integrated with Acquisition Wizard (6.5)
Basic curve fitting	With AcuityXpress software	With AcuityXpress software	Available with MetaXpress software (6.5)
Peak analysis	N/A	N/A	Available with MetaXpress software (6.5)
Windows compatibility	Windows XP	Windows 7 acquisition and analysis	<ul style="list-style-type: none"> <li>• Windows 7 acquisition and analysis</li> <li>• Windows 8 analysis</li> <li>• Windows 10 acquisition and analysis (6.5 and above)*</li> </ul>
<b>MetaXpress® PowerCore™ Software</b>			
Scale analysis to be faster than acquisition	Single time point and plane	Including time-lapse	Including time-lapse, multiple z planes, and custom modules
Run multi-core analysis on acquisition computer	No	No	Yes

\*Windows 10 acquisition for ImageXpress Micro XLS, ImageXpress Micro 4 and ImageXpress Confocal Systems. Windows 10 acquisition computer must be provided by Molecular Devices LLC.

## Application modules

We address hundreds of image-based analysis scenarios. Our turnkey application modules tackle your basic tasks. With MetaXpress Custom Module Editor, you are empowered to tailor image analysis to your research needs.

Application modules	Part number	Application modules	Part number
Custom Module Editor	9500-0158	Mitotic index	9500-0036
3D analysis (requires Custom Module Editor)	5047663	Micronuclei	9500-0045
Curve fitting	5058772	Monopole detection	9500-0039
Transmitted light segmentation (requires Custom Module Editors)	5058774	Multi-wavelength cell scoring	9500-0038
Angiogenesis tube formation	9500-0016	Multi-wavelength translocation	9500-0044
Cell cycle	9500-0040	Neurite outgrowth	9500-0015
Cell health	9500-0035	Transfluor® assay	9500-0033
Cell scoring	9500-0037	Translocation (includes Translocation-Enhanced)	9500-0014
Count nuclei	9500-0017	Transfluor HT assay	9500-0043
Granularity	9500-0032	Nuclear translocation HT	9500-0041
Live/dead	9500-0034	Cell proliferation HT	9500-0042
		All MetaXpress software application modules	5062036

### Contact Us

Phone: [+1.800.635.5577](tel:+18006355577)  
 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.