

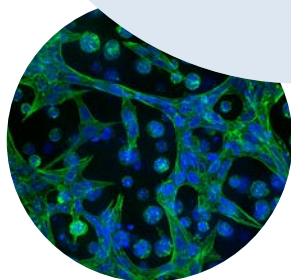


MetaXpress High-Content Image Acquisition and Analysis Software



BENEFITS

- Meet high throughput requirements with a scalable, streamlined workflow
- Adapt your analysis tools to tackle your toughest problems, including 3D analysis
- Run routine modules and custom analyses on parallel processors
- Schedule automatic data transfer between third-party hardware sources and secure database



Accelerate your discoveries

Quickly identify and quantify heterogeneous, sub-cellular and multi-parametric responses in diverse biological systems

Capture high-resolution images, recognize and segment objects, calculate parameters, and then translate your information into meaningful data. At Molecular Devices, we're inspired by your creativity to solve research challenges. Our goal is to provide you with the tools that let you build a successful High-Content Screening (HCS) infrastructure. Our MetaXpress® High-Content Image Acquisition and Analysis Software is the cornerstone of this infrastructure. It is built on our industry-leading software that has been providing scientists with insights into the world of cellular function for over 30 years. We took our extensive experience with cellular imaging and analysis techniques and put it all in one package, delivering an HCS solution as unique as your research.

An end-to-end solution for screening your most complex biological questions

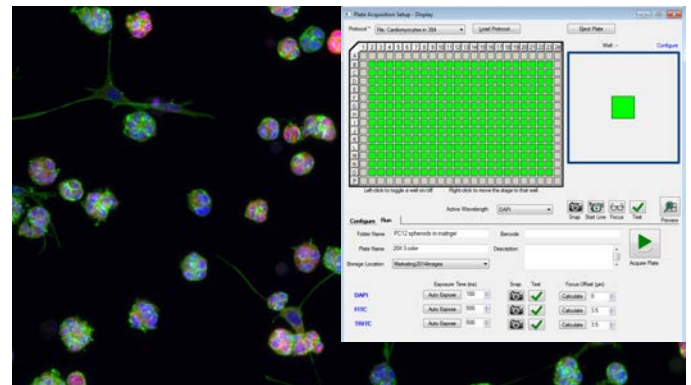
Creating an intuitive HCS workflow requires many components working in concert. Our MetaXpress software ties together every step seamlessly, from acquisition to data management.



Acquire

MetaXpress software offers the precise control you need to acquire images with the best possible quality.

- Capture crisp, high-resolution images using a range of objectives from 1x to 100x at various numerical apertures
- Maintain robust focus across a range of sample types with our easily configurable laser-based or image-based auto-focusing system
- Easily integrate live cell acquisition and quantify 3D measurements of thick samples or kinetic assays using the acquisition and analysis wizard
- Halve your exposure time, or increase assay quality by 30% using the digital confocal option in the acquisition interface



Analyze

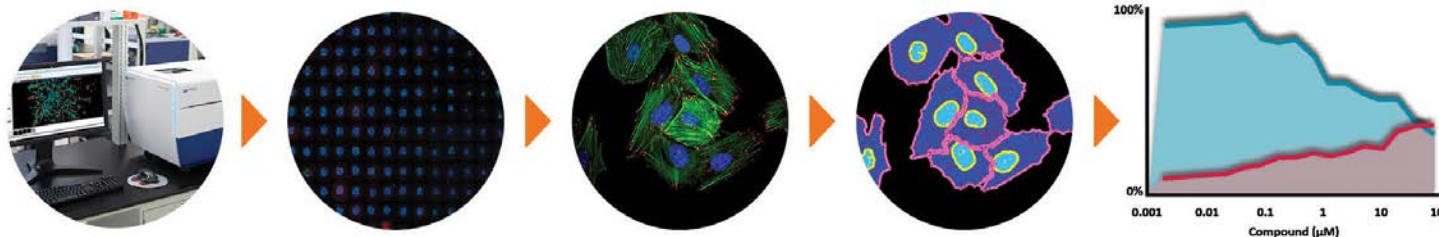
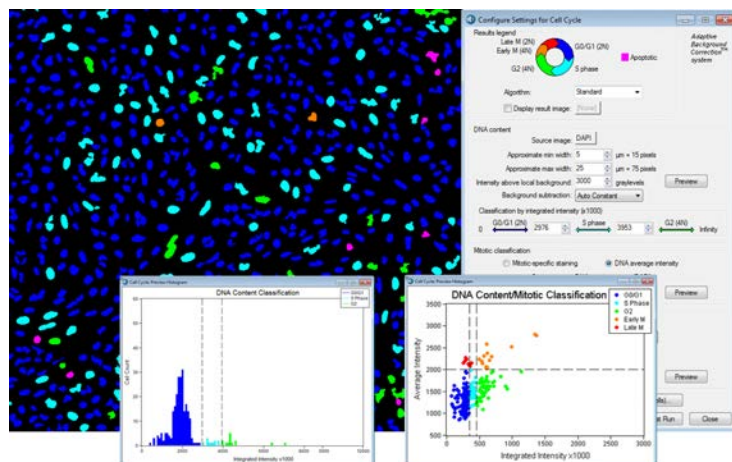
Expedite image analysis and data processing using our MetaXpress software with application modules that allow you to quickly and easily analyze your data.

Modular toolbox

Set up hundreds of routinely used HCS assays whether you're a novice or an expert. MetaXpress software application modules give you an interactive environment to quantify 2D and 3D measurements, including size, intensities, and distance.

Intelligent segmentation

All MetaXpress software application modules are equipped with a proprietary adaptive background correction algorithm, which adjusts image segmentation to the local intensity ranges and shape features within and between cells. This capability gives you the most robust segmentation available in an image-based screening system and improves detection on both phase contrast and fluorescently labeled images. Save analysis results at the cell-by-cell and image-by-image levels with segmentation overlays and review them interactively by selecting a cell and visualizing the corresponding analysis data or vice versa.



Store

MDCStore Data Management Solution and MDCStore Xchange Service provide a seamless solution for storing, organizing, and sharing your valuable data. No extra steps are required to transfer data to MDCStore after acquisition, so you can re-analyze your data without accessing the acquisition system, maximizing the availability of the instrument for further imaging. Save experiment protocols and share them with any ImageXpress or MetaXpress system connected to the MDCStore Data Management Solution to standardize routines in your laboratory or across your organization.

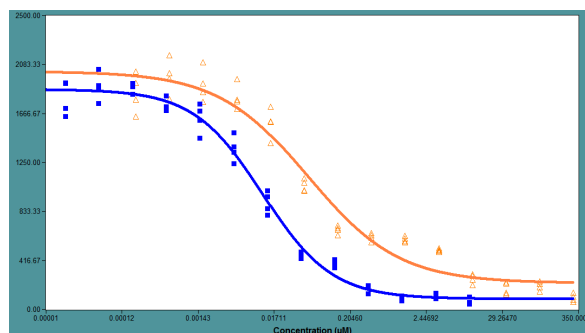
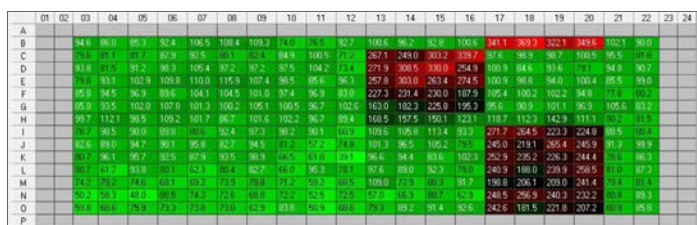
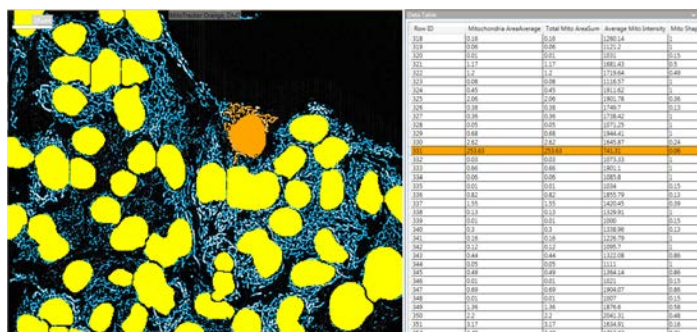
Schedule automatic data transfer between third-party hardware sources and your secure database, and export data to 3rd party and OME compatible files. MDCStore Xchange Service can also be set up to automatically transfer images across two designated locations, making it easy to share data in your file format of choice.

Visualize

Review your results with ease, minimizing the need to switch between applications by simultaneously viewing image thumbnails and plate heatmaps. Analyze, interpret, and mine data further with AcuityXpress™ High-Content Informatics software.

Key features include:

- Integration into MDCStore Data Management tool for easy access to needed files
- Bi-directional image viewer that facilitates data quality control by connecting images to numerical data
- Multi-parametric analysis and visualization
- Dose response analysis and curve fitting with over 30 different predefined functions



Customize

Our solution is scalable and adaptable to fit your ever changing and unique research needs.

Custom module editor

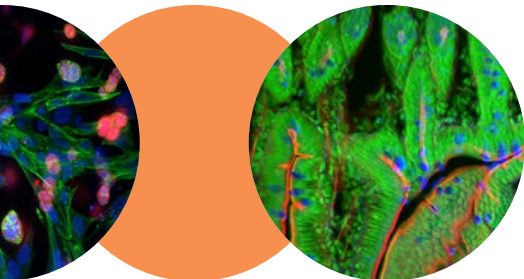
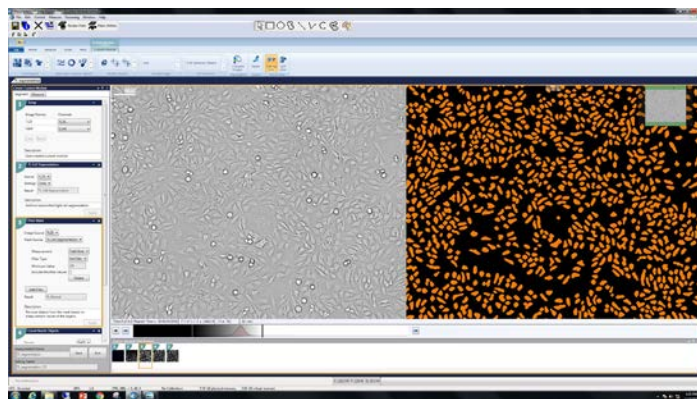
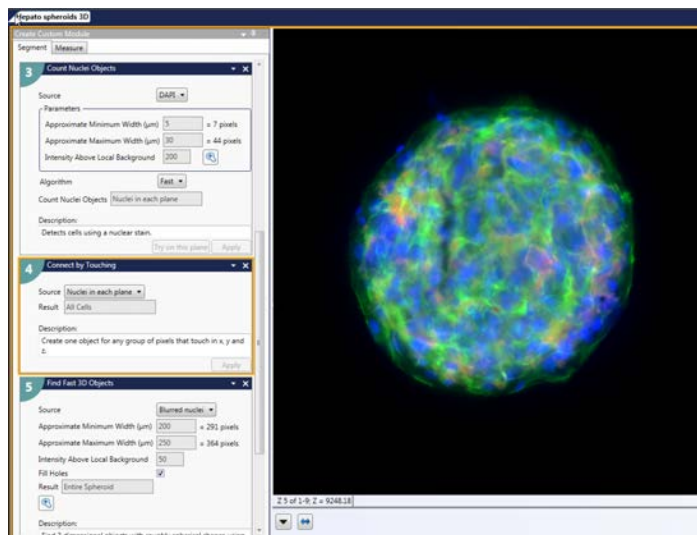
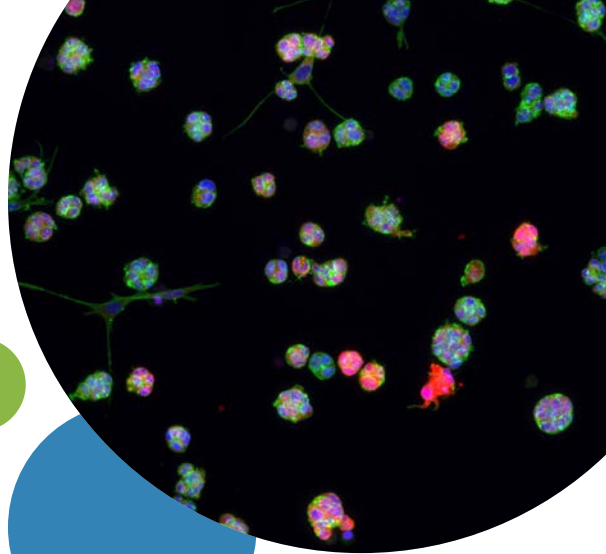
Build specific analysis capability by combining a series of subroutines. With our step-by-step, interactive editor, you can create challenging, multi-step algorithms such as counting punctae in neurite outgrowths, and then running them in minutes using MetaXpress® PowerCore™ software.

3D analysis

Increase the biological relevance of your screening assays with an integrated and seamless acquisition and analysis of 3D data. Simplify quantification of 3D structures with volume, intensity, and distance measurements.

Journals

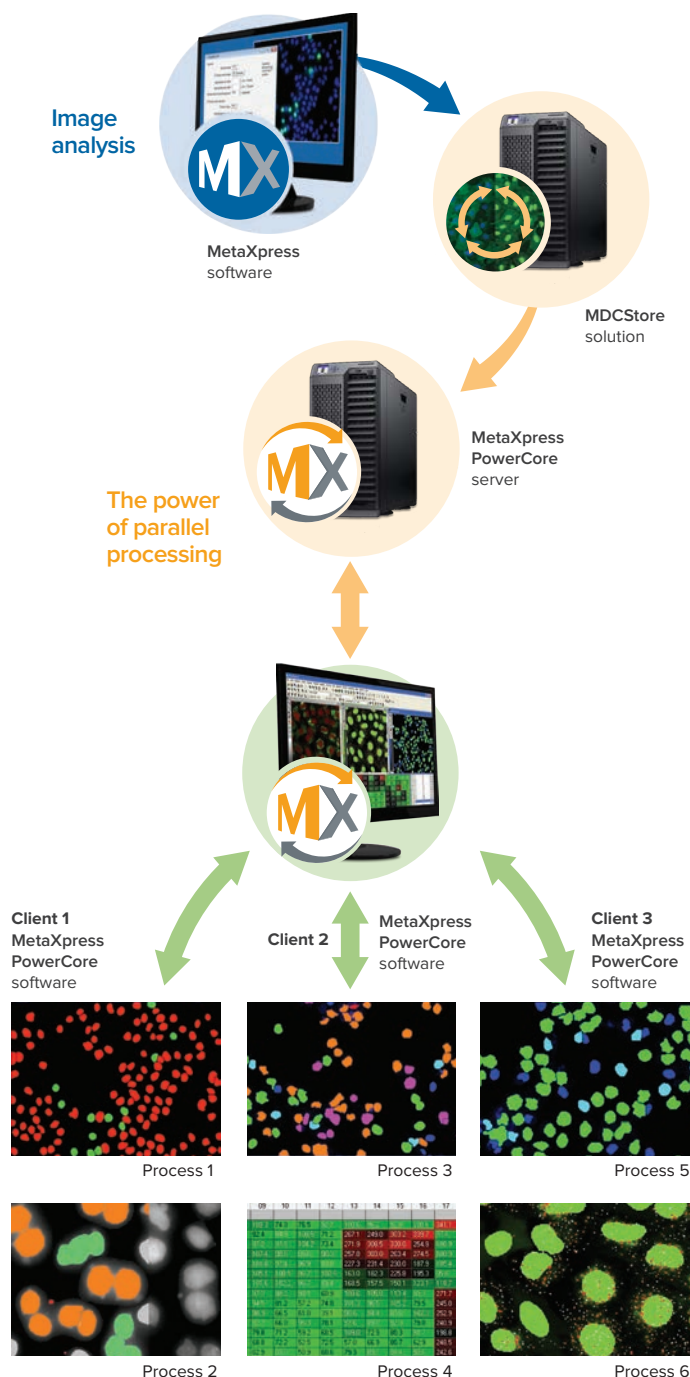
For ultimate customization of acquisition and analysis, journals enable multi-step routines such as whole-well scanning at low magnification and localizing objects of interest. Develop custom journal macros including automatic integration of intelligent hardware control based on your image analysis results.



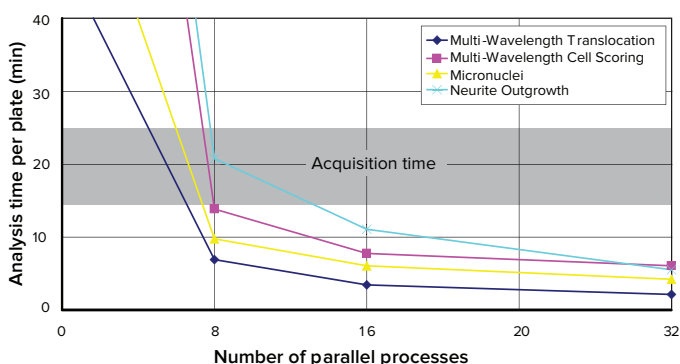
Accelerate

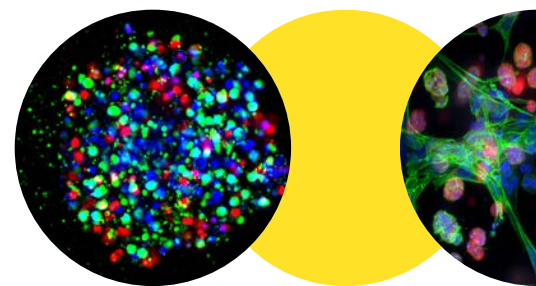
The MetaXpress PowerCore software is a distributed analysis platform with a server that divides plate analysis into small processes and distributes them to multiple clients. Each MetaXpress PowerCore software client, running on multiprocessor computers, executes analysis and relays the results to the server. Built-in flexibility accommodates time and 3D acquired images as well as single plane data. The scalable design accelerates your image analysis at no cost to segmentation quality, overlays, or the large number of analysis outputs. When the analysis is complete, the server efficiently saves all data into the MDCStore database, so results are accessible within your laboratory or across your organization.

Put your HCS capability into overdrive with MetaXpress PowerCore software. Typical basic research or HCS configurations will expedite MetaXpress software application module analysis times by 5x-10x—as fast as acquisition. Using our scalable solution, increase speed further by adding more processors. Core facilities or high-throughput imaging environments can accelerate analysis speeds 10x-40x to support multiple ImageXpress systems running in parallel at full capacity. Similarly, scale up allows re-analysis of archived images at speeds up to 25-fold faster than acquisition.



Analysis faster than acquisition





Ordering information

Our experienced scientific teams are available whenever you have support inquiries, issues, and requests for applications or technical information. When you're ready to take your research to the next level, our applications scientists, trained in the latest technological advancements and assay development techniques, will help get you there.

MetaXpress High-Content Image Acquisition and Analysis Software:

- MDCStore High Content Data Management Solution (included)
- Standard licensing is transferable to the computer of choice
- Network licensing allows multiple users to simultaneously connect to software
- AcuityXpress™ Informatics Software, data visualization, mining, and hit selection are ready to use upon system installation

MetaXpress PowerCore High-Content Distributed Image Analysis software for high-throughput image analysis:

- Various licensing is available to satisfy multiple parallel processes

MetaXpress software application modules:

- Optional modules can be added to the MetaXpress software license

| Product | Part number |
|---|-------------|
| Hardware control options | |
| Digital Confocal option for ImageXpress Micro systems | 5016604 |
| External control for screening acquisition | 9500-0024 |
| File transfer and scheduling | |
| MDCStore Xchange Service | 9500-0159 |
| Application modules | |
| Custom Module Editor | 9500-0158 |
| 3D Analysis (requires Custom Module Editor) | 5047663 |
| Curve fitting | 5058772 |
| Transmitted light segmentation (requires Custom Module Editors) | 5058774 |
| Angiogenesis tube formation | 9500-0016 |
| Cell cycle | 9500-0040 |
| Cell health | 9500-0035 |
| Cell scoring | 9500-0037 |
| Count nuclei | 9500-0017 |
| Granularity | 9500-0032 |
| Live/dead | 9500-0034 |
| Mitotic index | 9500-0036 |
| Micronuclei | 9500-0045 |
| Monopole detection | 9500-0039 |
| Multi-Wavelength cell scoring | 9500-0038 |
| Multi-wavelength translocation | 9500-0044 |
| Neurite outgrowth | 9500-0015 |
| Transfluor® Assay | 9500-0033 |
| Translocation* (includes Translocation-Enhanced*) | 9500-0014 |
| Transfluor HT Assay | 9500-0043 |
| Nuclear translocation HT | 9500-0041 |
| Cell proliferation HT | 9500-0042 |
| All MetaXpress software application modules | 5062036 |

Contact Us

Phone: +1.800.635.5577
 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
 United Kingdom
 +44.118.944.8000
 Europe*
 00800.665.32860

China (Beijing)
 +86.10.6410.8669
 China (Shanghai)
 +86.21.3372.1088
 Hong Kong
 +852.2248.6000

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