



Molecular Devices Imaging – Perfectly Right for Your Research

System Specifications	SpectraMax® i3x Multi-Mode Microplate Reader with MiniMax™ Imaging Cytometer	CloneSelect™ Imager	ImageXpress® Nano Automated Imaging System	ImageXpress® Micro 4 High-Content Imaging System	ImageXpress® Micro Confocal High-Content Imaging System
Measures cell confluency		•			
3 slides or 1 SBS plate/run	•	•	•	•	•
Label free cell counting	•	•			
Workflow for monoclonality verification		•			
4x imaging	•	•	•	•	•
2-color (red/green) + BF imaging	•		•	•	•
Plate reader + imaging in 1 system	•				
2x – 60x imaging			•	•	•
1x – 100x imaging				•	•
Tablet friendly			•		
Timelapse imaging				•	•
Z-stacking				•	•
Environmental control (CO ₂ , O ₂ , humidity, temperature)			•	•	•
UV-VIS widefield fluorescence				•	•
UV-NIR widefield fluorescence				•	•
Automated 3D analysis				•	•
AgileOptix confocal imaging					•
264 slides or 88 SBS plates/run				•	•
On-board liquid handling with simultaneous imaging and dispensing				•	•



Contact your Molecular Devices Technical Sales Specialist and get the perfectly right imager for your lab.

Contact Us モレキュラーデバイスジャパン株式会社 Phone : 0120-993-656 Web : www.moleculardevices.co.jp Email : info.japan@moldev.com	Regional Offices USA and Canada +1.800.635.5577 United Kingdom +44.118.944.8000 Europe* 00800.665.32860 <small>*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland</small>	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	Brazil +55.11.3616.6607
---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------	-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------	-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------	-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------	------------------------------------------------------------

Change to easier imaging for faster discoveries

Powerful and inspirational, our high-content imaging solutions transform your brilliance and intuition into elegant workflows, crystalline images, and compelling data.



Faster discoveries – More relevant data – Less time to image



SpectraMax® i3x Multi-Mode Microplate Reader with MiniMax™ Imaging Cytometer

Multi-Mode Microplate Reader + Basic Imaging

Features

- Softmax Pro software for capture and analysis of plate reader and imaging data
- Customer upgradeable to include reagent injection and western blot detection capabilities

Key Specifications

- 4X Objective Imaging
- 2 Color + Brightfield



CloneSelect™ Imager

Rapid confirmation of monoclonality and subsequent monitoring of cell growth are important for a number of applications including antibody discovery and cell line development.

Features

- Acquire and analyze plates for cell confluency in <90 sec.
- Compatible with 6- to 384-well plates
- Successfully used on suspension and adherent cell lines such as CHO, HEK, hybridomas, Per.C6®, Jurkat, SupT1, H-4-II-E, MCF-7, JT29, DLD1, and KB3
- Can be customized to include fluorescence detection for expanded monoclonality verification and cell based assay workflow options



ImageXpress® Nano Automated Imaging System

Automate Your Cell or Tissue-based Assays

The NANO features Widefield Fluorescence and Label-Free Automated Imaging with user-friendly software that all lab members can use.

Features

- CellReporterXpress™ Software is very easy to use for all lab members
- 20+ pre-built analysis routines
- Access data from a browser – anytime, anywhere
- Tablet friendly

Key Specifications

- Up to 3 Slides or 1 SBS plate per run
- 2X–60X Objective imaging
- UV-VIS Fluorescence
- CO₂, humidity, temperature control



ImageXpress® Micro 4 High-Content Imaging System

High Content Imager featuring Widefield, Fluorescence and Label-Free Automated Imaging Assay Applications

Features

- Software is very flexible, and simple to use, yet powerful
- High-content imaging workhorse that can be highly configured to fit any widefield assay needs
- Automated 3D imaging analysis

Key Specifications

- Up to 264 Slides or 88 SBS plates per run
- 1X–100X objective imaging
- Multi-threaded parallel processing for faster analysis speeds
- Optional on-board liquid handling with simultaneous imaging and compound addition
- UV-NIR fluorescence (customized light sources available)
- Optional CO₂/O₂/N₂, humidity and temperature control
- Integrate other lab components such as incubators, liquid handlers, and plate readers for a for fully automated work cell



ImageXpress® Micro Confocal High-Content Imaging System

Most sophisticated Molecular Devices High Content Imager featuring Widefield + Confocal + Fluorescence + Label-Free Automated Imaging Assay Applications

Features

- Software is very flexible, simple to use, yet powerful
- High content imaging workhorse that can be highly configured to fit any widefield and confocal assay needs
- Automated 3D imaging analysis

Key Specifications

- Up to 264 Slides or 88 SBS plates per run
- 1X–100X objective imaging
- Multiple disk geometries for AgileOptix confocal imaging – flexible to your assay!
- Multi-threaded parallel processing for faster analysis speeds
- Optional on-board liquid handling with simultaneous imaging and compound addition
- UV-NIR fluorescence (customized light sources available)
- Optional CO₂/O₂/N₂, humidity and temperature control
- Integrate other lab components such as incubators, liquid handlers, and plate readers for a for fully automated work cell

SEE HOW MIMETAS, THE ORGAN-ON-A-CHIP COMPANY, USES THE IMAGEXPRESS MICRO CONFOCAL

Mimetas uses their OrganoPlate® to develop disease, toxicology, and transport models for research, development and drug screening. For development and application of tissue models, they use a range of image-based readouts, such as phase-contrast time lapse, fluorescence microscopy, confocal microscopy, in combination with automated 3D image analysis. The team use the ImageXpress Micro Confocal High-Content Imaging System on a daily basis, serving as their standard high-throughput imager for most in-chip assays.

See this story and more at [MolecularDevices.com/Customer-Stories](https://www.moleculardevices.com/customer-stories).

