

Imaging solutions

to answer your biological questions

Stunning images | High throughput | Powerful analysis









ImageXpress HCS.ai

High-Content Screening System



The ImageXpress® HCS.ai systems feature a modular design and a wide array of configuration options, allowing for customization to meet your current needs and on-site upgrades as your requirements evolve.

- Upgraded MetaXpress® Acquire Software for easier, faster image acquisition
- AI-powered IN Carta® Image Analysis Software seamlessly integrated into your workflow
- Exceptional sensitivity with a 95% peak quantum efficiency sCMOS camera
- Easy switching between widefield mode and up to four confocal disc options with AgileOptix™ Spinning Disk Technology
- Six objective turret alongside optional magnification changer giving up to 12 effective levels of magnification
- Configurable system with a simple upgrade pathway to match your current and future research needs

ImageXpress HCS.ai WIDEFIELD	ImageXpress HCS.ai CONFOCAL	ImageXpress HCS.ai ADVANCED
	Widefield capability	
	2X–60X high NA and 20X–60X ELWD objective options plus a magnification changer option (1X and 1.5X)	
	LED illumination, 5 color LED (up to 220 mW per channel)	Laser illumination, 7 color laser (up to 800 mW per channel)
	DAPI, FITC, TRITC, TxRed, Cy5 (+ CFP, YFP as optional)	DAPI, CFP, FITC, YFP, TRITC, TxRed, Cy5, Cy7
		Standard spinning disk confocal 60 µm pinhole (single), 60 µm pinhole + 50 µm slit (dual)
		 Deep tissue spinning disk 50 µm pinhole high resolution and high-sensitivity disk (dual)
	Water immersion option (20X–60X)	
	Environmental control option (temperature, humidity, CO ₂ , O ₂)	

Upgradable modular design



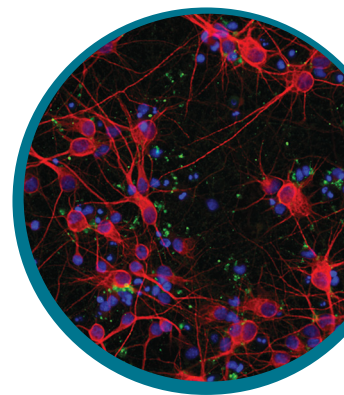
***Available models:**

ImageXpress Micro 4,
ImageXpress Micro Confocal,
ImageXpress Confocal HT.ai

ImageXpress High-Content Imaging Systems

High-content imaging systems with automated compound addition

- Integrated fluidics addition with tip-based liquid addition
- 2x sample plate capacity with temperature control
- Fast image acquisition for 2D, 3D, and kinetic assays
- AgileOptix Spinning Disk Confocal Technology
- Water immersion, laser illumination, and environmental control options available
- Analyze images with AI-powered IN Carta Image Analysis Software

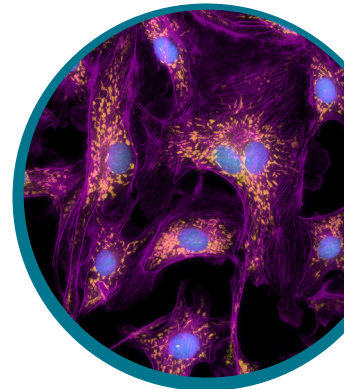


*ImageXpress Pico Digital Confocal uses AutoQuant 2D Real-Time Deconvolution

ImageXpress Pico Automated Cell Imaging System

Compact system that allows users to go from samples to results in minutes

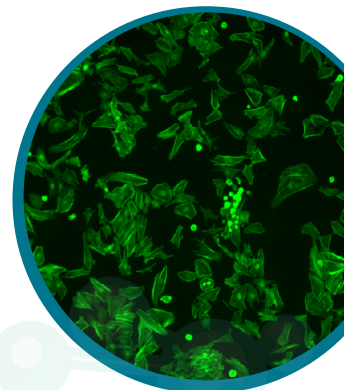
- Ideal for cell counting, transfection efficiency, and cell health assays
- 25+ preconfigured application protocols
- 3D z-stack acquisition
- On-the-fly analysis
- Environmental control for live cell assays
- Access data from a browser—anytime, anywhere
- Optional Digital Confocal on-the-fly 2D deconvolution*



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

Integrated multi-mode microplate reader and imaging system

- Ideal for cell counting, label-free imaging, and cell proliferation
- Western blot, imaging, and injectors on one configurable microplate reader
- Stain free, automated cell counting
- Preset plate reader and imaging protocol library
- Advanced curve fitting and statistical analysis



System specifications



SpectraMax i3x
Multi-Mode
Microplate Reader
with MiniMax
Imaging Cytometer



ImageXpress Pico
Automated Cell
Imaging System



ImageXpress HCS.ai
High-Content Imaging Systems

Product features			Widefield	Confocal	Advanced
Microplate types	6 to 1536 wells for plate reader 96- and 384-wells for imaging	Up to 384 wells	Up to 1536 wells		
Number of fluorescent channels	2	5	5–7	5–7	8
Camera technology	CCD	CMOS	Back thinned scientific CMOS		
Light source	1 LED	5 LED	5 LED	5 LED	7 LASER
Brightfield transmitted light	•	•	•	•	•
Magnification range*	4X	4–63X	2–90X		
Objective positions	1	4	6		
Water immersion option			•	•	•
Environmental control for CO ₂ , humidity and temperature control option		•	•	•	•
Z-stack acquisition		•	•	•	•
Confocal capability		Digital*	Digital*	•	•
Deep tissue spinning disk					•
Automated 3D analysis with volumetric output				•	•
Acquisition software	SoftMax® Pro	CellReporterXpress®	MetaXpress Acquire		
Analysis software	SoftMax Pro	CellReporterXpress (+export to IN Carta)	IN Carta		

*with automated magnification changer option

Please note: not all models are shown. For specific requirements including automated compound addition, please contact a member of our team.

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	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.120.993.656
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586	India	+1.800.266.5338

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom

