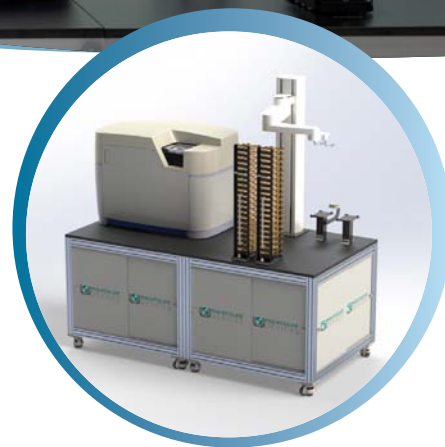


Deeper insights
for every lab



Extend the power of the ImageXpress HCS.ai system with scalable walkaway automation

The ImageXpress® HCS.ai High-Content Screening System with automation advances high-throughput imaging by integrating robotic plate handling and other flexible configurations. Enhance your screening by reducing manual workload, boosting reproducibility and throughput, and enabling you to perform new assays. Capture real-time cellular dynamics and perform fully automated time-course studies with improved accuracy.



The ImageXpress HCS.ai system with a dedicated, scalable automation workstation for expanded capacity.

Key capabilities



Image more plates in less time – Process 40 microtiter plates (96-well format) in just 2 hours—or 80 plates in 4 hours—with complete hands-off operation.*



Unlock new assays – Perform fully automated high-throughput screens or long-term, live cell assays



Streamline high-throughput imaging – Enhance the ImageXpress HCS.ai High-Content Screening System by reducing manual workload, improving reproducibility, and freeing up hands-on time.



Boost confidence in your data – Minimise human errors and get the reliable data you need through consistent, precise and reproducible imaging.



Simplify system installation – Reduce training time and screen from day one. Our team handles all aspects of system installation and integration – complete with the protocols you will need for your research.



Flexible software integration – Our system is compatible with three widely used software packages, allowing your lab to choose the software environment that best matches your experience, requirements, and budget.








Optimize the use of lab space – Fit seamlessly into your existing lab benches or workflows, minimizing disruption and maximizing efficiency.







Future-proof your investment – Start with a compact setup and expand your system as your scientific discovery evolves.

*Imaging speed an entire 96-well plate with two channels in under 90 seconds with 2D acquisition: 96-well, 1 FOV, 1 plane, DAPI (14ms), FITC (10ms).

Choose the right configuration for your workflow

Automated Set Up	Parts Included	Storage Capacity	System Dimensions
	ImageXpress HCS.ai FlexiLab <ul style="list-style-type: none"> • PF400 Robot Arm • 2 Universal Hotels (40 plate capacity) • Barcode Reader with Barcode Reader Station • Lid Station 	40 Microtiter plates at room temperature	36" x 84" 92cm x 214cm
	ImageXpress HCS.ai StoreMax <ul style="list-style-type: none"> • PF400 Robot Arm • 4 Universal Hotels (80 plate capacity) • Barcode Reader with Barcode Reader Station • Lid Station 	80 Microtiter plates at room temperature	36" x 84" 92cm x 214cm
	ImageXpress HCS.ai EnviroLab <ul style="list-style-type: none"> • PF400 Robot Arm • 4 Universal Hotels (80 plate capacity) • Barcode Reader with Barcode Reader Station • Lid Station • STX44BT – Automated Incubator (44 Plates Capacity) 	<ul style="list-style-type: none"> • 80 Microtiter plates at room temperature • 44 Microtiter plates at 37°C 5%, 95% RH, CO₂ 	36" x 84" 92cm x 214cm
	ImageXpress HCS.ai BaseStoreMax <ul style="list-style-type: none"> • PF400 Robot Arm • 4 Portrait Hotels (80 plate capacity) • Barcode Reader with Barcode Reader Station • Lid Station 	80 Microtiter plates at room temperature	24" x 73" 60cm x 186cm
	ImageXpress HCS.ai BaseEnviroLab <ul style="list-style-type: none"> • PF400 Robot Arm • 3 Portrait Hotels (60 plate capacity) • Barcode Reader with Barcode Reader Station • Lid Station • STX44BT – Automated Incubator (44 Plates Capacity) 	<ul style="list-style-type: none"> • 60 Microtiter plates at room temperature • 44 Microtiter plates at 37°C 5%, 95% RH, CO₂ 	24" x 96" 60cm x 243cm

Hardware

	PreciseFlex™*** PF400 <ul style="list-style-type: none"> • PF400 – 750 mm Z, 730 mm reach • PF400 SCARA is ISO-015066 compliant
	Ambient Storage Universal Plate Hotels <ul style="list-style-type: none"> • Universal Plate hotels for standard height SBS plates • Each hotel has 20 plate positions • Ambient hotel provides random access storage for microtiter plates. • The Universal hotel features a docking station, allowing for easy movement and secure placement of the hotel rack when needed. • Compatible with 6-well plates to 1536-well plates (max height 28mm) • The hotel is designed for easy removal of alternate nest locations, accommodating a deep well
	Ambient Storage Portrait Hotel <ul style="list-style-type: none"> • Portrait ambi hotel provides random access storage for microtiter plates. • Rear / Front Plate Feed • Each hotel has 20 plate positions • Compatible with 6-well plates to 1536-well plates (max height 28mm) • The hotel is designed for easy removal of alternate nest locations, accommodating a deep well
	Liconic™**** Automated STX44 – CO₂ Incubator <ul style="list-style-type: none"> • Benchtop model: 33°C to 50°C • 95% relative humidity, high speed handling • Plate hotels can be configured as follows: <ul style="list-style-type: none"> • 22 plate capacity (plate height up to 17mm), quantity 2. Max incubator capacity of 44 plates. • 17 plate capacity (plate height up to 23mm), quantity 2. Max incubator capacity of 34 plates.



Discover how the **ImageXpress HCS.ai High-Content Screening System** can accelerate your research today. Contact our expert team to learn more.
moleculardevices.com/imagepress-hcs-ai

Contact Us

Phone: +1.800.635.5577
Web: www.moleculardevices.com
Email: info@moldev.com

Regional Offices

USA and Canada	+1.800.635.5577	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
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

Check our website for a current listing of worldwide distributors.

