



## Automation solutions for high-throughput plate-based assays

Explore our fully-integrated, automated workflow solutions for cell and biochemical assays

As timelines compress and the amount of data required for IND submissions and publications increases, laboratories are pressed to do more with less. The introduction of automated work cells can help relieve this burden by allowing scientists to optimize their time in the lab and obtain data more efficiently.

We can build flexible, scalable, future-proof workflows backed by decades of consultative expertise to help you design and implement work cells that incorporate industry-leading technology to achieve your goals.

### ELISA Workcell

Walk away solution for ELISA's, cell based work and more

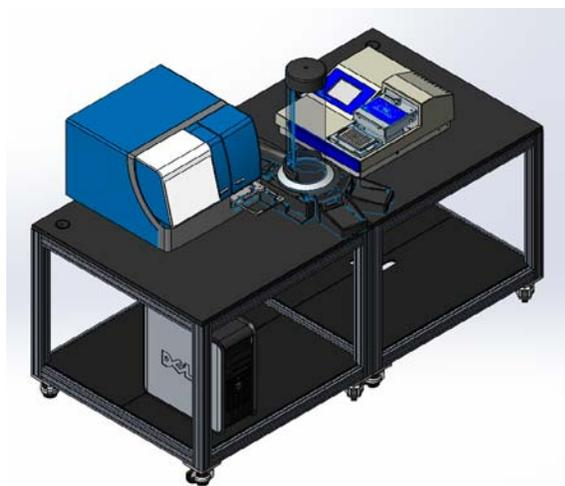


Plate Reader	SpectraMax iD5 multi-mode system
Plate Washer	AquaMax 4000 Microplate Washer
Robot	Peak Robotics S-LAB Plate Handler
Plate Storage	Plate Hotels – 40 microtiter plates
Controlling Software	Peak Robotics S-LAB Software

[Learn more](#)

### ELISA Advanced Workcell

Compact and flexible solution for a wide range of HTS work

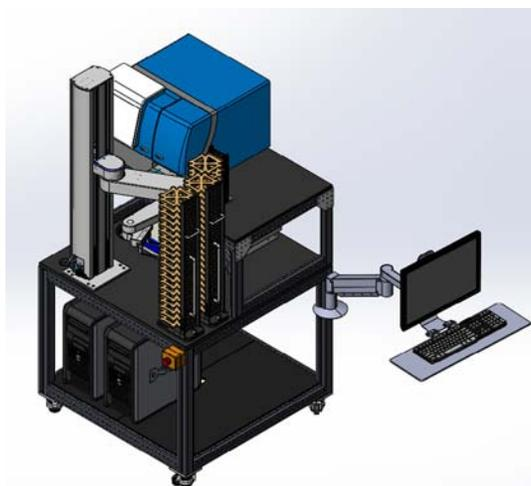


Plate Reader	SpectraMax iD5 multi-mode system
Plate Washer	AquaMax 4000 Microplate Washer
Robot	PreciseFlex 400 Sample Handler
Plate Storage	Plate Hotels – 40 microtiter plates
Controlling Software	Genera Control Software

[Learn more](#)

## ELISA Advanced EC Workcell

Incubator included to expand your capability

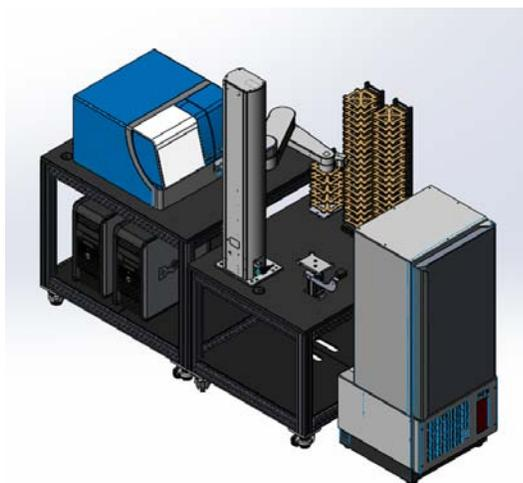


Plate Reader	SpectraMax iD5 multi-mode system
Plate Washer	AquaMax 4000 Microplate Washer
Robot	PreciseFlex 400 Sample Handler
Plate Storage	Plate Hotels – 40 microtiter plates
Controlling Software	Genera Control Software
Incubation	LiCONiC STX44 Automated CO <sub>2</sub> Incubator

[Learn more](#)

## ELISA HTS EC Workcell

Great for HTS assay work covering sample prep to analysis

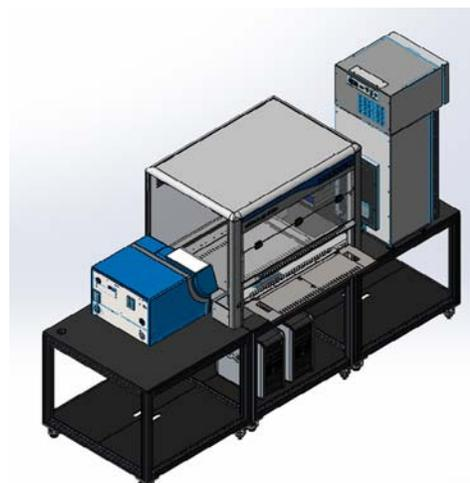


Plate Reader	SpectraMax iD5 multi-mode system
Plate Washer	AquaMax 4000 Microplate Washer
Incubation	LiCONiC STX44 Automated CO <sub>2</sub> Incubator
Liquid Handler	Microlab STAR Automated Liquid Handling

[Learn more](#)

## ELISA HTSx Workcell

Flexible and fast. No limit to throughput or capability

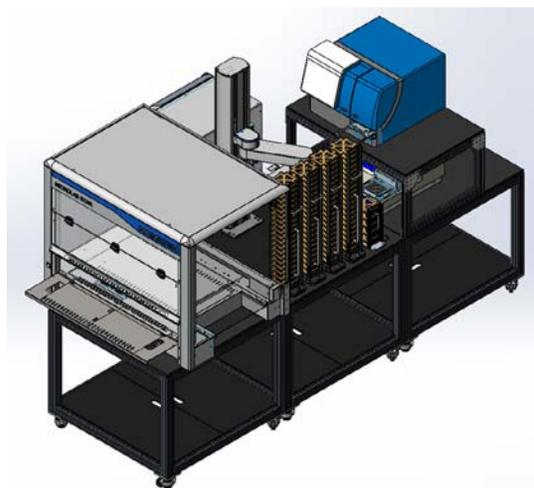


Plate Reader	SpectraMax iD5 multi-mode system
Plate Washer	AquaMax 4000 Microplate Washer
Robot	PreciseFlex 400 Sample Handler
Plate Storage	Plate Hotels – 40 microtiter plates
Controlling Software	Genera Control Software
Incubation	LiCONiC STX44 Automated CO <sub>2</sub> Incubator
Liquid Handler	Microlab STAR Automated Liquid Handling

[Learn more](#)

### Contact Us

Phone: +1.800.635.5577  
 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.

### Regional Offices

USA and Canada +1.800.635.5577  
 United Kingdom +44.118.944.8000  
 Europe\* 00800.665.32860  
 China +86.4008203586  
 Taiwan/Hong Kong +886.2.2656.7585  
 Japan +81.3.6362.9109  
 South Korea +82.2.3471.9531  
 India +91.73.8661.1198

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