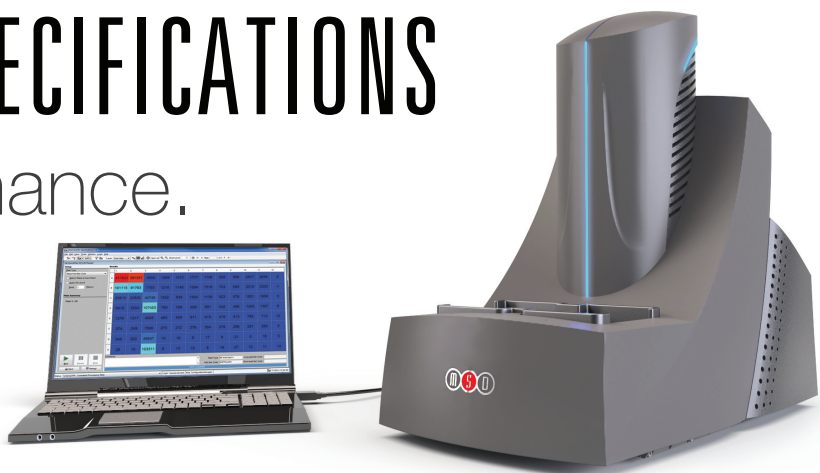




MESO[®] QuickPlex SQ 120

FEATURES AND SPECIFICATIONS

Affordable Performance.
More Results.



The QuickPlex[®] instrument offers affordable access to high-performance electrochemiluminescence immunoassays. This compact system has been engineered for reliability, ease of use, and low cost of ownership. The combination of rapid read times and the ability to perform multiple, simultaneous tests on a single sample increases productivity, saves sample, and delivers results quickly. The QuickPlex instrument has a wide menu of commercially available assay kits and a full line of components and reagents for developing your own assays.

Affordability

- MSD's most affordable ECL imager
- Special pricing for academic customers – please inquire
- Low cost of ownership

Assay Menu

- Wide menu of commercially available kits
- Full line of components for user-developed assays
- Fully compatible with all SECTOR[®] products

Validated

- Developed under design control
- Microsoft Windows compatible
- Simple, fast, and secure data analysis with DISCOVERY WORKBENCH[®] software

Performance

- ECL detection for superior sensitivity and dynamic range
- Single and multiplex assays
- Single-well read mode for enhanced performance

Compact Size

- Small footprint – ¼ the size of MSD's SECTOR instruments
- Simple set up, no special requirements
- Full-featured laptop

Convenience and Reliability

- Rapid read times (as fast as 90 seconds per plate)
- No fluidics – nothing to clog or clean
- No daily or monthly user maintenance





MESO QuickPlex SQ 120 Specifications

Instrument Name	MESO QuickPlex SQ 120	
Instrument Catalog No.	A10AA-0	
Detection Technology	Electrochemiluminescence (ECL)	
Detection System	Cooled scientific-grade CCD camera and precision custom lens assembly	
Multiplexing Capability	Up to 10 assays/well	
Plate Compatibility	MSD® 96-well SECTOR plates MSD QUICKPLEX® plates	
Sensitivity	Limits of detection as low as 0.05 to 1 pg/mL for many assays	
Dynamic Range	Greater than 10 ⁵	
Read Volume (96-well plates)	150 µL	
Plate Read Time	SECTOR plates	1 min 30 sec
(Singleplex to 10-spot multiplex plates)	QUICKPLEX plates	2 min 45 sec
Throughput	SECTOR plates	40 plates/hr
	QUICKPLEX plates	22 plates/hr
Plate Stacker Capability	5 plates input stack / 5 plates output stack	
Operating Temperature Range	68°F to 78°F (20°C to 26°C)	
Operating Humidity Range	10% to 80% relative humidity (non-condensing)	
Maximum Altitude	6,500 ft. (1,900 meters)	
Storage Conditions	0°F (-18°C) @ 50% relative humidity to 122°F (50°C) @ 85% relative humidity (non-condensing)	
Measured Sound Level	59 dbA	
Weight	44.5 lbs. (20.2 kg)	
Size (W x D x H)	13 in. x 23 in. x 21 in. (33 cm x 58 cm x 53 cm)	
Power	100–240 VAC, 50–60 Hz / 170W	
Regulatory	For research use only. Not for use in diagnostic procedures. Meets UL, CSA, FCC and applicable IEC safety and EMC standards and carries the CE and cSGSus marks in addition to the Australian regulatory compliance mark. Applicable China RoHS information is also found on the instrument label and in the user's manual. Please contact MSD for more information on regulatory certifications.	
User Maintenance	None required. Annual maintenance contracts are available for purchase and recommended for ensuring long-term value of the instrument	
Warranty	One year limited warranty on parts and labor. Service plans for extended warranty are available for separate purchase.	

Phone: 1-240-314-2795 . Fax: 1-301-990-2776 . Email: customerservice@mesoscale.com

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, MSD GOLD, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, R-PLEX, S-PLEX, U-PLEX, V-PLEX, STREPTAVIDIN GOLD, MESO, www.mesoscale.com, SMALL SPOT (design), 96 WELL 1, 4, 7, 9, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design), R-PLEX (design), S-PLEX (design), U-PLEX (design), V-PLEX (design), It's All About U, and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. All other trademarks or service marks are the property of their respective owners.

©2013, 2017 Meso Scale Diagnostics, LLC. All rights reserved.



For Research Use Only. Not for use in diagnostic procedures.

MK-DS-185-v3-2017Jul