



MESO[®] QuickPlex SQ 120MM

FEATURES AND SPECIFICATIONS

Greater Flexibility.
More Results.



The MESO QuickPlex[®] SQ 120MM is MSD's most versatile instrument. 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.

Broad Assay Menu

- Large menu of commercially available kits
- Full line of components for user-developed assays
- Access to all product lines for MSD[®] consumables

Flexibility

- Single and multiplex assays
- Compatible with all SECTOR[™] and QuickPlex[®] products
- Special pricing for academic customers

Validated

- Developed under design control
- Microsoft Windows 10 compatible
- Simple, fast, and secure data analysis with DISCOVERY WORKBENCH[®] software

Performance

- ECL detection for superior sensitivity and dynamic range
- Rapid read times (as fast as 90 seconds per plate)
- Single-well read mode for enhanced performance

Compact Size

- Small footprint— $\frac{1}{4}$ the size of MSD's SECTOR[®] instruments
- Simple set up, no special requirements
- Full-featured laptop

Convenience and Reliability

- No fluidics—nothing to clog or clean
- No daily or monthly user maintenance
- IQ/OQ/PQ, Preventative Maintenance, and Service Packages available





MESO QuickPlex SQ 120MM Specifications

Instrument Name	MESO QuickPlex SQ 120MM	
Instrument Catalog Number	AI1AA-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	
Plate Read Time	SECTOR plates	1 minute 30 seconds
(Singleplex to 10-spot multiplex plates)	QuickPlex plates	2 minutes 45 seconds
Throughput	SECTOR plates	40 plates/hour
	QuickPlex plates	22 plates/hour
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 (2,000 m)	
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 lb (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, 155 W max.	
Network	A wired or wireless network connection to the instrument laptop is required. Access to the *.methodicalmind.com™ and *.amazonaws.com websites through port 443 must be allowed through the company firewall.	
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 safety guide. Please contact MSD for more information on regulatory certifications.	
Warranty	One year limited warranty on parts and labor. Service plans for extended coverage are available for separate purchase.	

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

MESO SCALE DIAGNOSTICS, MSD, DISCOVERY WORKBENCH, QuickPlex, SECTOR, MESO, www.mesoscale.com, methodicalmind.com, 96 WELL 4-SPOT (design), MSD (design), 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, 2019 Meso Scale Diagnostics, LLC. All rights reserved.



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

MK-DS-910-v2-2020Jan

