



MESO[®] SECTOR S 600MM

FEATURES AND SPECIFICATIONS

Ultra-Sensitive Plate Imager
With Industrial Capacity



The MESO SECTOR[®] S 600MM is our flagship instrument. Its high-speed, ultra-low noise CCD camera provides exceptionally high sensitivity, wide dynamic range, and rapid read times. Developed under rigorous design control and MSD's certified Quality Management System, the MESO SECTOR S 600MM delivers the industrial-grade throughput and reliability required in the most demanding research environments.

Performance

- ECL detection for superior analytical performance
- Six logs of dynamic range
- Highly sensitive, large format imaging system
- 96-well and 384-well plate capability

Validated

- Developed under design control
- Microsoft Windows 10 compatible
- Simple, fast, and secure data analysis with DISCOVERY WORKBENCH[®] software
- Service, maintenance, and IQ/OQ/PQ plans available

Broad Assay Menu

- Extensive menu of off-the-shelf kits
- Singleplex and multiplex assays
- Access to all product lines for SECTOR[™] consumables
- Expert in-lab support for user-developed assays

Convenience and Reliability

- No fluidics—nothing to clean or clog
- High capacity stackers for walk-away use
- Rapid read times: 70 seconds per plate
- No daily or monthly user maintenance



MESO SECTOR S 600MM Specifications

Instrument Name	MESO SECTOR S 600MM
Instrument Catalog Number	IC1AA-0
Detection Technology	Electrochemiluminescence
Detection System	Cooled scientific-grade CCD camera and precision custom lens assembly
Multiplexing Capability	Up to 10 assays/well
Plate Compatibility	MSD® 96- and 384-well SECTOR plates
Sensitivity	Limits of detection as low as 0.05 to 1 pg/mL for many assays
Dynamic Range	Greater than 10 ⁶
Plate Read Time	70 sec (all singleplex and multiplex plates)
Throughput	50 plates/hr
Plate Stacker Capacity	Tall stack tube: 50 (96-well) or 75 (384-well) plates; short stack tube: 20 or 35 plates
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)
Weight	124 lb (56 kg)
Size (W x D x H)	27 in x 31 in x 30 in (69 cm x 79 cm x 76 cm)
Power	100–240 VAC, 50–60 Hz, 126 W max.
Network	A wired or wireless network connection to the instrument computer 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, SECTOR, MESO, www.mesoscale.com, methodicalmind.com, V-PLEX, 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.
©2014, 2017, 2019 Meso Scale Diagnostics, LLC. All rights reserved.



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

MK-DS-912-v2-2020Jan

