MESO® SECTOR S 600

FEATURES AND SPECIFICATIONS Ultra-Sensitive Plate Imager With Industrial Capacity

- Singleplex and multiplex formats
- Six-log dynamic range
- 96-well and 384-well plate capability

The MESO SECTOR[®] S 600 reader 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 600 delivers the industrial-grade throughput and reliability required in the most demanding research environments.

Performance

- ECL detection for superior analytical performance
- Rapid read times: 70 seconds per plate, up to 50 plates per hour
- Highly sensitive, large format imaging system
- 384-well multiplexing for high throughput and efficiency

Assay Menu

- Extensive menu of off-the-shelf kits
- Singleplex and multiplex assays
- V-PLEX[®] validated biomarker assays
- Expert in-lab support for user-developed assays

Validated

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

000

R

- Microsoft Windows compatible
- IQ/OQ/PQ plans available

Convenience and Reliability

- No fluidics—nothing to clean or clog
- High capacity stackers for walk-away use
- Multiplexing without compromise—equivalent performance with singleplex or multiplex assays
- No daily or monthly user maintenance



MESO SECTOR S 600 Specifications

Instrument Name	MESO SECTOR S 600
Instrument Catalog No.	ICOAA-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
Read Volume (96-well plates)	150 μL
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	155 lb (70 kg) for the installed instrument with computer and accessories
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.
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, DISCOVERY WORKBENCH, SECTOR, MESO, www.mesoscale.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. ©2014, 2017, 2019 Meso Scale Diagnostics, LLC. All rights reserved.



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

MK-DS-215-v3-2019Jul

