



# 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 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 ISO 9001:2008 certified Quality Management System, the 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

### Validated

- Developed under design control
- Simple, fast, and secure data analysis with DISCOVERY WORKBENCH® software
- Windows 7 compatible
- IQ/OQ/PQ plans available

### Assay Menu

- Extensive menu of off-the-shelf kits
- Singleplex and multiplex assays
- Validated with V-PLEX™ biomarkers
- 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
- 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.	IC0AA-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 <sup>6</sup>
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 (1900 m)
Storage Conditions	0°F (-18°C) @ 50% relative humidity to 122°F (50°C) @ 85% relative humidity (non-condensing)
Weight	155 lbs (70 kg) for the installed instrument with computer and accessories
Size (W x D x H)	26.8 in x 30.5 in x 29.5 in (68.1 cm x 77.5 cm x 74 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 cTUVus marks. 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.

MSD's Quality Management System is certified under ISO 9001:2008 for the design, development, manufacture, distribution and service of instruments, plates, reagents, assays, and kits for research purposes.

**Phone: 1-240-314-2795 . Fax: 1-301-990-2776 . Email: [customerservice@mesoscale.com](mailto:customerservice@mesoscale.com)**

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, V-PLEX, STREPTAVIDIN GOLD, MESO, [www.mesoscale.com](http://www.mesoscale.com), SMALL SPOT (design), 96 WELL 1, 4, 7, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

©2014 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-215-v1-2014Feb