



Serious Tools for Serious Science

# MESO® QuickPlex Q 60MM

The MESO Quickplex™ Q 60 MM instrument delivers 5 logs of dynamic range, attomolar sensitivity, multiplexing capability and access to a menu of over 600 assays.

Ask about our academic program, with special pricing and terms to help make the Q 60MM accessible to academic researchers.



## Over 600 Assays with Multiplexing Capabilities

### R-PLEX®

Fuel your research with singleplex assays with better results than traditional ELISAs

### U-PLEX®

Flexibility to customize multiplex assays and preconfigured application-specific sets

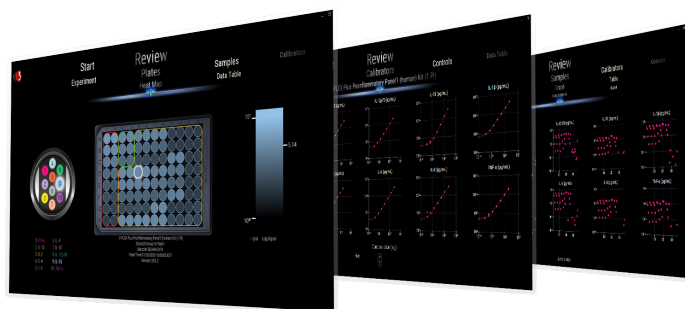
### V-PLEX®

Validated assays with consistent and reliable quantification every time

### S-PLEX®

Ultrasensitive assays deliver fg/mL detection of trace biomarkers or small sample volumes

## The Methodical Mind Experience



- Seamless integration with MSD® instruments for a streamlined workflow
- Cloud-based connectivity enables collaboration with easy access and sharing of data, assay protocols, and analysis templates
- Convenient and secure tool for managing users and teams. Built in audit trails for compliance.



www.mesoscale.com

# Features and Benefits

## High Performance Technology

- Industry leading sensitivity – 5 logs of dynamic range for precise quantification of even the most difficult sample types
- QuickPlex Ultra technology ensures fast read times - get your results in 83 seconds

## Integrated Workflow

- Built-in compatibility among consumables, hardware, and software
- Automatic integration of product information with data analysis makes reporting fast and easy

## Flexible Assays

- Large menu of pre-configured and customizable assays, kits, and multiplex panels
- Assay development components and tools to build your own assays

## Responsive and Expert Support

- Global team of field application scientists and service engineers for on-site support
- Highly skilled technical support team

## Robust, Worry-Free Instrumentation

- No daily or monthly calibration cycles
- Requires an internet connection for operation

## Instrument Service

- Service plan options to suit every budget
- No microfluidics means minimal maintenance

Instrument Name	MESO QuickPlex Q 60MM
Instrument Catalog Number	IB0AA-1
Detection Technology	Electrochemiluminescence (ECL) using cooled scientific-grade CCD camera and precision custom lens assembly
Plate Compatibility	MSD 96-well QuickPlex and QuickPlex Ultra™ plates
Sensitivity	Limits of detection as low as 0.05 to 1 pg/mL for many assays
Dynamic Range	10 <sup>5</sup> ECL counts
Plate Read Time	QuickPlex Ultra plates: 1 minute 23 seconds
Plate Stacker Capability	5 plates on the input stack and 5 plates on the output stack
Operating Environment	68°F to 78°F (20°C to 26°C) at 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	36 lb (16 kg)
Size (W x D x H)	13 in x 23 in x 14 in (33 cm x 58 cm x 36 cm)
Power	100–240 VAC, 50/60 Hz, 150 W max.
Network	The instrument's computer must be connected to the network using either the wireless or wired network adapter. The Methodical Mind® software requires an internet connection to the methodicalmind.com™ domain via HTTPS on TCP port 443.
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, UKCA, and cSGSUS marks in addition to the Australian regulatory compliance mark (RCM) and crossed out wheelie bin symbol with bar (WEEE). Applicable China RoHS information is also found on the instrument label and in the safety guide. Complies with FDA performance standards for laser products except for conformance with IEC 60825-1 Ed. 3., as described in Laser Notice No. 56, dated May 8, 2019. 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](mailto:customerservice@mesoscale.com)

MSD, [www.mesoscale.com](http://www.mesoscale.com), [methodicalmind.com](http://methodicalmind.com), MESO, Methodical Mind, QuickPlex, QuickPlex Ultra, R-PLEX, S-PLEX, U-PLEX, V-PLEX, and MSD (design) are all trademarks and/or service marks owned by or licensed to Meso Scale Diagnostics, LLC. All other trademarks and service marks are property of their respective owners.  
©2025 Meso Scale Diagnostics, LLC. All rights reserved.

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

D0000-3996-A