



Site Requirements for the MESO[®] SECTOR S 600MM, MESO QuickPlex[®] SQ 120MM, and MESO QuickPlex Q 60MM

This document outlines the site requirements for the installation of the MESO SECTOR[®] S 600MM, MESO QuickPlex SQ 120MM, and MESO QuickPlex Q 60MM instruments. These site requirements also apply to MESO SECTOR S 600 and MESO QuickPlex SQ 120 instruments running DISCOVERY WORKBENCH[®] software that will be upgraded to Windows 10.

Space Requirements

All systems include the instrument, an uninterruptible power supply (UPS), cables, and a computer. The MESO SECTOR S 600MM is supplied with a desktop computer and a touch screen monitor. The MESO QuickPlex SQ 120MM and MESO QuickPlex Q 60MM instruments are supplied with a laptop computer.

The instrument must be installed in a location where enough room is available around the instrument for adequate airflow for heat dissipation. The MESO SECTOR S 600MM requires 3 inches of space all of the way around the instrument. The MESO QuickPlex SQ 120MM and the MESO QuickPlex Q 60MM require 2 inches of space behind and 1 inch on each side of the instrument. The installation location must not be exposed to intense light such as direct sunlight, and it must not be in the direct path of heating and cooling ducts. The following table lists the space needed for the instrument as well as the overall system (instrument, UPS, and computer).

Instrument Name	Instrument Size (width x depth x height)	System Size (width x depth x height)
MESO SECTOR S 600MM	27 in x 31 in x 30 in (69 cm x 79 cm x 76 cm)	62 in x 34 in x 36 in (158 cm x 86 cm x 91 cm)
MESO QuickPlex SQ 120MM	13 in x 23 in x 21 in (33 cm x 58 cm x 53 cm)	30 in x 30 in x 21 in (76 cm x 76 cm x 53 cm)
MESO QuickPlex Q 60MM	13 in x 23 in x 14 in (33 cm x 58 cm x 36 cm)	30 in x 30 in x 14 in (76 cm x 76 cm x 36 cm)

Power Requirements

MSD[®] instruments are compatible with single-phase power supplied at 100–240 volts of alternating current and a frequency of 50 or 60 Hz. Ensure the location where the system will be installed has a power outlet within 4 ft (1.2 m) of this location. The table below indicates the peak power consumption for each system (instrument, UPS, and computer).

Instrument Name	Peak Power Consumption
MESO SECTOR S 600MM	231 watts
MESO QuickPlex SQ 120MM	350 watts
MESO QuickPlex Q 60MM	120 watts

Software Requirements

The Methodical Mind[™] software requires internet connectivity to the methodicalmind.com[™] and amazonaws.com websites through the company firewall. Enable the *.methodicalmind.com and *.amazonaws.com domains through TCP port 443.