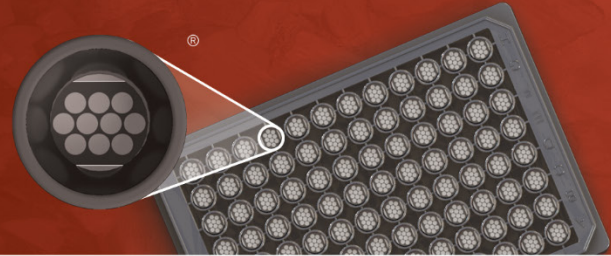


MSD™ Serology Control Pack 1



Ordering Information

MSD Customer Service
Phone: 1-240-314-2795
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com
www.mesoscale.com/support

Scientific Support

Phone: 1-240-314-2798
Email: ScientificSupport@mesoscale.com
www.mesoscale.com/support

Company Address

MESO SCALE DISCOVERY®
A division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173
USA

www.mesoscale.com®

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

Catalog Number	C4381-1
Contents	Serology Control 1.1 (1mL frozen liquid per vial) Serology Control 1.2 (1mL frozen liquid per vial) Serology Control 1.3 (1mL frozen liquid per vial)
Storage	≤-70 °C

Summary and Intended Use

The Serology Control Pack 1 consists of three levels of serum-based controls for use in the V-PLEX® COVID-19 Serology Kits. Each control contains assigned concentrations of human IgG, IgM, and IgA against antigens in the V-PLEX COVID-19 Serology Kits. For assigned values in MSD arbitrary units (AU/mL) and WHO international units in (BAU/mL), refer to the V-PLEX COVID-19 Serology Kits product insert or Serology Control Pack 1 lot-specific certificate of analysis (COA).

Storage and Handling

To maximize the consistency of measured values across vials, the controls must be stored at the recommended temperature. To use, first thaw the frozen control vials on ice and equilibrate to room temperature before loading into the plate. Vortex briefly and spin down before loading controls into the plate. Each control is supplied at the working concentration. Do not dilute prior to use.

Refer to the assay-specific product insert for control preparation and instruction for use.

Note: Stock control is stable for 5 years from the date of manufacture when stored at ≤-70 °C. Thawed control is stable through five freeze-thaw cycles.

Safety

Serology Control Pack 1 contains human serum and is a biosafety level 2 (BSL-2) product. Donors have been tested and found negative for HBsAg, HIV-1 and HIV-2 antibodies, and HCV. Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling controls. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines. Additional product-specific safety information is available in the safety data sheet (SDS), which can be obtained from MSD Customer Service or at www.mesoscale.com.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, www.mesoscale.com, MSD (design), V-PLEX, 96 WELL 10-SPOT (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

©2020, 2021 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-1006-v2-2021Mar