

MSD[®] Proinflammatory Panel 1 (human) Control Pack



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 #	C4049-1
Contents	Proinflammatory Panel 1 (human) Control 1 (5 vials,* lyophilized) Proinflammatory Panel 1 (human) Control 2 (5 vials,* lyophilized) Proinflammatory Panel 1 (human) Control 3 (5 vials,* lyophilized) <small>*If supplied as part of a V-PLEX[®] Plus kit, quantity will match number of plates ordered.</small>
Storage	2–8°C

Summary and Intended Use

The use of controls to monitor the analytical performance and reliability of test methods is an essential component of good laboratory practice.

The Proinflammatory Panel 1 (human) Control Pack consists of 3 levels of controls, each containing known concentrations of all human cytokines that are detected by the Proinflammatory Panel 1 (human) Kit. Proinflammatory Panel 1 (human) Controls 1, 2, and 3 are prepared by spiking recombinant calibrators into a non-human serum matrix. The controls are supplied as lyophilized powder.

Storage and Handling

To maximize the consistency of measured values across vials, the controls must be stored at the temperature recommended above. Once reconstituted in 250 µL of Diluent 2, the controls can be stored up to 10 days at 2–8°C. For long term storage, reconstituted controls must be stored frozen at ≤-70°C. Reconstituted controls can go through 3 freeze–thaw cycles without significantly changing analytes levels. Discard unused reconstituted material after the third freeze–thaw cycle.

To use, first reconstitute each vial of Proinflammatory Panel 1 (human) controls in 250 µL of Diluent 2, and then dilute to match the sample dilution (2-fold is the recommended sample dilution for this kit). Add diluted control solutions directly to the MSD Proinflammatory Panel 1 (human) plate, and assay as unknown samples. Discard excess diluted control material after use.

Safety

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, which can be obtained from MSD Customer Service.

Assignment of Control Values

The controls are provided to assess reproducibility of assay performance, and precision CVs are expected to be less than 25%. The certificate of analysis contains the concentrations of the controls measured at MSD across three lots. Even with good laboratory practices, site-to-site differences may occur; therefore, to establish accuracy specifications, it is recommended that each lab should establish its own nominal values and acceptance range for the controls concentrations.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, V-PLEX, www.mesoscale.com, MSD (design), 96 WELL 10-SPOT (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.
©2013, 2021 Meso Scale Diagnostics, LLC. All rights reserved.