

MSD[®] TH17 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 No.	C4085-1
Contents	TH17 Panel 1 (human) Control 1 (5 vials, lyophilized*) TH17 Panel 1 (human) Control 2 (5 vials, lyophilized*) TH17 Panel 1 (human) Control 3 (5 vials, lyophilized*) *Supplied as part of a V-PLEX [®] Plus kit. Quantity will match number of plates ordered.
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 TH17 Panel 1 (human) Control Pack consists of three levels of controls, each containing known concentrations of all human cytokines that are detected by the V-PLEX Plus and V-PLEX TH17 Panel 1 (human) Kits (catalog numbers K15085G and K15085D). TH17 Panel 1 (human) Controls 1, 2, and 3 are prepared by spiking recombinant calibrators into a nonmatrix-based diluent. The controls are supplied as a lyophilized powder.

Storage and Handling

To maximize the consistency of measured values across vials, the controls must be stored at the recommended temperature. Once reconstituted in 250 µL of Diluent 43 (catalog number R50AG), the controls may be stored frozen at ≤−70 °C. Reconstituted controls can go through three freeze-thaw cycles without significantly changing analyte levels. Discard unused reconstituted material after the third freeze-thaw cycle.

To use, first reconstitute each vial of TH17 Panel 1 (human) controls in 250 µL of Diluent 43 and then dilute to match the sample dilution (4-fold is the recommended sample dilution for this kit). Add diluted control solutions directly to the MSD TH17 Panel 1 (human) plate, and assay as unknown samples. Discard excess diluted control material after use.

Assignment of Control Values

The controls are provided to assess the reproducibility of assay performance, and precision CVs are expected to be less than 25%. For the lot-specific concentration of each analyte in the control pack, refer to the certificate of analysis supplied with the kit or available at www.mesoscale.com. 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 control concentrations.

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 (SDS), which can be obtained from MSD Customer Service or at www.mesoscale.com.

MESO SCALE DISCOVERY, 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.
©2020 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-424-v2-2020May