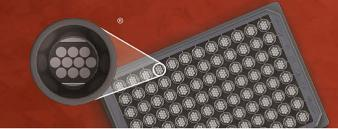
MSD® TH17 Panel 1 (mouse) Calibrator Blend



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 #	C0246-2
Contents	TH17 Panel 1 (mouse) Calibrator Rlend (1 vial: Ivonhilized)

Storage 2-8°C

Summary and Intended Use

The TH17 Panel 1 (mouse) Calibrator Blend is a lyophilized blend of ten recombinant mouse proteins: IL-16, IL-17A, IL-17C, IL-17E/IL-25, IL-17F, IL-21, IL-22, IL-23, IL-31, and MIP-3 α . TH17 Panel 1 (mouse) Calibrator Blend was developed for use as an assay calibrator for MSD assays in the V-PLEX® TH17 Panel 1 (mouse) and the V-PLEX TH17 Panel 1 (mouse) Plus Kits (catalog # K15246G). The utility of this product has not been evaluated in other applications.

Reagent

TH17 Panel 1 (mouse) Calibrator Blend includes full-length recombinant proteins expressed by Sf21 insect cells (IL-23) and E.coli (IL-16, IL-17A, IL-17C, IL-17F, IL-25, IL-17F, IL-21, IL-22, IL-31, and MIP-3 α). TH17 Panel 1 (mouse) Calibrator Blend is supplied as a lyophilized powder that yields top-of-curve calibrator concentration upon reconstitution in 1 mL of assay diluent (Diluent 41; catalog # R50AH).

Storage and Handling

In order to maximize consistency in calibrator values across assay runs, the calibrator must be stored at the recommended temperature and handled per the instructions provided in this document. Once reconstituted in 1 mL of Diluent 41, the calibrator blend is stable at \leq -70°C and can tolerate three freeze—thaw cycles.

TH17 Panel 1 (mouse) calibration curves are constructed based on the concentration of calibrator in each well. Refer to the assay-specific product insert for calibrator preparation instructions. Excess diluted calibrator should be discarded after use.

For the lot-specific concentration of each calibrator in the blend, refer to the certificate of analysis (COA) supplied with the kit or available at www.mesoscale.com.

Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling calibrators. 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, 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.

©2018 Meso Scale Diagnostics, LLC. All rights reserved.



