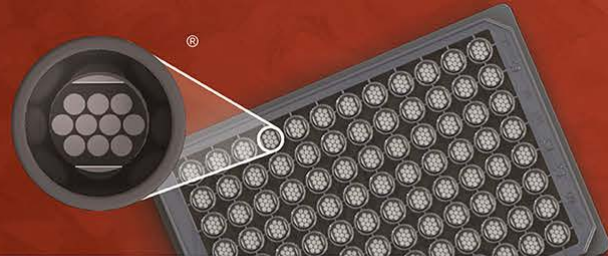


# MSD<sup>®</sup> TH17 Panel 1 (human) Calibrator Blend



## Ordering Information

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

## Scientific Support

Phone: 1-240-314-2798  
Email: ScientificSupport@mesoscale.com  
[www.mesoscale.com/support](http://www.mesoscale.com/support)

## Company Address

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

[www.mesoscale.com](http://www.mesoscale.com)<sup>®</sup>

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

<b>Catalog #</b>	C0085-2
<b>Contents</b>	TH17 Panel 1 (human) Calibrator Blend (1 vial; lyophilized)
<b>Storage</b>	2-8°C

## Summary and Intended Use

The TH17 Panel 1 (human) Calibrator Blend is a lyophilized blend of seven recombinant human proteins: IL-17A, IL-21, IL-22, IL-23, IL-27, IL-31, and MIP-3 $\alpha$ . TH17 Panel 1 (human) Calibrator Blend was developed for use as an assay calibrator for MSD assays in the V-PLEX<sup>®</sup> Plus and the V-PLEX TH17 Panel 1 (human) Kits (catalog # K15085G, K15085D). The utility of this product has not been evaluated in other applications.

## Reagent

TH17 Panel 1 (human) Calibrator Blend includes full-length recombinant proteins expressed by *Sf21* insect cells (IL-23); NSO cells (IL-27); and *E.coli* (IL-17A, IL-21, IL-22, IL-31, and MIP-3 $\alpha$ ). TH17 Panel 1 (human) Calibrator Blend is supplied as a lyophilized powder that yields top-of-curve calibrator concentration upon reconstitution in 1 mL of assay diluent (Diluent 43; catalog # R50AG).

## 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 assay diluent (catalog # R50AG), the calibrator blend is stable at  $\leq -70^{\circ}\text{C}$  and can tolerate three freeze-thaw cycles.

TH17 Panel 1 (human) 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](http://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](http://www.mesoscale.com).

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, V-PLEX, [www.mesoscale.com](http://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. ©2017 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-423-v1-2017Feb