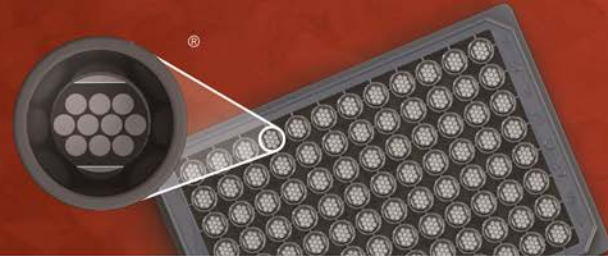


MSD® Proinflammatory Panel 1 (mouse)

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 No	C0048-2
Contents	Proinflammatory Panel 1 (mouse) Calibrator Blend (1 vial; lyophilized)
Storage	2–8°C

Summary and Intended Use

The Proinflammatory Panel 1 (mouse) Calibrator Blend is a lyophilized blend of 10 recombinant mouse proteins: IFN- γ , IL-1 β , IL-2, IL-4, IL-5, IL-6, KC/GRO, IL-10, IL-12p70, and TNF- α . Proinflammatory Panel 1 (mouse) Calibrator Blend was developed for use as an assay calibrator for MSD assays in the V-PLEX™ Proinflammatory Panel 1 (mouse) Kit (catalog # K15048). The utility of this product has not been evaluated in other applications.

Reagent

Proinflammatory Panel 1 (mouse) Calibrator Blend includes full-length *E. coli* recombinant proteins (IFN- γ , IL-1 β , IL-2, IL-4, IL-6, KC/GRO, IL-10, and TNF- α), as well as full-length Sf21 insect cell-expressed recombinant proteins (IL-5 and IL-12p70). Proinflammatory 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 (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 assay diluent (catalog # R50AH), the calibrator blend is stable for 30 days at 2–8°C and can tolerate 5 freeze–thaw cycles.

Proinflammatory 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.

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. Refer to the material safety data sheet, included in the product package and available from MSD Customer Service, for additional product safety and handling information.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, 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 Meso Scale Diagnostics, LLC. All rights reserved.