

MSD® Proinflammatory Panel 1 (rat) 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 #	C0044-2
Contents	Proinflammatory Panel 1 (rat) Calibrator Blend (1 vial; lyophilized)
Storage	2-8°C

Summary and Intended Use

The Proinflammatory Panel 1 (rat) Calibrator Blend is a lyophilized blend of nine recombinant rat proteins: IFN- γ , IL-1 β , IL-4, IL-5, IL-6, KC/GRO, IL-10, IL-13, and TNF- α . Proinflammatory Panel 1 (rat) Calibrator Blend was developed for use as an assay calibrator for MSD assays in the V-PLEX® and V-PLEX Plus Proinflammatory Panel 2 (rat) Kits (catalog # K15059D, K15059G). The utility of this product has not been evaluated in other applications.

Reagent

Proinflammatory Panel 1 (rat) Calibrator Blend includes full-length *E. coli* recombinant proteins (IFN- γ , IL-1 β , IL-4, IL-5, KC/GRO, IL-10, IL-13, and TNF- α) as well as full-length insect cell-expressed recombinant protein (IL-6). Proinflammatory Panel 1 (rat) Calibrator Blend is supplied as a lyophilized powder that yields top-of-curve calibrator concentration upon reconstitution in 1 mL of assay diluent (Diluent 42; catalog # R50AK).

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 42, the calibrator blend is stable at $\leq -70^{\circ}\text{C}$ and can tolerate five freeze-thaw cycles.

Proinflammatory Panel 1 (rat) 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.