

Chemokine Panel 1 (human) Gen. B Calibrator Blend



www.mesoscale.com[®]

Catalog No.	C0705-2
Contents	Chemokine Panel 1 (human) Gen. B Calibrator Blend (1 vial; lyophilized)
Storage	2–8 °C

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: 1-301-947-2025
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

Summary and Intended Use

The Chemokine Panel 1 (human) Gen. B Calibrator Blend is a lyophilized blend of nine recombinant human proteins: Eotaxin, Eotaxin-3, IP-10, MCP-1, MCP-4, MDC, MIP-1 α , MIP-1 β , and TARC. Chemokine Panel 1 (human) Gen. B Calibrator Blend was developed for use as an assay calibrator for assays in the V-PLEX[®] and V-PLEX Plus Chemokine Panel 1 (human) Gen. B (Catalog Nos. K15705D and K15705G), V-PLEX and V-PLEX Plus Chemokine Panel 1 (NHP) Gen. B (Catalog Nos. K15712D and K15712G), and V-PLEX and V-PLEX Plus Gen. B singleplex assays (Eotaxin, Eotaxin-3, IP-10, MCP-1, MCP-4, MDC, MIP-1 α , MIP-1 β , and TARC) that are part of the Chemokine Panel 1 (human) and (NHP) Gen. B Kits. The utility of this product has not been evaluated in other products or applications.

Reagent

Chemokine Panel 1 (human) Gen. B Calibrator Blend is supplied as a lyophilized powder that yields top-of-curve calibrator concentration upon reconstitution in 1 mL of assay diluent (Diluent 57, Catalog No. R50BZ).

Storage and Handling

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 assay diluent, the calibrator blend is stable at ≤ -70 °C and can tolerate 3 freeze-thaw cycles.

Chemokine Panel 1 (human) Gen. B 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. Additional product-specific safety information is available in the safety data sheet (SDS), which can be obtained from MSD Customer Service.

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

MESO SCALE DISCOVERY, Meso Scale Diagnostics, MSD, www.mesoscale.com, MSD (design), V-PLEX, 96 WELL 10-SPOT (design), and Spot the Difference are trademarks and/or service marks of Meso Scale Diagnostics, LLC.
©2023 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-1317-v1-2024Feb