

Chemokine Panel 1 (human) Gen. B Control Pack



www.mesoscale.com®

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

Catalog No.	C4705-1
Contents	Chemokine Panel 1 (human) Gen. B Control 1* (5 vial; lyophilized) Chemokine Panel 1 (human) Gen. B Control 2* (5 vial; lyophilized) Chemokine Panel 1 (human) Gen. B Control 3* (5 vial; lyophilized) *Supplied as part of a V-PLEX Plus kit, the quantity will match the number of plates ordered.
Storage	2–8 °C

Summary and Intended Use

The use of controls to monitor the analytical performance and reliability of test methods is an essential component of good laboratory practice.

The Chemokine Panel 1 (human) Gen. B Control Pack consists of 3 levels of controls, each containing known concentrations of all human chemokine proteins that are detected by 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 Gen. B (NHP) (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 controls span the linear ranges and are designed to test the analytical range of the assays.

Reagent

Chemokine Panel 1 (human) Gen. B Controls 1, 2, and 3 are prepared by spiking recombinant calibrators into a nonmatrix-based diluent. The controls are supplied as a lyophilized powder.

To use, first reconstitute each vial of Chemokine Panel 1 (human) Gen. B controls in 250 μ L of assay diluent (Diluent 57, Catalog No. R50BZ), and then dilute to match the sample dilution (4-fold is the recommended sample dilution for this kit). Add diluted control solutions directly to the MSD® Chemokine Panel 1 (human) Gen. B plate and assay as unknown samples. Discard unused diluted control material.

Storage and Handling

To maximize the consistency of measured values across vials, the controls must be stored at the recommended temperature. Once reconstituted, the controls must be stored frozen at ≤ -70 °C. Reconstituted controls can go through 3 freeze-thaw cycles without significantly changing the level of the analyte. Discard unused reconstituted material after the third freeze-thaw cycle.

Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling controls. 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, which can be obtained from MSD Customer Service.

Assignment of Control Values

The controls are provided to assess reproducibility of assay performance, and precision CVs are expected to be less than 25%. The certificate of analysis contains the concentrations of the controls measured at MSD across three lots. Even with good laboratory practices, site-to-site differences may occur; therefore, to establish accuracy specifications, it is recommended that each lab should establish its own nominal values and acceptance range for the control concentrations.

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-1318-v1-2024Feb