

**CERTIFICATE OF ANALYSIS**

**Product Description:** Chemokine Panel 1 (human) Gen. B Calibrator Blend  
**Catalog Number:** C0705-2  
**Lot Number:** A0080372  
**Expiration Date:** 31 JUL 2028  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1450	pg/mL
MIP-1 $\beta$	594	pg/mL
Eotaxin-3	16900	pg/mL
TARC/CCL17	721	pg/mL
IP-10	1940	pg/mL
MIP-1 $\alpha$	618	pg/mL
MCP-1	397	pg/mL
MDC	4650	pg/mL
MCP-4	377	pg/mL

**Reconstitution Volume:** 1000  $\mu$ L

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		30 AUG 2023