



## CERTIFICATE OF ANALYSIS

**Product Description:** Chemokine Panel 1 (human) Gen. B Calibrator Blend  
**Catalog Number:** C0705-2  
**Lot Number:** A0080389  
**Expiration Date:** 30 APR 2029  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1480	pg/mL
MIP-1 $\beta$	612	pg/mL
Eotaxin-3	15700	pg/mL
TARC/CCL17	778	pg/mL
IP-10	1800	pg/mL
MIP-1 $\alpha$	633	pg/mL
MCP-1	387	pg/mL
MDC	4710	pg/mL
MCP-4	374	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		23 MAY 2024