

**CERTIFICATE OF ANALYSIS**


**Product Description:** Chemokine Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0047-2  
**Lot Number:** A0080418  
**Expiration Date:** 28 FEB 2030  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1810	pg/mL
MIP-1 $\beta$	954	pg/mL
Eotaxin-3	5770	pg/mL
TARC/CCL17	1730	pg/mL
IP-10	2470	pg/mL
MIP-1 $\alpha$	1180	pg/mL
IL-8 (HA)	73600	pg/mL
MCP-1	546	pg/mL
MDC	5630	pg/mL
MCP-4	748	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		22 APR 2025