

**CERTIFICATE OF ANALYSIS**


Product Description: Chemokine Panel 1 (human) Calibrator Blend
Catalog Number: C0047-2
Lot Number: A0080456
Expiration Date: 31 DEC 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1860	pg/mL
MIP-1 β	889	pg/mL
Eotaxin-3	5620	pg/mL
TARC/CCL17	1780	pg/mL
IP-10	2310	pg/mL
MIP-1 α	1180	pg/mL
IL-8 (HA)	75200	pg/mL
MCP-1	580	pg/mL
MDC	5560	pg/mL
MCP-4	744	pg/mL

Reconstitution Volume: 1000 μ 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
Review/Approval	Karen Hamilla	Quality		03 FEB 2026