



CERTIFICATE OF ANALYSIS

Product Description: Chemokine Panel 1 (human) Calibrator Blend
Catalog Number: C0047-2
Lot Number: A0080432
Expiration Date: 30 APR 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1890	pg/mL
MIP-1β	890	pg/mL
Eotaxin-3	6330	pg/mL
TARC/CCL17	1780	pg/mL
IP-10	2520	pg/mL
MIP-1α	1180	pg/mL
IL-8 (HA)	76100	pg/mL
MCP-1	519	pg/mL
MDC	5600	pg/mL
MCP-4	746	pg/mL

Reconstitution Volume: 1000 µL

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		21 MAY 2025