

**CERTIFICATE OF ANALYSIS**

Product Description: Chemokine Panel 1 (human) Calibrator Blend
Catalog Number: C0047-2
Lot Number: A0080347
Expiration Date: 31 OCT 2027
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1820	pg/mL
MIP-1 β	1160	pg/mL
Eotaxin-3	6120	pg/mL
TARC/CCL17	1820	pg/mL
IP-10	2870	pg/mL
MIP-1 α	1160	pg/mL
IL-8 (HA)	72000	pg/mL
MCP-1	545	pg/mL
MDC	6030	pg/mL
MCP-4	703	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		01 DEC 2022