



CERTIFICATE OF ANALYSIS

Product Description: Chemokine Panel 1 (human) Calibrator Blend
Catalog Number: C0047-2
Lot Number: A0080388
Expiration Date: 31 MAR 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1820	pg/mL
MIP-1 β	1170	pg/mL
Eotaxin-3	5790	pg/mL
TARC/CCL17	1750	pg/mL
IP-10	2320	pg/mL
MIP-1 α	1170	pg/mL
IL-8 (HA)	70900	pg/mL
MCP-1	553	pg/mL
MDC	5870	pg/mL
MCP-4	758	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		04 JUN 2024