



CERTIFICATE OF ANALYSIS

Product Description: Chemokine Panel 1 (human) Gen. B Calibrator Blend
Catalog Number: C0705-2
Lot Number: A0080416
Expiration Date: 31 DEC 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1450	pg/mL
MIP-1 β	583	pg/mL
Eotaxin-3	21500	pg/mL
TARC/CCL17	744	pg/mL
IP-10	1860	pg/mL
MIP-1 α	618	pg/mL
MCP-1	379	pg/mL
MDC	4620	pg/mL
MCP-4	358	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		31 JAN 2025