



CERTIFICATE OF ANALYSIS

Product Description: Chemokine Panel 1 (human) Gen. B Calibrator Blend
Catalog Number: C0705-2
Lot Number: A0080462
Expiration Date: 31 MAR 2031
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1500	pg/mL
MIP-1 β	668	pg/mL
Eotaxin-3	17800	pg/mL
TARC/CCL17	784	pg/mL
IP-10	1970	pg/mL
MIP-1 α	683	pg/mL
MCP-1	413	pg/mL
MDC	4740	pg/mL
MCP-4	365	pg/mL

Reconstitution Volume:

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		24 APR 2026