



CERTIFICATE OF ANALYSIS

Product Description: Chemokine Panel 1 (human) Gen. B Calibrator Blend
Catalog Number: C0705-2
Lot Number: A0080377
Expiration Date: 31 AUG 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1480	pg/mL
MIP-1 β	606	pg/mL
Eotaxin-3	17900	pg/mL
TARC/CCL17	744	pg/mL
IP-10	1920	pg/mL
MIP-1 α	614	pg/mL
MCP-1	395	pg/mL
MDC	4710	pg/mL
MCP-4	383	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		27 SEP 2023