

**CERTIFICATE OF ANALYSIS**


Product Description: Chemokine Panel 1 (human) Gen. B Calibrator Blend
Catalog Number: C0705-2
Lot Number: A0080431
Expiration Date: 31 MAR 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1430	pg/mL
MIP-1 β	633	pg/mL
Eotaxin-3	17400	pg/mL
TARC/CCL17	730	pg/mL
IP-10	1980	pg/mL
MIP-1 α	590	pg/mL
MCP-1	370	pg/mL
MDC	4500	pg/mL
MCP-4	353	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		29 MAY 2025