



CERTIFICATE OF ANALYSIS

Product Description: Chemokine Panel 1 (human) Gen. B Calibrator Blend
Catalog Number: C0705-2
Lot Number: A0080370
Expiration Date: 30 JUN 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1420	pg/mL
MIP-1β	578	pg/mL
Eotaxin-3	16500	pg/mL
TARC/CCL17	746	pg/mL
IP-10	1780	pg/mL
MIP-1α	602	pg/mL
MCP-1	398	pg/mL
MDC	4700	pg/mL
MCP-4	392	pg/mL

Reconstitution Volume: 1000 µL

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	<i>Karen Hamilla</i>	28 AUG 2023