



CERTIFICATE OF ANALYSIS

Product Description: Chemokine Panel 1 (human) Calibrator Blend
Catalog Number: C0047-2
Lot Number: A0080338
Expiration Date: 31 AUG 2027
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1770	pg/mL
MIP-1β	1010	pg/mL
Eotaxin-3	6760	pg/mL
TARC/CCL17	1750	pg/mL
IP-10	2850	pg/mL
MIP-1α	1180	pg/mL
IL-8 (HA)	75400	pg/mL
MCP-1	564	pg/mL
MDC	5740	pg/mL
MCP-4	720	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		08 NOV 2022