



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Chemokine Panel 1 (human) Calibrator Blend

Catalog Number: C0047-2

Lot Number: A0080266

Expiration Date: 31 AUG 2025

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1885	pg/mL
MIP-1 β	1190	pg/mL
Eotaxin-3	5405	pg/mL
TARC	1845	pg/mL
IP-10	2360	pg/mL
MIP-1 α	1160	pg/mL
IL-8	76850	pg/mL
MCP-1	551.5	pg/mL
MDC	10150	pg/mL
MCP-4	738	pg/mL

Reconstitution Volume: 1000 μ L

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		05 NOV 2020