



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Chemokine Panel 1 (human) Calibrator Blend

Catalog Number: C0047-2

Lot Number: A0080280

Expiration Date: 28 FEB 2026

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1890	pg/mL
MIP-1 β	905	pg/mL
Eotaxin-3	5550	pg/mL
TARC	1860	pg/mL
IP-10	2520	pg/mL
MIP-1 α	1090	pg/mL
IL-8	68900	pg/mL
MCP-1	544	pg/mL
MDC	6080	pg/mL
MCP-4	691	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		28 APR 2021