



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Chemokine Panel 1 (human) Calibrator Blend

Catalog Number: C0047-2

Lot Number: A0080320

Expiration Date: 31 MAR 2027

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1960	pg/mL
MIP-1 β	1150	pg/mL
Eotaxin-3	6460	pg/mL
TARC	1990	pg/mL
IP-10	2510	pg/mL
MIP-1 α	1160	pg/mL
IL-8	70600	pg/mL
MCP-1	558	pg/mL
MDC	5880	pg/mL
MCP-4	734	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		24 JUN 2022