



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Chemokine Panel 1 (human) Calibrator Blend

Catalog Number: C0047-2

Lot Number: A0080259

Expiration Date: 31 MAY 2025

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1880	pg/mL
MIP-1 β	1270	pg/mL
Eotaxin-3	4990	pg/mL
TARC	1800	pg/mL
IP-10	2340	pg/mL
MIP-1 α	1100	pg/mL
IL-8	78600	pg/mL
MCP-1	553	pg/mL
MDC	9730	pg/mL
MCP-4	705	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		22 JUN 2020