



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Chemokine Panel 1 (human) Calibrator Blend

**Catalog Number:** C0047-2

**Lot Number:** A0080306

**Expiration Date:** 30 SEP 2026

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1945	pg/mL
MIP-1 $\beta$	1145	pg/mL
Eotaxin-3	5320	pg/mL
TARC	2080	pg/mL
IP-10	2555	pg/mL
MIP-1 $\alpha$	1110	pg/mL
IL-8	72900	pg/mL
MCP-1	546	pg/mL
MDC	6025	pg/mL
MCP-4	688	pg/mL

**Reconstitution Volume:** 1000  $\mu$ L

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		15 NOV 2021