



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0047-2

**Lot Number:** A0080228

**Expiration Date:** 31 AUG 2024

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	1870	pg/mL
MIP-1 $\beta$	1330	pg/mL
Eotaxin-3	5970	pg/mL
TARC	1830	pg/mL
IP-10	2360	pg/mL
MIP-1 $\alpha$	1060	pg/mL
IL-8	64500	pg/mL
MCP-1	552	pg/mL
MDC	10200	pg/mL
MCP-4	724	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
Review/Approval	Karen Hamilla	Quality		20 NOV 2019