



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0050-2

**Lot Number:** A0080331

**Expiration Date:** 31 JUL 2027

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	1180	pg/mL
IL-1 $\alpha$	422	pg/mL
IL-5	888	pg/mL
IL-7	913	pg/mL
IL-12/IL-23p40	3450	pg/mL
IL-15	840	pg/mL
IL-16	2620	pg/mL
IL-17A	6010	pg/mL
TNF- $\beta$	717	pg/mL
VEGF	1280	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		26 SEP 2022