



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0050-2

**Lot Number:** A0080240

**Expiration Date:** 30 NOV 2024

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	1260	pg/mL
IL-1 $\alpha$	320	pg/mL
IL-5	857	pg/mL
IL-7	807	pg/mL
IL-12/IL-23p40	3380	pg/mL
IL-15	941	pg/mL
IL-16	2910	pg/mL
IL-17A	6040	pg/mL
TNF- $\beta$	623	pg/mL
VEGF	1350	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>	Carissa Hubbard	Quality		23 DEC 2019