



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (human) Calibrator Blend

Catalog Number: C0050-2

Lot Number: A0080253

Expiration Date: 30 APR 2025

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	1280	pg/mL
IL-1 α	464	pg/mL
IL-5	906	pg/mL
IL-7	826	pg/mL
IL-12/IL-23p40	3650	pg/mL
IL-15	863	pg/mL
IL-16	3150	pg/mL
IL-17A	5340	pg/mL
TNF- β	705	pg/mL
VEGF	1520	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		01 JUL 2020