



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (human) Calibrator Blend

Catalog Number: C0050-2

Lot Number: A0080283

Expiration Date: 31 MAR 2026

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	1200	pg/mL
IL-1 α	438	pg/mL
IL-5	853	pg/mL
IL-7	766	pg/mL
IL-12/IL-23p40	3270	pg/mL
IL-15	714	pg/mL
IL-16	2930	pg/mL
IL-17A	6710	pg/mL
TNF- β	792	pg/mL
VEGF	1150	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		05 MAY 2021