



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (human) Calibrator Blend

Catalog Number: C0050-2

Lot Number: A0080307

Expiration Date: 30 SEP 2026

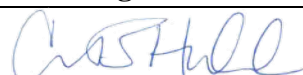
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	1250	pg/mL
IL-1 α	448	pg/mL
IL-5	849	pg/mL
IL-7	829	pg/mL
IL-12/IL-23p40	3220	pg/mL
IL-15	810	pg/mL
IL-16	2700	pg/mL
IL-17A	6970	pg/mL
TNF- β	761	pg/mL
VEGF	1140	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	Carissa B. Hubbard	Quality		21 DEC 2021