



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0235

Expiration Date: 30 SEP 2026

Storage Temperature: 2-8°C

Intended Assay Species: Human

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	36750	pg	147000	pg/mL
IL-1 β	5738	pg	22950	pg/mL
IL-2	2600	pg	10400	pg/mL
IL-4	2838	pg	11350	pg/mL
IL-6	2625	pg	10500	pg/mL
IL-8	2988	pg	11950	pg/mL
IL-10	5413	pg	21650	pg/mL
IL-12p70	7825	pg	31300	pg/mL
IL-13	2375	pg	9500	pg/mL
TNF- α	4188	pg	16750	pg/mL
IL-5	5825	pg	23300	pg/mL
GM-CSF	13750	pg	55000	pg/mL
IL-17A	31375	pg	125500	pg/mL
VEGF-A	4138	pg	16550	pg/mL

Reconstitution Volume: 250 μ L

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 DEC 2021