



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060 -2

Lot Number: A00U0227

Expiration Date: 31 AUG 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	33750	pg	135000	pg/mL
IL-1 β	5538	pg	22150	pg/mL
IL-2	2488	pg	9950	pg/mL
IL-4	2713	pg	10850	pg/mL
IL-6	2563	pg	10250	pg/mL
IL-8	2863	pg	11450	pg/mL
IL-10	5150	pg	20600	pg/mL
IL-12p70	7500	pg	30000	pg/mL
IL-13	2350	pg	9400	pg/mL
TNF- α	4275	pg	17100	pg/mL
IL-5	5638	pg	22550	pg/mL
GM-CSF	13500	pg	54000	pg/mL
IL-17A	31250	pg	125000	pg/mL
VEGF-A	4125	pg	16500	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		28 OCT 2021