



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0252

Expiration Date: 28 FEB 2027

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	32250	pg	129000	pg/mL
IL-1 β	5638	pg	22550	pg/mL
IL-2	2538	pg	10150	pg/mL
IL-4	2263	pg	9050	pg/mL
IL-6	2375	pg	9500	pg/mL
IL-8	2763	pg	11050	pg/mL
IL-10	4925	pg	19700	pg/mL
IL-12p70	7213	pg	28850	pg/mL
IL-13	2088	pg	8350	pg/mL
TNF- α	3625	pg	14500	pg/mL
IL-5	5525	pg	22100	pg/mL
GM-CSF	13500	pg	54000	pg/mL
IL-17A	33250	pg	133000	pg/mL
VEGF-A	4038	pg	16150	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		19 APR 2022