



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0274

Expiration Date: 31 JUL 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	38375	pg	153500	pg/mL
IL-1 β	5913	pg	23650	pg/mL
IL-2	2513	pg	10050	pg/mL
IL-4	2188	pg	8750	pg/mL
IL-6	2463	pg	9850	pg/mL
IL-8	3125	pg	12500	pg/mL
IL-10	4975	pg	19900	pg/mL
IL-12p70	7150	pg	28600	pg/mL
IL-13	2188	pg	8750	pg/mL
TNF- α	3900	pg	15600	pg/mL
IL-5	5650	pg	22600	pg/mL
GM-CSF	13500	pg	54000	pg/mL
IL-17A	36375	pg	145500	pg/mL
VEGF-A	3163	pg	12650	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		30 SEP 2022