



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0206

Expiration Date: 31 MAR 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	34250	pg	137000	pg/mL
IL-1 β	5650	pg	22600	pg/mL
IL-2	2613	pg	10450	pg/mL
IL-4	2500	pg	10000	pg/mL
IL-6	2575	pg	10300	pg/mL
IL-8	2938	pg	11750	pg/mL
IL-10	5350	pg	21400	pg/mL
IL-12p70	7875	pg	31500	pg/mL
IL-13	2613	pg	10450	pg/mL
TNF- α	4600	pg	18400	pg/mL
IL-5	5750	pg	23000	pg/mL
GM-CSF	13875	pg	55500	pg/mL
IL-17A	34625	pg	138500	pg/mL
VEGF-A	4425	pg	17700	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		03 JUN 2021