



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0268

Expiration Date: 30 APR 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	39250	pg	157000	pg/mL
IL-1 β	5663	pg	22650	pg/mL
IL-2	2550	pg	10200	pg/mL
IL-4	2200	pg	8800	pg/mL
IL-6	2313	pg	9250	pg/mL
IL-8	2688	pg	10750	pg/mL
IL-10	4775	pg	19100	pg/mL
IL-12p70	6788	pg	27150	pg/mL
IL-13	2013	pg	8050	pg/mL
TNF- α	3625	pg	14500	pg/mL
IL-5	5388	pg	21550	pg/mL
GM-CSF	13250	pg	53000	pg/mL
IL-17A	36875	pg	147500	pg/mL
VEGF-A	4075	pg	16300	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		15 JUN 2022