



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0276

Expiration Date: 31 AUG 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	35000	pg	140000	pg/mL
IL-1 β	5700	pg	22800	pg/mL
IL-2	2438	pg	9750	pg/mL
IL-4	2488	pg	9950	pg/mL
IL-6	2300	pg	9200	pg/mL
IL-8	2925	pg	11700	pg/mL
IL-10	4663	pg	18650	pg/mL
IL-12p70	6825	pg	27300	pg/mL
IL-13	2175	pg	8700	pg/mL
TNF- α	3600	pg	14400	pg/mL
IL-5	5388	pg	21550	pg/mL
GM-CSF	11913	pg	47650	pg/mL
IL-17A	38375	pg	153500	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