



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0368
Expiration Date: 31 MAR 2029
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	37500	pg	150000	pg/mL
IL-1 β	5425	pg	21700	pg/mL
IL-2	2575	pg	10300	pg/mL
IL-4	2150	pg	8600	pg/mL
IL-6	2600	pg	10400	pg/mL
IL-8	2825	pg	11300	pg/mL
IL-10	5012.5	pg	20050	pg/mL
IL-12p70	7450	pg	29800	pg/mL
IL-13	2312.5	pg	9250	pg/mL
TNF- α	3737.5	pg	14950	pg/mL
IL-5	5500	pg	22000	pg/mL
GM-CSF	12875	pg	51500	pg/mL
IL-17A	25250	pg	101000	pg/mL
VEGF-A	4062.5	pg	16250	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 MAY 2024