



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0462
Expiration Date: 31 DEC 2030
Storage Temperature: 2-8°C
Intended Assay Species: Human, NHP

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	23550	pg	94200	pg/mL
IL-1 β	4925	pg	19700	pg/mL
IL-2	2700	pg	10800	pg/mL
IL-4	2485	pg	9940	pg/mL
IL-6	2342.5	pg	9370	pg/mL
IL-8	2500	pg	10000	pg/mL
IL-10	5400	pg	21600	pg/mL
IL-12p70	7375	pg	29500	pg/mL
IL-13	2387.5	pg	9550	pg/mL
TNF- α	4475	pg	17900	pg/mL
IL-5	5225	pg	20900	pg/mL
GM-CSF	13075	pg	52300	pg/mL
IL-17A	25250	pg	101000	pg/mL
VEGF-A	4675	pg	18700	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		19 FEB 2026