



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0410
Expiration Date: 28 FEB 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	26250	pg	105000	pg/mL
IL-1 β	5012.5	pg	20050	pg/mL
IL-2	2387.5	pg	9550	pg/mL
IL-4	2712.5	pg	10850	pg/mL
IL-6	2450	pg	9800	pg/mL
IL-8	2575	pg	10300	pg/mL
IL-10	5225	pg	20900	pg/mL
IL-12p70	7750	pg	31000	pg/mL
IL-13	2625	pg	10500	pg/mL
TNF- α	4375	pg	17500	pg/mL
IL-5	5437.5	pg	21750	pg/mL
GM-CSF	11050	pg	44200	pg/mL
IL-17A	27750	pg	111000	pg/mL
VEGF-A	5087.5	pg	20350	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		08 APR 2025