



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0411
Expiration Date: 30 APR 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	20625	pg	82500	pg/mL
IL-1 β	4925	pg	19700	pg/mL
IL-2	2687.5	pg	10750	pg/mL
IL-4	2500	pg	10000	pg/mL
IL-6	2337.5	pg	9350	pg/mL
IL-8	2387.5	pg	9550	pg/mL
IL-10	4925	pg	19700	pg/mL
IL-12p70	7325	pg	29300	pg/mL
IL-13	2550	pg	10200	pg/mL
TNF- α	4250	pg	17000	pg/mL
IL-5	5012.5	pg	20050	pg/mL
GM-CSF	10887.5	pg	43550	pg/mL
IL-17A	25125	pg	100500	pg/mL
VEGF-A	4700	pg	18800	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		13 JUN 2025