



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0447
Expiration Date: 31 JUL 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	24125	pg	96500	pg/mL
IL-1 β	4875	pg	19500	pg/mL
IL-2	2712.5	pg	10850	pg/mL
IL-4	2512.5	pg	10050	pg/mL
IL-6	2362.5	pg	9450	pg/mL
IL-8	2437.5	pg	9750	pg/mL
IL-10	4900	pg	19600	pg/mL
IL-12p70	7250	pg	29000	pg/mL
IL-13	2337.5	pg	9350	pg/mL
TNF- α	4112.5	pg	16450	pg/mL
IL-5	5050	pg	20200	pg/mL
GM-CSF	10950	pg	43800	pg/mL
IL-17A	27375	pg	109500	pg/mL
VEGF-A	4900	pg	19600	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		17 SEP 2025