



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0453
Expiration Date: 31 AUG 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	23875	pg	95500	pg/mL
IL-1 β	4675	pg	18700	pg/mL
IL-2	2612.5	pg	10450	pg/mL
IL-4	2437.5	pg	9750	pg/mL
IL-6	2337.5	pg	9350	pg/mL
IL-8	2325	pg	9300	pg/mL
IL-10	4737.5	pg	18950	pg/mL
IL-12p70	7112.5	pg	28450	pg/mL
IL-13	2175	pg	8700	pg/mL
TNF- α	4287.5	pg	17150	pg/mL
IL-5	5062.5	pg	20250	pg/mL
GM-CSF	10950	pg	43800	pg/mL
IL-17A	25125	pg	100500	pg/mL
VEGF-A	4725	pg	18900	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		22 SEP 2025