



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0474
Expiration Date: 30 APR 2031
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	25500	pg	102000	pg/mL
IL-1 β	5050	pg	20200	pg/mL
IL-2	2875	pg	11500	pg/mL
IL-4	2385	pg	9540	pg/mL
IL-6	2395	pg	9580	pg/mL
IL-8	2525	pg	10100	pg/mL
IL-10	5350	pg	21400	pg/mL
IL-12p70	7350	pg	29400	pg/mL
IL-13	2425	pg	9700	pg/mL
TNF- α	4425	pg	17700	pg/mL
IL-5	5200	pg	20800	pg/mL
GM-CSF	13325	pg	53300	pg/mL
IL-17A	24425	pg	97700	pg/mL
VEGF-A	4650	pg	18600	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		14 MAY 2026