



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0477
Expiration Date: 31 MAY 2031
Storage Temperature: 2-8°C
Intended Assay Species: Human, NHP

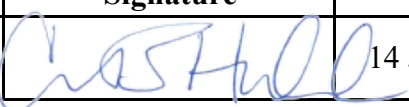
Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	25250	pg	101000	pg/mL
IL-1 β	4725	pg	18900	pg/mL
IL-2	2347.5	pg	9390	pg/mL
IL-4	2387.5	pg	9550	pg/mL
IL-6	2312.5	pg	9250	pg/mL
IL-8	2482.5	pg	9930	pg/mL
IL-10	5350	pg	21400	pg/mL
IL-12p70	7225	pg	28900	pg/mL
IL-13	2287.5	pg	9150	pg/mL
TNF- α	4375	pg	17500	pg/mL
IL-5	5125	pg	20500	pg/mL
GM-CSF	12925	pg	51700	pg/mL
IL-17A	23650	pg	94600	pg/mL
VEGF-A	4550	pg	18200	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	Carissa B. Hubbard	Quality		14 JUL 2026