



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0423
Expiration Date: 30 JUN 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	28750	pg	115000	pg/mL
IL-1 β	5825	pg	23300	pg/mL
IL-2	3037.5	pg	12150	pg/mL
IL-4	3075	pg	12300	pg/mL
IL-6	2662.5	pg	10650	pg/mL
IL-8	2925	pg	11700	pg/mL
IL-10	5637.5	pg	22550	pg/mL
IL-12p70	8187.5	pg	32750	pg/mL
IL-13	2637.5	pg	10550	pg/mL
TNF- α	4825	pg	19300	pg/mL
IL-5	6000	pg	24000	pg/mL
GM-CSF	11925	pg	47700	pg/mL
IL-17A	31625	pg	126500	pg/mL
VEGF-A	5462.5	pg	21850	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		23 JUL 2025