



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0448
Expiration Date: 31 OCT 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	25750	pg	103000	pg/mL
IL-1 β	5075	pg	20300	pg/mL
IL-2	2925	pg	11700	pg/mL
IL-4	2625	pg	10500	pg/mL
IL-6	2550	pg	10200	pg/mL
IL-8	2525	pg	10100	pg/mL
IL-10	5175	pg	20700	pg/mL
IL-12p70	7600	pg	30400	pg/mL
IL-13	2625	pg	10500	pg/mL
TNF- α	4400	pg	17600	pg/mL
IL-5	4775	pg	19100	pg/mL
GM-CSF	10850	pg	43400	pg/mL
IL-17A	24800	pg	99200	pg/mL
VEGF-A	4250	pg	17000	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		06 JAN 2026