



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0376
Expiration Date: 30 JUN 2029
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	40125	pg	160500	pg/mL
IL-1 β	5550	pg	22200	pg/mL
IL-2	2362.5	pg	9450	pg/mL
IL-4	2062.5	pg	8250	pg/mL
IL-6	2600	pg	10400	pg/mL
IL-8	2812.5	pg	11250	pg/mL
IL-10	5037.5	pg	20150	pg/mL
IL-12p70	7212.5	pg	28850	pg/mL
IL-13	2237.5	pg	8950	pg/mL
TNF- α	3387.5	pg	13550	pg/mL
IL-5	5350	pg	21400	pg/mL
GM-CSF	11550	pg	46200	pg/mL
IL-17A	22875	pg	91500	pg/mL
VEGF-A	3237.5	pg	12950	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		17 SEP 2024