



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0356
Expiration Date: 31 DEC 2028
Storage Temperature: 2-8°C
Intended Assay Species: Human

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	39500	pg	158000	pg/mL
IL-1 β	6100	pg	24400	pg/mL
IL-2	2825	pg	11300	pg/mL
IL-4	3062.5	pg	12250	pg/mL
IL-6	3300	pg	13200	pg/mL
IL-8	3250	pg	13000	pg/mL
IL-10	6062.5	pg	24250	pg/mL
IL-12p70	8825	pg	35300	pg/mL
IL-13	2750	pg	11000	pg/mL
TNF- α	4900	pg	19600	pg/mL
IL-5	6300	pg	25200	pg/mL
GM-CSF	14500	pg	58000	pg/mL
IL-17A	26250	pg	105000	pg/mL
VEGF-A	4800	pg	19200	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		02 FEB 2024