



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0363
Expiration Date: 28 FEB 2029
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	39375	pg	157500	pg/mL
IL-1 β	6287.5	pg	25150	pg/mL
IL-2	2650	pg	10600	pg/mL
IL-4	2200	pg	8800	pg/mL
IL-6	2725	pg	10900	pg/mL
IL-8	2762.5	pg	11050	pg/mL
IL-10	5062.5	pg	20250	pg/mL
IL-12p70	7112.5	pg	28450	pg/mL
IL-13	2287.5	pg	9150	pg/mL
TNF- α	3625	pg	14500	pg/mL
IL-5	5462.5	pg	21850	pg/mL
GM-CSF	13125	pg	52500	pg/mL
IL-17A	25000	pg	100000	pg/mL
VEGF-A	3950	pg	15800	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		12 MAR 2024