



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0397
Expiration Date: 31 OCT 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	42625	pg	170500	pg/mL
IL-1 β	5700	pg	22800	pg/mL
IL-2	2362.5	pg	9450	pg/mL
IL-4	2875	pg	11500	pg/mL
IL-6	2612.5	pg	10450	pg/mL
IL-8	3037.5	pg	12150	pg/mL
IL-10	5287.5	pg	21150	pg/mL
IL-12p70	7875	pg	31500	pg/mL
IL-13	2375	pg	9500	pg/mL
TNF- α	5025	pg	20100	pg/mL
IL-5	5825	pg	23300	pg/mL
GM-CSF	11900	pg	47600	pg/mL
IL-17A	28250	pg	113000	pg/mL
VEGF-A	5425	pg	21700	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 Hamila	Quality		04 DEC 2024