



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0149
Expiration Date: 31 MAR 2025
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	35750	pg	143000	pg/mL
IL-1 β	5400	pg	21600	pg/mL
IL-2	2538	pg	10150	pg/mL
IL-4	2788	pg	11150	pg/mL
IL-6	2525	pg	10100	pg/mL
IL-8	2738	pg	10950	pg/mL
IL-10	4625	pg	18500	pg/mL
IL-12p70	7813	pg	31250	pg/mL
IL-13	3175	pg	12700	pg/mL
TNF- α	4525	pg	18100	pg/mL
IL-5	5613	pg	22450	pg/mL
GM-CSF	15500	pg	62000	pg/mL
IL-17A	26250	pg	105000	pg/mL
VEGF-A	8463	pg	33850	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		20 APR 2020