



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0161
Expiration Date: 31 JUL 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	34250	pg	137000	pg/mL
IL-1 β	5263	pg	21050	pg/mL
IL-2	2625	pg	10500	pg/mL
IL-4	2563	pg	10250	pg/mL
IL-6	2638	pg	10550	pg/mL
IL-8	2888	pg	11550	pg/mL
IL-10	4713	pg	18850	pg/mL
IL-12p70	7850	pg	31400	pg/mL
IL-13	3213	pg	12850	pg/mL
TNF- α	4725	pg	18900	pg/mL
IL-5	5588	pg	22350	pg/mL
GM-CSF	15875	pg	63500	pg/mL
IL-17A	30875	pg	123500	pg/mL
VEGF-A	6288	pg	25150	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 AUG 2020