



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0169

Expiration Date: 31 AUG 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	35625	pg	142500	pg/mL
IL-1 β	5050	pg	20200	pg/mL
IL-2	2600	pg	10400	pg/mL
IL-4	2538	pg	10150	pg/mL
IL-6	2338	pg	9350	pg/mL
IL-8	2575	pg	10300	pg/mL
IL-10	4675	pg	18700	pg/mL
IL-12p70	7613	pg	30450	pg/mL
IL-13	2950	pg	11800	pg/mL
TNF- α	4325	pg	17300	pg/mL
IL-5	5488	pg	21950	pg/mL
GM-CSF	16000	pg	64000	pg/mL
IL-17A	31375	pg	125500	pg/mL
VEGF-A	6438	pg	25750	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 OCT 2020