



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0175

Expiration Date: 31 OCT 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	29750	pg	119000	pg/mL
IL-1 β	5013	pg	20050	pg/mL
IL-2	2538	pg	10150	pg/mL
IL-4	2363	pg	9450	pg/mL
IL-6	2538	pg	10150	pg/mL
IL-8	2575	pg	10300	pg/mL
IL-10	4825	pg	19300	pg/mL
IL-12p70	7763	pg	31050	pg/mL
IL-13	2975	pg	11900	pg/mL
TNF- α	4575	pg	18300	pg/mL
IL-5	5163	pg	20650	pg/mL
GM-CSF	14375	pg	57500	pg/mL
IL-17A	30250	pg	121000	pg/mL
VEGF-A	6900	pg	27600	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		14 JAN 2021