



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0181

Expiration Date: 31 DEC 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	31000	pg	124000	pg/mL
IL-1 β	5325	pg	21300	pg/mL
IL-2	2563	pg	10250	pg/mL
IL-4	2375	pg	9500	pg/mL
IL-6	2550	pg	10200	pg/mL
IL-8	2788	pg	11150	pg/mL
IL-10	4988	pg	19950	pg/mL
IL-12p70	7925	pg	31700	pg/mL
IL-13	2800	pg	11200	pg/mL
TNF- α	4600	pg	18400	pg/mL
IL-5	5438	pg	21750	pg/mL
GM-CSF	14750	pg	59000	pg/mL
IL-17A	30625	pg	122500	pg/mL
VEGF-A	6800	pg	27200	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		25 JAN 2021