



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1

**Catalog Number:** C0060-2

**Lot Number:** A00U0213

**Expiration Date:** 31 MAY 2026

**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- $\gamma$	35875	pg	143500	pg/mL
IL-1 $\beta$	5850	pg	23400	pg/mL
IL-2	2650	pg	10600	pg/mL
IL-4	2313	pg	9250	pg/mL
IL-6	2625	pg	10500	pg/mL
IL-8	2988	pg	11950	pg/mL
IL-10	5438	pg	21750	pg/mL
IL-12p70	7850	pg	31400	pg/mL
IL-13	2525	pg	10100	pg/mL
TNF- $\alpha$	4425	pg	17700	pg/mL
IL-5	5663	pg	22650	pg/mL
GM-CSF	13500	pg	54000	pg/mL
IL-17A	34375	pg	137500	pg/mL
VEGF-A	4363	pg	17450	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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		31 AUG 2021