



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1

**Catalog Number:** C0060-2

**Lot Number:** A00U0260

**Expiration Date:** 31 MAR 2027

**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- $\gamma$	31000	pg	124000	pg/mL
IL-1 $\beta$	5613	pg	22450	pg/mL
IL-2	2463	pg	9850	pg/mL
IL-4	2175	pg	8700	pg/mL
IL-6	2388	pg	9550	pg/mL
IL-8	2788	pg	11150	pg/mL
IL-10	4875	pg	19500	pg/mL
IL-12p70	7100	pg	28400	pg/mL
IL-13	2150	pg	8600	pg/mL
TNF- $\alpha$	3613	pg	14450	pg/mL
IL-5	5425	pg	21700	pg/mL
GM-CSF	13250	pg	53000	pg/mL
IL-17A	32875	pg	131500	pg/mL
VEGF-A	3900	pg	15600	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
<b>Review/Approval</b>	Karen Hamilla	Quality		03 MAY 2022