



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1

**Catalog Number:** C0060-2

**Lot Number:** A00U0291

**Expiration Date:** 30 SEP 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- $\gamma$	37125	pg	148500	pg/mL
IL-1 $\beta$	6038	pg	24150	pg/mL
IL-2	2525	pg	10100	pg/mL
IL-4	2475	pg	9900	pg/mL
IL-6	2413	pg	9650	pg/mL
IL-8	3088	pg	12350	pg/mL
IL-10	4950	pg	19800	pg/mL
IL-12p70	7300	pg	29200	pg/mL
IL-13	2163	pg	8650	pg/mL
TNF- $\alpha$	3750	pg	15000	pg/mL
IL-5	5700	pg	22800	pg/mL
GM-CSF	12750	pg	51000	pg/mL
IL-17A	36500	pg	146000	pg/mL
VEGF-A	3338	pg	13350	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		09 NOV 2022