



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1

**Catalog Number:** C0060-2

**Lot Number:** A00U0240

**Expiration Date:** 28 FEB 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- $\gamma$	31250	pg	125000	pg/mL
IL-1 $\beta$	5288	pg	21150	pg/mL
IL-2	2488	pg	9950	pg/mL
IL-4	2225	pg	8900	pg/mL
IL-6	2338	pg	9350	pg/mL
IL-8	2725	pg	10900	pg/mL
IL-10	4713	pg	18850	pg/mL
IL-12p70	6775	pg	27100	pg/mL
IL-13	2075	pg	8300	pg/mL
TNF- $\alpha$	3675	pg	14700	pg/mL
IL-5	5213	pg	20850	pg/mL
GM-CSF	12750	pg	51000	pg/mL
IL-17A	32750	pg	131000	pg/mL
VEGF-A	3363	pg	13450	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		15 MAR 2022