



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1

**Catalog Number:** C0060-2

**Lot Number:** A00U0138

**Expiration Date:** 31 DEC 2024

**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- $\gamma$	37875	pg	151500	pg/mL
IL-1 $\beta$	5463	pg	21850	pg/mL
IL-2	2588	pg	10350	pg/mL
IL-4	2925	pg	11700	pg/mL
IL-6	2713	pg	10850	pg/mL
IL-8	2600	pg	10400	pg/mL
IL-10	4750	pg	19000	pg/mL
IL-12p70	7963	pg	31850	pg/mL
IL-13	3300	pg	13200	pg/mL
TNF- $\alpha$	5313	pg	21250	pg/mL
IL-5	5638	pg	22550	pg/mL
GM-CSF	16000	pg	64000	pg/mL
IL-17A	28500	pg	114000	pg/mL
VEGF-A	8600	pg	34400	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		06 JAN 2020