



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1

**Catalog Number:** C0060-2

**Lot Number:** A00U0204

**Expiration Date:** 28 FEB 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- $\gamma$	32875	pg	131500	pg/mL
IL-1 $\beta$	5163	pg	20650	pg/mL
IL-2	2525	pg	10100	pg/mL
IL-4	2400	pg	9600	pg/mL
IL-6	2538	pg	10150	pg/mL
IL-8	2863	pg	11450	pg/mL
IL-10	4950	pg	19800	pg/mL
IL-12p70	7663	pg	30650	pg/mL
IL-13	2650	pg	10600	pg/mL
TNF- $\alpha$	4363	pg	17450	pg/mL
IL-5	5463	pg	21850	pg/mL
GM-CSF	14625	pg	58500	pg/mL
IL-17A	30625	pg	122500	pg/mL
VEGF-A	4213	pg	16850	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		26 MAR 2021