



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 1  
**Catalog Number:** C0060-2  
**Lot Number:** A00U0343  
**Expiration Date:** 31 OCT 2028  
**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- $\gamma$	37250	pg	149000	pg/mL
IL-1 $\beta$	6050	pg	24200	pg/mL
IL-2	2575	pg	10300	pg/mL
IL-4	2275	pg	9100	pg/mL
IL-6	2337.5	pg	9350	pg/mL
IL-8	2700	pg	10800	pg/mL
IL-10	4962.5	pg	19850	pg/mL
IL-12p70	7037.5	pg	28150	pg/mL
IL-13	2300	pg	9200	pg/mL
TNF- $\alpha$	3287.5	pg	13150	pg/mL
IL-5	5275	pg	21100	pg/mL
GM-CSF	12625	pg	50500	pg/mL
IL-17A	24375	pg	97500	pg/mL
VEGF-A	3950	pg	15800	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		08 DEC 2023