



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 1  
**Catalog Number:** C0060-2  
**Lot Number:** A00U0409  
**Expiration Date:** 31 JAN 2030  
**Storage Temperature:** 2-8°C  
**Intended Assay Species:** Human, NHP

**Concentration Assignment:**

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	25125	pg	100500	pg/mL
IL-1 $\beta$	4812.5	pg	19250	pg/mL
IL-2	2212.5	pg	8850	pg/mL
IL-4	2662.5	pg	10650	pg/mL
IL-6	2350	pg	9400	pg/mL
IL-8	2487.5	pg	9950	pg/mL
IL-10	5475	pg	21900	pg/mL
IL-12p70	7512.5	pg	30050	pg/mL
IL-13	2575	pg	10300	pg/mL
TNF- $\alpha$	5862.5	pg	23450	pg/mL
IL-5	5125	pg	20500	pg/mL
GM-CSF	10750	pg	43000	pg/mL
IL-17A	26875	pg	107500	pg/mL
VEGF-A	6400	pg	25600	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		04 MAR 2025