



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 1  
**Catalog Number:** C0060-2  
**Lot Number:** A00U0401  
**Expiration Date:** 30 NOV 2029  
**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- $\gamma$	25500	pg	102000	pg/mL
IL-1 $\beta$	4875	pg	19500	pg/mL
IL-2	2437.5	pg	9750	pg/mL
IL-4	2637.5	pg	10550	pg/mL
IL-6	2425	pg	9700	pg/mL
IL-8	2612.5	pg	10450	pg/mL
IL-10	5175	pg	20700	pg/mL
IL-12p70	7637.5	pg	30550	pg/mL
IL-13	2675	pg	10700	pg/mL
TNF- $\alpha$	5875	pg	23500	pg/mL
IL-5	5150	pg	20600	pg/mL
GM-CSF	10700	pg	42800	pg/mL
IL-17A	27000	pg	108000	pg/mL
VEGF-A	6650	pg	26600	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		12 FEB 2025