



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 1  
**Catalog Number:** C0060-2  
**Lot Number:** A00U0419  
**Expiration Date:** 31 MAR 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- $\gamma$	20625	pg	82500	pg/mL
IL-1 $\beta$	5087.5	pg	20350	pg/mL
IL-2	2700	pg	10800	pg/mL
IL-4	2612.5	pg	10450	pg/mL
IL-6	2425	pg	9700	pg/mL
IL-8	2562.5	pg	10250	pg/mL
IL-10	5087.5	pg	20350	pg/mL
IL-12p70	7750	pg	31000	pg/mL
IL-13	2575	pg	10300	pg/mL
TNF- $\alpha$	4425	pg	17700	pg/mL
IL-5	5487.5	pg	21950	pg/mL
GM-CSF	11325	pg	45300	pg/mL
IL-17A	26375	pg	105500	pg/mL
VEGF-A	5087.5	pg	20350	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 MAY 2025