



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 3  
**Catalog Number:** C0062-2  
**Lot Number:** A00U0418  
**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
IL-1 $\alpha$	7575	pg	30300	pg/mL
IL-7	6462.5	pg	25850	pg/mL
IL-12/IL-23p40	33250	pg	133000	pg/mL
IL-15	3687.5	pg	14750	pg/mL
IL-16	25125	pg	100500	pg/mL
TNF- $\beta$	4462.5	pg	17850	pg/mL
G-CSF	13250	pg	53000	pg/mL
IFN- $\alpha$ 2a	48875	pg	195500	pg/mL
IL-18	16375	pg	65500	pg/mL
TPO	50750	pg	203000	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		16 MAY 2025