



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 3  
**Catalog Number:** C0062-2  
**Lot Number:** A00U0459  
**Expiration Date:** 31 OCT 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 $\alpha$	7225	pg	28900	pg/mL
IL-7	6000	pg	24000	pg/mL
IL-12/IL-23p40	31750	pg	127000	pg/mL
IL-15	3550	pg	14200	pg/mL
IL-16	32500	pg	130000	pg/mL
TNF- $\beta$	4425	pg	17700	pg/mL
G-CSF	12675	pg	50700	pg/mL
IFN- $\alpha$ 2a	47000	pg	188000	pg/mL
IL-18	15975	pg	63900	pg/mL
TPO	48750	pg	195000	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		11 DEC 2025