



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 3  
**Catalog Number:** C0062-2  
**Lot Number:** A00U0378  
**Expiration Date:** 31 MAY 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
IL-1 $\alpha$	7437.5	pg	29750	pg/mL
IL-7	6562.5	pg	26250	pg/mL
IL-12/IL-23p40	36125	pg	144500	pg/mL
IL-15	3850	pg	15400	pg/mL
IL-16	26750	pg	107000	pg/mL
TNF- $\beta$	4512.5	pg	18050	pg/mL
G-CSF	12500	pg	50000	pg/mL
IFN- $\alpha$ 2a	50625	pg	202500	pg/mL
IL-18	17125	pg	68500	pg/mL
TPO	53125	pg	212500	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		13 JUN 2024