



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 3
Catalog Number: C0062-2
Lot Number: A00U0443
Expiration Date: 30 APR 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 α	7912.5	pg	31650	pg/mL
IL-7	6700	pg	26800	pg/mL
IL-12/IL-23p40	34875	pg	139500	pg/mL
IL-15	3887.5	pg	15550	pg/mL
IL-16	34625	pg	138500	pg/mL
TNF- β	4700	pg	18800	pg/mL
G-CSF	12387.5	pg	49550	pg/mL
IFN- α 2a	50250	pg	201000	pg/mL
IL-18	16750	pg	67000	pg/mL
TPO	50750	pg	203000	pg/mL

Reconstitution Volume: 250 μ 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
Review/Approval	Karen Hamilla	Quality		29 MAY 2025