



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 3
Catalog Number: C0062-2
Lot Number: A00U0478
Expiration Date: 31 MAY 2031
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	7525	pg	30100	pg/mL
IL-7	5825	pg	23300	pg/mL
IL-12/IL-23p40	34500	pg	138000	pg/mL
IL-15	3700	pg	14800	pg/mL
IL-16	34500	pg	138000	pg/mL
TNF- β	4575	pg	18300	pg/mL
G-CSF	12650	pg	50600	pg/mL
IFN- α 2a	48000	pg	192000	pg/mL
IL-18	16425	pg	65700	pg/mL
TPO	47750	pg	191000	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		17 JUL 2026