



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0215

Expiration Date: 31 AUG 2026

Storage Temperature: 2-8°C

Intended Assay Species: Human

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-1 α	8275	pg	33100	pg/mL
IL-7	8213	pg	32850	pg/mL
IL-12/IL-23p40	30250	pg	121000	pg/mL
IL-15	4000	pg	16000	pg/mL
IL-16	31375	pg	125500	pg/mL
TNF- β	4925	pg	19700	pg/mL
G-CSF	12750	pg	51000	pg/mL
IFN- α 2a	48000	pg	192000	pg/mL
IL-18	17625	pg	70500	pg/mL
TPO	35875	pg	143500	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		01 SEP 2021