



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0168

Expiration Date: 31 AUG 2025

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	7100	pg	28400	pg/mL
IL-7	7988	pg	31950	pg/mL
IL-12/IL-23p40	27000	pg	108000	pg/mL
IL-15	3813	pg	15250	pg/mL
IL-16	27875	pg	111500	pg/mL
TNF- β	5213	pg	20850	pg/mL
IFN- α 2a	44375	pg	177500	pg/mL
G-CSF	23000	pg	92000	pg/mL
IL-18	46750	pg	187000	pg/mL
TPO	41875	pg	167500	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		04 SEP 2020