



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0245

Expiration Date: 31 JAN 2027

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	7400	pg	29600	pg/mL
IL-7	8888	pg	35550	pg/mL
IL-12/IL-23p40	32375	pg	129500	pg/mL
IL-15	3838	pg	15350	pg/mL
IL-16	24750	pg	99000	pg/mL
TNF- β	4488	pg	17950	pg/mL
G-CSF	12750	pg	51000	pg/mL
IFN- α 2a	47875	pg	191500	pg/mL
IL-18	16500	pg	66000	pg/mL
TPO	50250	pg	201000	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		23 MAR 2022