



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0257

Expiration Date: 30 JUN 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	7063	pg	28250	pg/mL
IL-7	6163	pg	24650	pg/mL
IL-12/IL-23p40	30625	pg	122500	pg/mL
IL-15	3713	pg	14850	pg/mL
IL-16	26375	pg	105500	pg/mL
TNF- β	4013	pg	16050	pg/mL
G-CSF	11550	pg	46200	pg/mL
IFN- α 2a	45250	pg	181000	pg/mL
IL-18	15875	pg	63500	pg/mL
TPO	45250	pg	181000	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		11 AUG 2022