



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0239

Expiration Date: 31 OCT 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	7850	pg	31400	pg/mL
IL-7	8088	pg	32350	pg/mL
IL-12/IL-23p40	30000	pg	120000	pg/mL
IL-15	3750	pg	15000	pg/mL
IL-16	25250	pg	101000	pg/mL
TNF- β	4725	pg	18900	pg/mL
G-CSF	12875	pg	51500	pg/mL
IFN- α 2a	48750	pg	195000	pg/mL
IL-18	18750	pg	75000	pg/mL
TPO	44375	pg	177500	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	Carissa B. Hubbard	Quality		24 NOV 2021