



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0228

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	7813	pg	31250	pg/mL
IL-7	8188	pg	32750	pg/mL
IL-12/IL-23p40	29125	pg	116500	pg/mL
IL-15	3900	pg	15600	pg/mL
IL-16	26500	pg	106000	pg/mL
TNF- β	5225	pg	20900	pg/mL
G-CSF	12450	pg	49800	pg/mL
IFN- α 2a	46375	pg	185500	pg/mL
IL-18	49000	pg	196000	pg/mL
TPO	39500	pg	158000	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	Xiomara Talavera	Quality		07 OCT 2021