



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3
Catalog Number: C0062-2
Lot Number: A00U0164
Expiration Date: 31 JUL 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	7125	pg	28500	pg/mL
IL-7	7788	pg	31150	pg/mL
IL-12/IL-23p40	27250	pg	109000	pg/mL
IL-15	4050	pg	16200	pg/mL
IL-16	28500	pg	114000	pg/mL
TNF- β	5338	pg	21350	pg/mL
G-CSF	23750	pg	95000	pg/mL
IFN- α 2a	46250	pg	185000	pg/mL
IL-18	47125	pg	188500	pg/mL
TPO	41250	pg	165000	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		31 JUL 2020