



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0106

Expiration Date: 30 APR 2024

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	6 638	pg	26 550	pg/mL
IL-7	7 175	pg	28 700	pg/mL
IL-12/IL-23p40	25 750	pg	103 000	pg/mL
IL-15	3 688	pg	14 750	pg/mL
IL-16	32 500	pg	130 000	pg/mL
TNF- β	4 888	pg	19 550	pg/mL
G-CSF	23 500	pg	94 000	pg/mL
IFN- α 2 α	46 375	pg	185 500	pg/mL
IL-18	53 500	pg	214 000	pg/mL
TPO	43 875	pg	175 500	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	Dawn Stollar	Quality		24 APR 2019