



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0134

Expiration Date: 30 NOV 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	4913	pg	19650	pg/mL
IL-7	7225	pg	28900	pg/mL
IL-12/IL-23p40	26625	pg	106500	pg/mL
IL-15	4375	pg	17500	pg/mL
IL-16	33500	pg	134000	pg/mL
TNF- β	3988	pg	15950	pg/mL
G-CSF	25500	pg	102000	pg/mL
IFN- α 2a	43375	pg	173500	pg/mL
IL-18	53250	pg	213000	pg/mL
TPO	41750	pg	167000	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		06 DEC 2019