



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 3
Catalog Number: C0062-2
Lot Number: A00U0402
Expiration Date: 30 NOV 2029
Storage Temperature: 2-8°C
Intended Assay Species: Human, NHP

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-1 α	7162.5	pg	28650	pg/mL
IL-7	6600	pg	26400	pg/mL
IL-12/IL-23p40	33750	pg	135000	pg/mL
IL-15	3787.5	pg	15150	pg/mL
IL-16	25625	pg	102500	pg/mL
TNF- β	4450	pg	17800	pg/mL
G-CSF	12625	pg	50500	pg/mL
IFN- α 2a	48750	pg	195000	pg/mL
IL-18	20125	pg	80500	pg/mL
TPO	49625	pg	198500	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		20 DEC 2024