



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 3
Catalog Number: C0062-2
Lot Number: A00U0379
Expiration Date: 30 APR 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	7375	pg	29500	pg/mL
IL-7	6287.5	pg	25150	pg/mL
IL-12/IL-23p40	34250	pg	137000	pg/mL
IL-15	3562.5	pg	14250	pg/mL
IL-16	27375	pg	109500	pg/mL
TNF- β	3887.5	pg	15550	pg/mL
G-CSF	12262.5	pg	49050	pg/mL
IFN- α 2a	47500	pg	190000	pg/mL
IL-18	23500	pg	94000	pg/mL
TPO	51000	pg	204000	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		18 JUN 2024