



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 6
Catalog Number: C0072-2
Lot Number: A00U0432
Expiration Date: 31 JUL 2030
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-17A/F	17250	pg	69000	pg/mL
IL-17E/IL-25	9175	pg	36700	pg/mL
IL-17F	140000	pg	560000	pg/mL
IL-21	16125	pg	64500	pg/mL
IL-22	3612.5	pg	14450	pg/mL
IL-23	34750	pg	139000	pg/mL
IL-27	51750	pg	207000	pg/mL
IL-29/IFN- λ 1	10575	pg	42300	pg/mL
IL-31	11750	pg	47000	pg/mL
IL-33	10925	pg	43700	pg/mL
TSLP	9337.5	pg	37350	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		27 AUG 2025