



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 6
Catalog Number: C0072-2
Lot Number: A00U0463
Expiration Date: 31 JAN 2031
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	17325	pg	69300	pg/mL
IL-17E/IL-25	8350	pg	33400	pg/mL
IL-17F	152750	pg	611000	pg/mL
IL-21	17125	pg	68500	pg/mL
IL-22	3600	pg	14400	pg/mL
IL-23	35000	pg	140000	pg/mL
IL-27	45500	pg	182000	pg/mL
IL-29/IFN- λ 1	10500	pg	42000	pg/mL
IL-31	11850	pg	47400	pg/mL
IL-33	11175	pg	44700	pg/mL
TSLP	9100	pg	36400	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 MAR 2026