



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 6

Catalog Number: C0072-2

Lot Number: A00U0237

Expiration Date: 31 OCT 2026

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-17A/F	18875	pg	75500	pg/mL
IL-17E/IL-25	9200	pg	36800	pg/mL
IL-17F	124375	pg	497500	pg/mL
IL-21	19875	pg	79500	pg/mL
IL-22	3325	pg	13300	pg/mL
IL-23	33125	pg	132500	pg/mL
IL-27	36750	pg	147000	pg/mL
IL-29/IFN-λ1	9650	pg	38600	pg/mL
IL-31	11388	pg	45550	pg/mL
IL-33	12325	pg	49300	pg/mL
TSLP	8613	pg	34450	pg/mL

Reconstitution Volume: 250 µL

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		21 APR 2022