



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 6

Catalog Number: C0072 -2

Lot Number: A00U0094

Expiration Date: 31 MAY 2024

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	18 750	pg	75 000	pg/mL
IL-17E/IL-25	9 375	pg	37 500	pg/mL
IL-17F	128 750	pg	515 000	pg/mL
IL-21	11 900	pg	47 600	pg/mL
IL-22	3 888	pg	15 550	pg/mL
IL-23	27 375	pg	109 500	pg/mL
IL-27	49 750	pg	199 000	pg/mL
IL-29/IFN- λ 1	10 688	pg	42 750	pg/mL
IL-31	12 188	pg	48 750	pg/mL
IL-33	11 738	pg	46 950	pg/mL
TSLP	9 713	pg	38 850	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	Dawn Stollar	Quality		28 MAY 2019