



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 6

Catalog Number: C0072-2

Lot Number: A00U0190

Expiration Date: 31 MAY 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	18125	pg	72500	pg/mL
IL-17E/IL-25	9075	pg	36300	pg/mL
IL-17F	107875	pg	431500	pg/mL
IL-21	19250	pg	77000	pg/mL
IL-22	3500	pg	14000	pg/mL
IL-23	31375	pg	125500	pg/mL
IL-27	38375	pg	153500	pg/mL
IL-29/IFN- λ 1	9400	pg	37600	pg/mL
IL-31	10638	pg	42550	pg/mL
IL-33	12188	pg	48750	pg/mL
TSLP	8475	pg	33900	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		01 JUL 2021