



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 6

Catalog Number: C0072-2

Lot Number: A00U0143

Expiration Date: 31 JAN 2025

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	16875	pg	67500	pg/mL
IL-17E/IL-25	8900	pg	35600	pg/mL
IL-17F	131250	pg	525000	pg/mL
IL-21	32750	pg	131000	pg/mL
IL-22	3613	pg	14450	pg/mL
IL-23	32125	pg	128500	pg/mL
IL-27	36000	pg	144000	pg/mL
IL-29/IFN- λ 1	11038	pg	44150	pg/mL
IL-31	11463	pg	45850	pg/mL
IL-33	11663	pg	46650	pg/mL
TSLP	10075	pg	40300	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		15 JUN 2020