



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 6
Catalog Number: C0072-2
Lot Number: A00U0295
Expiration Date: 31 JAN 2028
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	21375	pg	85500	pg/mL
IL-17E/IL-25	8525	pg	34100	pg/mL
IL-17F	127500	pg	510000	pg/mL
IL-21	18375	pg	73500	pg/mL
IL-22	3625	pg	14500	pg/mL
IL-23	34750	pg	139000	pg/mL
IL-27	37750	pg	151000	pg/mL
IL-29/IFN- λ 1	10675	pg	42700	pg/mL
IL-31	11375	pg	45500	pg/mL
IL-33	13000	pg	52000	pg/mL
TSLP	9125	pg	36500	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		22 FEB 2023