



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 6  
**Catalog Number:** C0072-2  
**Lot Number:** A00U0332  
**Expiration Date:** 30 JUN 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	20375	pg	81500	pg/mL
IL-17E/IL-25	9400	pg	37600	pg/mL
IL-17F	141250	pg	565000	pg/mL
IL-21	17000	pg	68000	pg/mL
IL-22	3387.5	pg	13550	pg/mL
IL-23	32875	pg	131500	pg/mL
IL-27	39750	pg	159000	pg/mL
IL-29/IFN-λ1	11200	pg	44800	pg/mL
IL-31	11437.5	pg	45750	pg/mL
IL-33	11675	pg	46700	pg/mL
TSLP	9687.5	pg	38750	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 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
<b>Review/Approval</b>	Karen Hamilla	Quality		08 AUG 2023