



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 6  
**Catalog Number:** C0072-2  
**Lot Number:** A00U0385  
**Expiration Date:** 30 SEP 2029  
**Storage Temperature:** 2-8°C  
**Intended Assay Species:** Human, NHP

**Concentration Assignment:**

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A/F	21000	pg	84000	pg/mL
IL-17E/IL-25	7875	pg	31500	pg/mL
IL-17F	151250	pg	605000	pg/mL
IL-21	20750	pg	83000	pg/mL
IL-22	3462.5	pg	13850	pg/mL
IL-23	36250	pg	145000	pg/mL
IL-27	47375	pg	189500	pg/mL
IL-29/IFN-λ1	11462.5	pg	45850	pg/mL
IL-31	11050	pg	44200	pg/mL
IL-33	11350	pg	45400	pg/mL
TSLP	8762.5	pg	35050	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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		01 NOV 2024