



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 6  
**Catalog Number:** C0072-2  
**Lot Number:** A00U0339  
**Expiration Date:** 30 NOV 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	21125	pg	84500	pg/mL
IL-17E/IL-25	8837.5	pg	35350	pg/mL
IL-17F	132500	pg	530000	pg/mL
IL-21	19250	pg	77000	pg/mL
IL-22	3400	pg	13600	pg/mL
IL-23	35625	pg	142500	pg/mL
IL-27	40250	pg	161000	pg/mL
IL-29/IFN-λ1	11062.5	pg	44250	pg/mL
IL-31	11425	pg	45700	pg/mL
IL-33	12412.5	pg	49650	pg/mL
TSLP	9612.5	pg	38450	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		17 JAN 2024