



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 6

**Catalog Number:** C0072-2

**Lot Number:** A00U0283

**Expiration Date:** 30 SEP 2027

**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	18000	pg	72000	pg/mL
IL-17E/IL-25	9450	pg	37800	pg/mL
IL-17F	150000	pg	600000	pg/mL
IL-21	20375	pg	81500	pg/mL
IL-22	3350	pg	13400	pg/mL
IL-23	32875	pg	131500	pg/mL
IL-27	35250	pg	141000	pg/mL
IL-29/IFN- $\lambda$ 1	10500	pg	42000	pg/mL
IL-31	11225	pg	44900	pg/mL
IL-33	12238	pg	48950	pg/mL
TSLP	9088	pg	36350	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		19 OCT 2022