



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 7

**Catalog Number:** C0073-2

**Lot Number:** A00U0249

**Expiration Date:** 28 FEB 2027

**Storage Temperature:** 2-8°C

**Intended Assay Species:** Mouse

### Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-16	7275	pg	29100	pg/mL
IL-17A	2775	pg	11100	pg/mL
IL-17C	48875	pg	195500	pg/mL
IL-17E/IL-25	16125	pg	64500	pg/mL
IL-21	42875	pg	171500	pg/mL
IL-22	2188	pg	8750	pg/mL
IL-23	24125	pg	96500	pg/mL
IL-12/IL-23p40	24125	pg	96500	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		26 APR 2022