



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 7  
**Catalog Number:** C0073-2  
**Lot Number:** A00U0108  
**Expiration Date:** 31 MAY 2024  
**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	7 713	pg	30 850	pg/mL
IL-17A	3 438	pg	13 750	pg/mL
IL-17C	55 875	pg	223 500	pg/mL
IL-17E/IL-25	20 875	pg	83 500	pg/mL
IL-21	37 500	pg	150 000	pg/mL
IL-22	2 263	pg	9 050	pg/mL
IL-23	23 875	pg	95 500	pg/mL
IL-12/IL-23p40	23 875	pg	95 500	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		26 AUG 2019