



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 7  
**Catalog Number:** C0073-2  
**Lot Number:** A00U0310  
**Expiration Date:** 31 MAR 2028  
**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	7487.5	pg	29950	pg/mL
IL-17A	2687.5	pg	10750	pg/mL
IL-17C	46625	pg	186500	pg/mL
IL-17E/IL-25	17750	pg	71000	pg/mL
IL-21	43875	pg	175500	pg/mL
IL-22	2125	pg	8500	pg/mL
IL-23	24750	pg	99000	pg/mL
IL-12/IL-23p40	24750	pg	99000	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. 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		27 APR 2023