



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 8  
**Catalog Number:** C0074-2  
**Lot Number:** A00U0315  
**Expiration Date:** 29 FEB 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-15	122500	pg	490000	pg/mL
IL-17F	67875	pg	271500	pg/mL
IL-31	86750	pg	347000	pg/mL
IL-33	44750	pg	179000	pg/mL
IL-27p28/IL-30	94625	pg	378500	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		16 JUN 2023