



CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 8  
**Catalog Number:** C0074-2  
**Lot Number:** A00U0340  
**Expiration Date:** 28 FEB 2029  
**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	132500	pg	530000	pg/mL
IL-17F	62375	pg	249500	pg/mL
IL-31	86375	pg	345500	pg/mL
IL-33	42875	pg	171500	pg/mL
IL-27p28/IL-30	90500	pg	362000	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		11 APR 2024