



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 8

**Catalog Number:** C0074-2

**Lot Number:** A00U0238

**Expiration Date:** 30 NOV 2026

**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	119500	pg	478000	pg/mL
IL-17F	61125	pg	244500	pg/mL
IL-31	83875	pg	335500	pg/mL
IL-33	46875	pg	187500	pg/mL
IL-27p28/IL-30	104625	pg	418500	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
Review/Approval	Karen Hamilla	Quality		07 APR 2022