



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 8

**Catalog Number:** C0074-2

**Lot Number:** A00U0133

**Expiration Date:** 30 NOV 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-15	132500	pg	530000	pg/mL
IL-17F	55125	pg	220500	pg/mL
IL-31	75625	pg	302500	pg/mL
IL-33	40375	pg	161500	pg/mL
IL-27p28/IL-30	98250	pg	393000	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		10 JAN 2020