



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 8
Catalog Number: C0074-2
Lot Number: A00U0372
Expiration Date: 30 JUN 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	137500	pg	550000	pg/mL
IL-17F	74250	pg	297000	pg/mL
IL-31	83250	pg	333000	pg/mL
IL-33	46500	pg	186000	pg/mL
IL-27p28/IL-30	86500	pg	346000	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		17 JUL 2024