



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 7
Catalog Number: C0073-2
Lot Number: A00U0392
Expiration Date: 31 JAN 2030
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-16	6412.5	pg	25650	pg/mL
IL-17A	2612.5	pg	10450	pg/mL
IL-17C	44875	pg	179500	pg/mL
IL-17E/IL-25	16250	pg	65000	pg/mL
IL-21	61125	pg	244500	pg/mL
IL-22	1937.5	pg	7750	pg/mL
IL-23	21875	pg	87500	pg/mL
IL-12/IL-23p40	21875	pg	87500	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		24 FEB 2025