



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 7
Catalog Number: C0073-2
Lot Number: A00U0361
Expiration Date: 31 MAR 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	Unit of Measure
IL-16	7012.5	pg	28050	pg/mL
IL-17A	2787.5	pg	11150	pg/mL
IL-17C	42625	pg	170500	pg/mL
IL-17E/IL-25	17000	pg	68000	pg/mL
IL-21	54500	pg	218000	pg/mL
IL-22	2125	pg	8500	pg/mL
IL-23	23000	pg	92000	pg/mL
IL-12/IL-23p40	23000	pg	92000	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		01 MAY 2024