



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 7

Catalog Number: C0073-2

Lot Number: A00U0219

Expiration Date: 30 JUN 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-16	7100	pg	28400	pg/mL
IL-17A	2563	pg	10250	pg/mL
IL-17C	47625	pg	190500	pg/mL
IL-17E/IL-25	15250	pg	61000	pg/mL
IL-21	30000	pg	120000	pg/mL
IL-22	2200	pg	8800	pg/mL
IL-23	23625	pg	94500	pg/mL
IL-12/IL-23p40	23625	pg	94500	pg/mL

Reconstitution Volume: 250 μ L

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		22 OCT 2021