



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 7

Catalog Number: C0073-2

Lot Number: A00U0191

Expiration Date: 31 JAN 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	7213	pg	28850	pg/mL
IL-17A	2638	pg	10550	pg/mL
IL-17C	55250	pg	221000	pg/mL
IL-17E/IL-25	16375	pg	65500	pg/mL
IL-21	36250	pg	145000	pg/mL
IL-22	2213	pg	8850	pg/mL
IL-23	26000	pg	104000	pg/mL
IL-12/IL-23p40	26000	pg	104000	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		15 MAR 2021