



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 7
Catalog Number: C0073-2
Lot Number: A00U0336
Expiration Date: 31 JUL 2028
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	6725	pg	26900	pg/mL
IL-17A	2562.5	pg	10250	pg/mL
IL-17C	43375	pg	173500	pg/mL
IL-17E/IL-25	16375	pg	65500	pg/mL
IL-21	37500	pg	150000	pg/mL
IL-22	2125	pg	8500	pg/mL
IL-23	22625	pg	90500	pg/mL
IL-12/IL-23p40	22625	pg	90500	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		06 SEP 2023