



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 8
Catalog Number: C0074-2
Lot Number: A00U0296
Expiration Date: 30 NOV 2027
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-15	138750	pg	555000	pg/mL
IL-17F	65375	pg	261500	pg/mL
IL-31	86625	pg	346500	pg/mL
IL-33	45375	pg	181500	pg/mL
IL-27p28/IL-30	100250	pg	401000	pg/mL

Reconstitution Volume: 250 µL 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		04 JAN 2023