



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 8

Catalog Number: C0074-2

Lot Number: A00U0059

Expiration Date: 31 AUG 2023

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
ms IL-15	137 500	pg	550 000	pg/mL
ms IL-17F	59 500	pg	238 000	pg/mL
ms IL-31	80 750	pg	323 000	pg/mL
ms IL-33	40 500	pg	162 000	pg/mL
ms IL-27p28/IL-30	102 125	pg	408 500	pg/mL

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.

Statement: The above calibrator lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	30 AUG 2018