



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 7

Catalog Number: C0073-2

Lot Number: A00U0065

Expiration Date: 30 JUN 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-16	7 300	pg	29 200	pg/mL
ms IL-17A	2 588	pg	10 350	pg/mL
ms IL-17C	49 500	pg	198 000	pg/mL
ms IL-17E	19 875	pg	79 500	pg/mL
ms IL-21	45 375	pg	181 500	pg/mL
ms IL-22	2 213	pg	8 850	pg/mL
ms IL-23	24 000	pg	96 000	pg/mL
ms IL-12/IL-23p40	24 000	pg	96 000	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>	29 AUG 2018