



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 7

Catalog Number: C0073-2

Lot Number: A00U0019

Expiration Date: 31 MAR 2021

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator (pg)	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL) (pg/mL)	Concentration of Standard 1 (after 5-fold Dilution) (pg/mL)
ms IL-12/IL-23p40	25 625	102 500	20 500
ms IL-16	7 675	30 700	6 140
ms IL-17A	2 675	10 700	2 140
ms IL-17C	78 750	315 000	63 000
ms IL-17E	23 375	93 500	18 700
ms IL-21	66 750	267 000	53 400
ms IL-22	2 338	9 350	1 870
ms IL-23	25 625	102 500	20 500

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>	21 JUL 2017