



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 7

Catalog Number: C0073-2

Lot Number: A00U0020

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	25 375	101 500	20 300
ms IL-16	7 675	30 700	6 140
ms IL-17A	2 650	10 600	2 120
ms IL-17C	57 250	229 000	45 800
ms IL-17E	23 625	94 500	18 900
ms IL-21	50 250	201 000	40 200
ms IL-22	2 288	9 150	1 830
ms IL-23	25 375	101 500	20 300

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