



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 7

Catalog Number: C0073-2

Lot Number: A00U0023

Expiration Date: 30 APR 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)
Mouse IL-12/IL-23p40	23 875	95 500	19 100
Mouse IL-16	7 888	31 550	6 310
Mouse IL-17A	2 650	10 600	2 120
Mouse IL-17C	62 625	250 500	50 100
Mouse IL-17E	22 750	91 000	18 200
Mouse IL-21	41 625	166 500	33 300
Mouse IL-22	2 275	9 100	1 820
Mouse IL-23	23 875	95 500	19 100

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>	14 NOV 2017