



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 6

Catalog Number: C0072-2

Lot Number: A00U0018

Expiration Date: 28 FEB 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)
Human IL-17AF	23 000	92 000	18 400
Human IL-17E/IL-25	11 563	46 250	9 250
Human IL-17F	140 000	560 000	112 000
Human IL-21	15 750	63 000	12 600
Human IL-22	4 275	17 100	3 420
Human IL-23	27 125	108 500	21 700
Human IL-27	63 375	253 500	50 700
Human IL-29	14 750	59 000	11 800
Human IL-31	14 000	56 000	11 200
Human IL-33	13 000	52 000	10 400
Human TSLP	12 625	50 500	10 100

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>	22 FEB 2017