



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0006

Expiration Date: 31 MAR 2020

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 IFN- γ	24 375	97 500	19 500
Human IL-1 β	4 963	19 850	3 970
Human IL-2	2 538	10 150	2 030
Human IL-4	2 925	11 700	2 340
Human IL-6	2 613	10 450	2 090
Human IL-8	2 900	11 600	2 320
Human IL-10	4 913	19 650	3 930
Human IL-12p70	7 138	28 550	5 710
Human IL-13	2 575	10 300	2 060
Human TNF- α	4 825	19 300	3 860
Human IL-5	5 525	22 100	4 420
Human GM-CSF	12 750	51 000	10 200
Human IL-17A	31 250	125 000	25 000
Human VEGF-A	6 163	24 650	4 930

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