



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0045

Expiration Date: 31 JAN 2023

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)
hu IFN- γ	33 250	133 000	26 600
hu IL-1 β	4 463	17 850	3 570
hu IL-2	2 475	9 900	1 980
hu IL-4	3 250	13 000	2 600
hu IL-6	2 550	10 200	2 040
hu IL-8	2 550	10 200	2 040
hu IL-10	4 163	16 650	3 330
hu IL-12p70	7 925	31 700	6 340
hu IL-13	3 075	12 300	2 460
hu TNF- α	4 438	17 750	3 550
hu IL-5	4 825	19 300	3 860
hu GM-CSF	12 025	48 100	9 620
hu IL-17A	22 000	88 000	17 600
hu VEGF-A	7 413	29 650	5 930

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>	19 JAN 2018