



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0025

Expiration Date: 31 JUL 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-1 α	6 675	26 700	5 340
Human IL-7	9 225	36 900	7 380
Human IL-12/23p40	28 250	113 000	22 600
Human IL-15	3 913	15 650	3 130
Human IL-16	28 000	112 000	22 400
Human TNF- β	5 950	23 800	4 760
Human G-CSF	27 625	110 500	22 100
Human IFN- α 2a	51 000	204 000	40 800
Human IL-18	56 125	224 500	44 900
Human TPO	51 500	206 000	41 200

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>	17 JUL 2017