



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0040

Expiration Date: 31 JUL 2022

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	6 475	25 900	5 180
Human IL-7	7 663	30 650	6 130
Human IL-12/IL-23p40	25 375	101 500	20 300
Human IL-15	3 675	14 700	2 940
Human IL-16	25 875	103 500	20 700
Human TNF- β	4 938	19 750	3 950
Human G-CSF	26 000	104 000	20 800
Human IFN- α 2a	50 375	201 500	40 300
Human IL-18	51 000	204 000	40 800
Human TPO	48 750	195 000	39 000

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>	01 AUG 2017