



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0038

Expiration Date: 31 MAY 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)
hu IL1α	7038	28 150	5630
hu IL-7	8263	33 050	6610
hu IL-12/IL-23p40	27375	109 500	21900
hu IL-15	3838	15 350	3070
hu IL-16	28000	112 000	22400
hu TNF-β	6313	25 250	5050
hu G-CSF	29125	116 500	23300
hu IFN-α2a	55750	223 000	44600
hu IL-18	97125	388 500	77700
hu TPO	51875	207 500	41500

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>	26 JUN 2017