



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 3

**Catalog Number:** C0062-2

**Lot Number:** A00U0036

**Expiration Date:** 31 MAR 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 $\alpha$	6 238	24 950	4 990
Human IL-7	8 538	34 150	6 830
Human IL-12/IL-23p40	24 875	99 500	19 900
Human IL-15	3 600	14 400	2 880
Human IL-16	26 375	105 500	21 100
Human TNF- $\beta$	5 513	22 050	4 410
Human G-CSF	26 125	104 500	20 900
Human IFN- $\alpha$ 2a	49 500	198 000	39 600
Human IL-18	47 625	190 500	38 100
Human TPO	49 750	199 000	39 800

**Statement:** The above calibrator lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	18 JUL 2017