



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1

**Catalog Number:** C0060-2

**Lot Number:** A00U0007

**Expiration Date:** 30 APR 2020

**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 IFN- $\gamma$	21 875	87 500	17 500
Human IL-1 $\beta$	4 850	19 400	3 880
Human IL-2	2 458	9 832	1 966
Human IL-4	2 450	9 800	1 960
Human IL-6	2 550	10 200	2 040
Human IL-8	2 875	11 500	2 300
Human IL-10	4775	19 100	3 820
Human IL-12p70	6 975	27 900	5 580
Human IL-13	2 470	9 880	1 976
Human TNF- $\alpha$	4 725	18 900	3 780
Human IL-5	5 125	20 500	4 100
Human GM-CSF	12 300	49 200	9 840
Human IL-17A	29 250	117 000	23 400
Human VEGF-A	7 100	28 400	5 680

**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>	06 APR 2017