



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1

**Catalog Number:** C0060-2

**Lot Number:** A00U0032

**Expiration Date:** 28 FEB 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 IFN-gamma	23 300	93 200	18 640
hu IL-1β	4 700	18 800	3 760
hu IL-2	2 300	9 200	1 840
hu IL-4	2 048	8 190	1 638
hu IL-6	2 525	10 100	2 020
hu IL-8	2 310	9 240	1 848
hu IL-10	4 675	18 700	3 740
hu IL-12p70	6 750	27 000	5 400
hu IL-13	2 300	9 200	1 840
hu TNF-α	4 375	17 500	3 500
hu IL-5	5 250	21 000	4 200
hu GM-CSF	12 625	50 500	10 100
hu IL-17A	22 200	88 800	17 760
hu VEGF-A	6 050	24 200	4 840

**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>	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	23MAR2017