



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 9

**Catalog Number:** C0090-2

**Lot Number:** A00U0046

**Expiration Date:** 30 NOV 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 EPO	25 250	101 000	20 200
Human FLT3L	8 025	32 100	6 420
Human IFN-β	123 750	495 000	99 000
Human IL-3	20 500	82 000	16 400
Human IL-9	1 450	5 800	1 160
Human IL-17B	4 188	16 750	3 350
Human IL-17C	26 625	106 500	21 300
Human IL-17D	35 750	143 000	28 600
Human IL-1RA	4 200	16 800	3 360
Human IL-2Rα	75 000	300 000	60 000

**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>	21 DEC 2017