



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0029

Expiration Date: 30 NOV 2021

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	26 375	105 500	21 100
Human FLT3L	8 250	33 000	6 600
Human IFN-β	125 000	500 000	100 000
Human IL-3	22 000	88 000	17 600
Human IL-9	1 538	6 150	1 230
Human IL-17B	4 275	17 100	3 420
Human IL-17C	27 500	110 000	22 000
Human IL-17D	35 125	140 500	28 100
Human IL-1RA	5 063	20 250	4 050
Human IL-2Rα	73 375	293 500	58 700

Statement: The above calibrator lot was tested using MSD procedures and conforms to MSD's specifications. It 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>	15 FEB 2017