



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 12

Catalog Number: C0092-2

Lot Number: A00U0033

Expiration Date: 31 DEC 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)
ms IL-9	11 388	45 550	9 110
ms IL-17 A/F	13 000	52 000	10 400
ms IP-10	6 575	26 300	5 260
ms MCP-1	1 625	6 500	1 300
ms MIP-1 α	2 488	9 950	1 990
ms MIP-1 β	33 750	135 000	27 000
ms MIP-2	2 550	10 200	2 040
ms MIP-3 α	3 525	14 100	2 820

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