



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 12

Catalog Number: C0092-2

Lot Number: A00U0048

Expiration Date: 31 DEC 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)
ms IL-9	12 013	48 050	9 610
ms IL-17 A/F	13 750	55 000	11 000
ms IP-10	7 275	29 100	5 820
ms MCP-1	1 688	6 750	1 350
ms MIP-1 α	2 488	9 950	1 990
ms MIP-1 β	34 500	138 000	27 600
ms MIP-2	2 600	10 400	2 080
ms MIP-3 α	3 713	14 850	2 970

Statement: The above calibrator lot 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>	23 JAN 2018