



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 12

Catalog Number: C0092-2

Lot Number: A00U0076

Expiration Date: 31 AUG 2023

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
ms IL-9	11 563	pg	46 250	pg/mL
ms IL-17A/F	12 875	pg	51 500	pg/mL
ms IP-10	7 213	pg	28 850	pg/mL
ms MCP-1	1 588	pg	6 350	pg/mL
ms MIP-1 α	2 438	pg	9 750	pg/mL
ms MIP-1 β	34 250	pg	137 000	pg/mL
ms MIP-2	2 550	pg	10 200	pg/mL
ms MIP-3 α	3 375	pg	13 500	pg/mL

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

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>	27 SEP 2018