



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 12

Catalog Number: C0092-2

Lot Number: A00U0056

Expiration Date: 31 MAY 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	11 413	pg	45 650	pg/mL
ms IL-17A/F	12 475	pg	49 900	pg/mL
ms IP-10	7 575	pg	30 300	pg/mL
ms MCP-1	1 575	pg	6 300	pg/mL
ms MIP-1 α	2 413	pg	9 650	pg/mL
ms MIP-1 β	33 500	pg	134 000	pg/mL
ms MIP-2	2 563	pg	10 250	pg/mL
ms MIP-3 α	3 438	pg	13 750	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>	29 AUG 2018