



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0051

Expiration Date: 28 FEB 2023

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
ms IFN- γ	3 600	pg	14 400	pg/mL
ms IL-1 β	15 750	pg	63 000	pg/mL
ms IL-2	13 500	pg	54 000	pg/mL
ms IL-4	12 200	pg	48 800	pg/mL
ms IL-5	3 125	pg	12 500	pg/mL
ms IL-6	19 375	pg	77 500	pg/mL
ms KC	2 650	pg	10 600	pg/mL
ms IL-10	27 500	pg	110 000	pg/mL
ms IL-12p70	102 875	pg	411 500	pg/mL
ms TNF- α	7 513	pg	30 050	pg/mL
ms GM-CSF	1 275	pg	5 100	pg/mL
ms IL-13	23 250	pg	93 000	pg/mL
ms EPO	14 625	pg	58 500	pg/mL
ms VEGF-A	15 500	pg	62 000	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