



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0074

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 IFN- γ	3 063	pg	12 250	pg/mL
ms IL-1 β	20 000	pg	80 000	pg/mL
ms IL-2	14 250	pg	57 000	pg/mL
ms IL-4	12 625	pg	50 500	pg/mL
ms IL-5	3 188	pg	12 750	pg/mL
ms IL-6	19 500	pg	78 000	pg/mL
ms KC	2 625	pg	10 500	pg/mL
ms IL-10	28 125	pg	112 500	pg/mL
ms IL-12p70	107 250	pg	429 000	pg/mL
ms TNF- α	8 750	pg	35 000	pg/mL
ms GM-CSF	1 338	pg	5 350	pg/mL
ms IL-13	23 750	pg	95 000	pg/mL
ms EPO	16 875	pg	67 500	pg/mL
ms VEGF-A	15 000	pg	60 000	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>	28 AUG 2018