



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0084

Expiration Date: 31 OCT 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	15 125	pg	60 500	pg/mL
ms IL-2	12 625	pg	50 500	pg/mL
ms IL-4	12 875	pg	51 500	pg/mL
ms IL-5	3 450	pg	13 800	pg/mL
ms IL-6	18 875	pg	75 500	pg/mL
ms KC	2 688	pg	10 750	pg/mL
ms IL-10	27 875	pg	111 500	pg/mL
ms IL-12p70	102 375	pg	409 500	pg/mL
ms TNF- α	8 163	pg	32 650	pg/mL
ms GM-CSF	1 300	pg	5 200	pg/mL
ms IL-13	22 500	pg	90 000	pg/mL
ms EPO	15 875	pg	63 500	pg/mL
ms VEGF-A	14 000	pg	56 000	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>	01 NOV 2018