



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0063

Expiration Date: 30 JUN 2023

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
hu IFN- γ	34 250	pg	137 000	pg/mL
hu IL-1 β	4 538	pg	18 150	pg/mL
hu IL-2	2 538	pg	10 150	pg/mL
hu IL-4	2 550	pg	10 200	pg/mL
hu IL-6	2 575	pg	10 300	pg/mL
hu IL-8	2 513	pg	10 050	pg/mL
hu IL-10	4 513	pg	18 050	pg/mL
hu IL-12p70	8 300	pg	33 200	pg/mL
hu IL-13	3 050	pg	12 200	pg/mL
hu TNF- α	4 563	pg	18 250	pg/mL
hu IL-5	4 888	pg	19 550	pg/mL
hu GM-CSF	11 963	pg	47 850	pg/mL
hu IL-17A	23 750	pg	95 000	pg/mL
hu VEGF-A	7 238	pg	28 950	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