



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0079

Expiration Date: 30 NOV 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	32 500	pg	130 000	pg/mL
hu IL-1 β	4 663	pg	18 650	pg/mL
hu IL-2	2 550	pg	10 200	pg/mL
hu IL-4	2 625	pg	10 500	pg/mL
hu IL-6	2 600	pg	10 400	pg/mL
hu IL-8	2 563	pg	10 250	pg/mL
hu IL-10	4 625	pg	18 500	pg/mL
hu IL-12p70	8 313	pg	33 250	pg/mL
hu IL-13	2 613	pg	10 450	pg/mL
hu TNF- α	4 563	pg	18 250	pg/mL
hu IL-5	5 075	pg	20 300	pg/mL
hu GM-CSF	12 150	pg	48 600	pg/mL
hu IL-17A	24 625	pg	98 500	pg/mL
hu VEGF-A	7 975	pg	31 900	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>	20 NOV 2018