



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0093

Expiration Date: 31 JAN 2024

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	34 875	pg	139 500	pg/mL
hu IL-1 β	4 788	pg	19 150	pg/mL
hu IL-2	2 688	pg	10 750	pg/mL
hu IL-4	2 700	pg	10 800	pg/mL
hu IL-6	2 650	pg	10 600	pg/mL
hu IL-8	2 575	pg	10 300	pg/mL
hu IL-10	4 625	pg	18 500	pg/mL
hu IL-12p70	9 063	pg	36 250	pg/mL
hu IL-13	3 238	pg	12 950	pg/mL
hu TNF- α	4 538	pg	18 150	pg/mL
hu IL-5	5 125	pg	20 500	pg/mL
hu GM-CSF	12 500	pg	50 000	pg/mL
hu IL-17A	47 375	pg	189 500	pg/mL
hu VEGF-A	7 813	pg	31 250	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>	04 FEB 2019