



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0085

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 EPO	25 250	pg	101 000	pg/mL
hu FLT-3L	7 700	pg	30 800	pg/mL
hu IFN-β	112 500	pg	450 000	pg/mL
hu IL-3	17 750	pg	71 000	pg/mL
hu IL-9	1 675	pg	6 700	pg/mL
hu IL-17B	3 988	pg	15 950	pg/mL
hu IL-17C	26 375	pg	105 500	pg/mL
hu IL-17D	57 875	pg	231 500	pg/mL
hu IL-1RA	5 613	pg	22 450	pg/mL
hu IL-2Rα	71 375	pg	285 500	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>	07 DEC 2018