



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0055

Expiration Date: 30 APR 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	28 375	pg	113 500	pg/mL
hu FLT3L	8 125	pg	32 500	pg/mL
hu IFN-β	131 250	pg	525 000	pg/mL
hu IL-3	21 375	pg	85 500	pg/mL
hu IL-9	2 088	pg	8 350	pg/mL
hu IL-17B	5 338	pg	21 350	pg/mL
hu IL-17C	30 500	pg	122 000	pg/mL
hu IL-17D	57 625	pg	230 500	pg/mL
hu IL-1RA	6 850	pg	27 400	pg/mL
hu IL-2Rα	70 875	pg	283 500	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>	29 AUG 2018