



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0064

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 EPO	63 625	pg	254 500	pg/mL
hu FLT-3L	7 638	pg	30 550	pg/mL
hu IFN-β	116 000	pg	464 000	pg/mL
hu IL-3	18 375	pg	73 500	pg/mL
hu IL-9	1 863	pg	7 450	pg/mL
hu IL-17B	4 475	pg	17 900	pg/mL
hu IL-17C	25 000	pg	100 000	pg/mL
hu IL-17D	57 375	pg	229 500	pg/mL
hu IL-1RA	5 788	pg	23 150	pg/mL
hu IL-2Rα	66 375	pg	265 500	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>	30 AUG 2018