



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 6

Catalog Number: C0072-2

Lot Number: A00U0087

Expiration Date: 31 OCT 2023

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
hu IL-17A/F	18 000	pg	72 000	pg/mL
hu-17E/IL-25	9 650	pg	38 600	pg/mL
hu IL-17F	123 500	pg	494 000	pg/mL
hu IL-21	14 250	pg	57 000	pg/mL
hu IL-22	3 675	pg	14 700	pg/mL
hu IL-23	26 125	pg	104 500	pg/mL
hu IL-27	50 125	pg	200 500	pg/mL
hu IL-29/IFN-λ1	11 225	pg	44 900	pg/mL
hu IL-31	11 838	pg	47 350	pg/mL
hu IL-33	11 288	pg	45 150	pg/mL
hu TSLP	10 050	pg	40 200	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>	16 NOV 2018