



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 6

Catalog Number: C0072-2

Lot Number: A00U0057

Expiration Date: 31 JUL 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	18 750	pg	75 000	pg/ml
hu IL-17E/IL-25	10 238	pg	40 950	pg/ml
hu IL-17F	128 750	pg	515 000	pg/ml
hu IL-21	16 500	pg	66 000	pg/ml
hu IL-22	6 538	pg	26 150	pg/ml
hu IL-23	25 125	pg	100 500	pg/ml
hu IL-27	59 625	pg	238 500	pg/ml
hu IL-29/IFN- λ1	14 250	pg	57 000	pg/ml
hu IL-31	12 625	pg	50 500	pg/ml
hu IL-33	11 513	pg	46 050	pg/ml
hu TSLP	9 163	pg	36 650	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>	17 AUG 2018