



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 6

**Catalog Number:** C0072-2

**Lot Number:** A00U0043

**Expiration Date:** 31 OCT 2022

**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	20 000	pg	80 000	pg/mL
hu-17E/IL-25	12 075	pg	48 300	pg/mL
hu IL-17F	115 250	pg	461 000	pg/mL
hu IL-21	16 375	pg	65 500	pg/mL
hu IL-22	4 488	pg	17 950	pg/mL
hu IL-23	26 250	pg	105 000	pg/mL
hu IL-27	66 875	pg	267 500	pg/mL
hu IL-29/IFN-λ1	14 000	pg	56 000	pg/mL
hu IL-31	12 413	pg	49 650	pg/mL
hu IL-33	13 250	pg	53 000	pg/mL
hu TSLP	9 900	pg	39 600	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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	29 AUG 2018