



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0044

Expiration Date: 30 NOV 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-1α	6 625	pg	26 500	pg/mL
hu IL-7	7 463	pg	29 850	pg/mL
hu IL-12/IL-23p40	26 125	pg	104 500	pg/mL
hu IL-15	3 613	pg	14 450	pg/mL
hu IL-16	25 125	pg	100 500	pg/mL
hu TNF-β	4 900	pg	19 600	pg/mL
hu G-CSF	27 750	pg	111 000	pg/mL
hu IFN-α2a	53 250	pg	213 000	pg/mL
hu IL-18	52 875	pg	211 500	pg/mL
hu TPO	50 000	pg	200 000	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