



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0035

Expiration Date: 28 FEB 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	6 863	pg	27 450	pg/mL
hu IL-7	9 288	pg	37 150	pg/mL
hu IL-12/IL-23p40	29 000	pg	116 000	pg/mL
hu IL-15	3 988	pg	15 950	pg/mL
hu IL-16	28 375	pg	113 500	pg/mL
hu TNF-β	6 200	pg	24 800	pg/mL
hu G-CSF	28 250	pg	113 000	pg/mL
hu IL-18	97 500	pg	390 000	pg/mL
hu TPO	53 000	pg	212 000	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