



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0083

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-1α	7 338	pg	29 350	pg/mL
hu IL-7	8 025	pg	32 100	pg/mL
hu IL-12/IL-23p40	28 500	pg	114 000	pg/mL
hu IL-15	3 863	pg	15 450	pg/mL
hu IL-16	26 375	pg	105 500	pg/mL
hu TNF-β	5 050	pg	20 200	pg/mL
hu G-CSF	26 875	pg	107 500	pg/mL
hu IFN-α2a	53 125	pg	212 500	pg/mL
hu IL-18	50 500	pg	202 000	pg/mL
hu TPO	51 500	pg	206 000	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>	19 OCT 2018