



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 3

**Catalog Number:** C0062-2

**Lot Number:** A00U0061

**Expiration Date:** 31 MAY 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	7 250	pg	29 000	pg/mL
hu IL-7	8 213	pg	32 850	pg/mL
hu IL-12/IL-23p40	25 750	pg	103 000	pg/mL
hu IL-15	3 750	pg	15 000	pg/mL
hu IL-16	26 375	pg	105 500	pg/mL
hu TNF-β	5 213	pg	20 850	pg/mL
hu G-CSF	27 875	pg	111 500	pg/mL
hu IFN-α2a	56 000	pg	224 000	pg/mL
hu IL-18	48 875	pg	195 500	pg/mL
hu TPO	56 000	pg	224 000	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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	29 AUG 2018