



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1

**Catalog Number:** C0060-2

**Lot Number:** A00U0053

**Expiration Date:** 31 MAR 2023

**Storage Condition:** 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
hu IFN- $\gamma$	39 000	pg	156 000	pg/mL
hu IL-1 $\beta$	4 613	pg	18 450	pg/mL
hu IL-2	2 563	pg	10 250	pg/mL
hu IL-4	2 825	pg	11 300	pg/mL
hu IL-6	2 525	pg	10 100	pg/mL
hu IL-8	2 575	pg	10 300	pg/mL
hu IL-10	4 288	pg	17 150	pg/mL
hu IL-12p70	8 363	pg	33 450	pg/mL
hu IL-13	2 913	pg	11 650	pg/mL
hu TNF- $\alpha$	4 538	pg	18 150	pg/mL
hu IL-5	6 688	pg	26 750	pg/mL
hu GM-CSF	8 725	pg	34 900	pg/mL
hu IL-17A	24 125	pg	96 500	pg/mL
hu VEGF-A	9 125	pg	36 500	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