



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 10

**Catalog Number:** C0091-2

**Lot Number:** A00U0054

**Expiration Date:** 30 APR 2023

**Storage Condition:** 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
hu Eotaxin-2	9 150	pg	36 600	pg/mL
hu GRO-α	3 613	pg	14 450	pg/mL
hu I-309	3 775	pg	15 100	pg/mL
hu MCP-2	3 150	pg	12 600	pg/mL
hu MCP-3	7 275	pg	29 100	pg/mL
hu M-CSF	2 900	pg	11 600	pg/mL
hu MIF	39 250	pg	157 000	pg/mL
hu MIP-5	43 125	pg	172 500	pg/mL
hu TRAIL	17 500	pg	70 000	pg/mL
hu YKL-40	7 013	pg	28 050	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