



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 10

**Catalog Number:** C0091-2

**Lot Number:** A00U0078

**Expiration Date:** 31 AUG 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	7 463	pg	29 850	pg/mL
hu GRO- $\alpha$	3 000	pg	12 000	pg/mL
hu I-309	2 900	pg	11 600	pg/mL
hu MCP-2	2 700	pg	10 800	pg/mL
hu MCP-3	5 825	pg	23 300	pg/mL
hu M-CSF	2 450	pg	9 800	pg/mL
hu MIF	33 875	pg	135 500	pg/mL
hu MIP-5	38 250	pg	153 000	pg/mL
hu TRAIL	14 125	pg	56 500	pg/mL
hu YKL-40	6 075	pg	24 300	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>	21 SEP 2018