



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 4

Catalog Number: C0063-2

Lot Number: A00U0058

Expiration Date: 30 JUN 2023

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
hu I-TAC	2 450	pg	9 800	pg/mL
hu MIP-3β	2 375	pg	9 500	pg/mL
hu SDF-1α	120 125	pg	480 500	pg/mL
hu CTACK	4 713	pg	18 850	pg/mL
hu ENA-78	5 088	pg	20 350	pg/mL
hu MIP-3α	23 750	pg	95 000	pg/mL
hu Fractalkine	118 500	pg	474 000	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>	29 AUG 2018