



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 4

Catalog Number: C0063-2

Lot Number: A00U0095

Expiration Date: 29 FEB 2024

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	2 425	pg	9 700	pg/mL
hu MIP-3β	2 250	pg	9 000	pg/mL
hu SDF-1α	114 500	pg	458 000	pg/mL
hu CTACK	4 388	pg	17 550	pg/mL
hu ENA-78	5 013	pg	20 050	pg/mL
hu MIP-3α	22 500	pg	90 000	pg/mL
hu Fractalkine	118 750	pg	475 000	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		11 APR 2019