



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 4

Catalog Number: C0063-2

Lot Number: A00U0080

Expiration Date: 30 SEP 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	2 600	pg	10 400	pg/mL
hu MIP-3β	2 575	pg	10 300	pg/mL
hu SDF-1α	130 000	pg	520 000	pg/mL
hu CTACK	5 050	pg	20 200	pg/mL
hu ENA-78	5 438	pg	21 750	pg/mL
hu MIP-3α	25 000	pg	100 000	pg/mL
hu Fractalkine	127 500	pg	510 000	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>	06 NOV 2018