



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 4

Catalog Number: C0063-2

Lot Number: A00U0041

Expiration Date: 31 OCT 2022

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator (pg)	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL) (pg/mL)	Concentration of Standard 1 (after 5-fold Dilution) (pg/mL)
Human I-TAC	2 725	10 900	2 180
Human MIP-3β	2 513	10 050	2 010
Human SDF-1α	125 000	500 000	100 000
Human CTACK	5 300	21 200	4 240
Human ENA-78	5 588	22 350	4 470
Human MIP-3α	24 875	99 500	19 900
Human Fractalkine	133 750	535 000	107 000

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>	25 OCT 2017