



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 4

Catalog Number: C0063-2

Lot Number: A00U0005

Expiration Date: 31 MAR 2020

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	2 525	10 100	2 020
Human MIP-3 α	26 000	104 000	20 800
Human MIP-3 β	2 525	10 100	2 020
Human SDF-1 α	128 750	515 000	103 000
Human Fractalkine	226 000	904 000	180 800
Human CTACK	5 250	21 000	4 200
Human ENA-78	4 925	19 700	3 940

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>	27 FEB 2017