



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 4

**Catalog Number:** C0063-2

**Lot Number:** A00U0039

**Expiration Date:** 30 JUN 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	2 650	10 600	2 120
Human MIP-3α	24 875	99 500	19 900
Human MIP-3β	2 563	10 250	2 050
Human SDF-1α	127 500	510 000	102 000
Human Fractalkine	131 250	525 000	105 000
Human CTACK	5 113	20 450	4 090
Human ENA-78	5 313	21 250	4 250

**Statement:** The above calibrator lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	30 JUN 2017