



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 4

**Catalog Number:** C0063-2

**Lot Number:** A00U0280

**Expiration Date:** 30 JUN 2027

**Storage Temperature:** 2-8°C

**Intended Assay Species:** Human

### Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
I-TAC	1875	pg	7500	pg/mL
MIP-3 $\beta$	2063	pg	8250	pg/mL
SDF-1 $\alpha$	121125	pg	484500	pg/mL
CTACK	5263	pg	21050	pg/mL
ENA-78	4550	pg	18200	pg/mL
MIP-3 $\alpha$	28250	pg	113000	pg/mL
Fractalkine	122875	pg	491500	pg/mL

**Reconstitution Volume:** 250  $\mu$ L

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.

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		08 AUG 2022