



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 4

**Catalog Number:** C0063-2

**Lot Number:** A00U0255

**Expiration Date:** 31 MAY 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	1900	pg	7600	pg/ml
MIP-3 $\beta$	2050	pg	8200	pg/ml
SDF-1 $\alpha$	120500	pg	482000	pg/ml
CTACK	4000	pg	16000	pg/ml
ENA-78	4800	pg	19200	pg/ml
MIP-3 $\alpha$	28375	pg	113500	pg/ml
Fractalkine	118500	pg	474000	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		29 JUN 2022