



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 4  
**Catalog Number:** C0063-2  
**Lot Number:** A00U0360  
**Expiration Date:** 31 JAN 2029  
**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          | 2087.5                          | pg              | 8350  | pg/mL           |
| MIP-3 $\beta$  | 2450                            | pg              | 9800  | pg/mL           |
| SDF-1 $\alpha$ | 120750                          | pg              | 483000  | pg/mL           |
| CTACK          | 5300                            | pg              | 21200   | pg/mL           |
| ENA-78         | 4962.5                          | pg              | 19850   | pg/mL           |
| MIP-3 $\alpha$ | 30875                           | pg              | 123500  | pg/mL           |
| Fractalkine    | 120375                          | pg              | 481500  | 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  |  | 12 FEB 2024 |