



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 4

Catalog Number: C0063-2

Lot Number: A00U0209

Expiration Date: 31 MAR 2026

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	2013	pg	8050	pg/mL
MIP-3 β	2238	pg	8950	pg/mL
SDF-1 α	124750	pg	499000	pg/mL
CTACK	4125	pg	16500	pg/mL
ENA-78	4825	pg	19300	pg/mL
MIP-3 α	29625	pg	118500	pg/mL
Fractalkine	123000	pg	492000	pg/mL

Reconstitution Volume: 250 μ 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
Review/Approval	Karen Hamilla	Quality		08 SEP 2021