



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 4

Catalog Number: C0063-2

Lot Number: A00U0177

Expiration Date: 31 OCT 2025

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	2363	pg	9450	pg/mL
MIP-3 β	2413	pg	9650	pg/mL
SDF-1 α	131250	pg	525000	pg/mL
CTACK	5350	pg	21400	pg/mL
ENA-78	4988	pg	19950	pg/mL
MIP-3 α	31250	pg	125000	pg/mL
Fractalkine	121125	pg	484500	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		02 DEC 2020