



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 4
Catalog Number: C0063-2
Lot Number: A00U0436
Expiration Date: 30 APR 2030
Storage Temperature: 2-8°C
Intended Assay Species: Human, NHP


Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
I-TAC	2137.5	pg	8550	pg/mL
MIP-3 β	2275	pg	9100	pg/mL
SDF-1 α	137500	pg	550000	pg/mL
CTACK	4350	pg	17400	pg/mL
ENA-78	4612.5	pg	18450	pg/mL
MIP-3 α	23500	pg	94000	pg/mL
Fractalkine	128750	pg	515000	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		22 MAY 2025