



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 4
Catalog Number: C0063-2
Lot Number: A00U0124
Expiration Date: 31 AUG 2024
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	2 488	pg	9 950	pg/mL
MIP-3 β	2 025	pg	8 100	pg/mL
SDF-1 α	118 625	pg	474 500	pg/mL
CTACK	4 363	pg	17 450	pg/mL
ENA-78	4 688	pg	18 750	pg/mL
MIP-3 α	22 375	pg	89 500	pg/mL
Fractalkine	122 625	pg	490 500	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	Xiomara Talavera	Quality		30 SEP 2019