



CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 4  
**Catalog Number:** C0063-2  
**Lot Number:** A00U0381  
**Expiration Date:** 31 DEC 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	2375	pg	9500	pg/mL
MIP-3β	2375	pg	9500	pg/mL
SDF-1α	122625	pg	490500	pg/mL
CTACK	4750	pg	19000	pg/mL
ENA-78	4725	pg	18900	pg/mL
MIP-3α	30250	pg	121000	pg/mL
Fractalkine	128750	pg	515000	pg/mL

Reconstitution Volume: 250 µL

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		24 JAN 2025