



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 4  
**Catalog Number:** C0063-2  
**Lot Number:** A00U0308  
**Expiration Date:** 31 JAN 2028  
**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	2300	pg	9200	pg/mL
MIP-3β	2275	pg	9100	pg/mL
SDF-1α	114500	pg	458000	pg/mL
CTACK	4937.5	pg	19750	pg/mL
ENA-78	4612.5	pg	18450	pg/mL
MIP-3α	27500	pg	110000	pg/mL
Fractalkine	131250	pg	525000	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Xiomara Talavera	Quality		25 MAY 2023