



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 4
Catalog Number: C0063-2
Lot Number: A00U0289
Expiration Date: 31 OCT 2027
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	2262.5	pg	9050	pg/mL
MIP-3β	2375	pg	9500	pg/mL
SDF-1α	126250	pg	505000	pg/mL
CTACK	5037.5	pg	20150	pg/mL
ENA-78	4625	pg	18500	pg/mL
MIP-3α	28625	pg	114500	pg/mL
Fractalkine	132500	pg	530000	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
Review/Approval	Xiomara Talavera	Quality	<i>Talavera</i>	20 DEC 2022