



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 10
Catalog Number: C0091-2
Lot Number: A00U0293
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
Eotaxin-2	5787.5	pg	23150	pg/mL
GRO- α	2975	pg	11900	pg/mL
I-309	2737.5	pg	10950	pg/mL
MCP-2	1812.5	pg	7250	pg/mL
MCP-3	5000	pg	20000	pg/mL
M-CSF	2187.5	pg	8750	pg/mL
MIF	26375	pg	105500	pg/mL
MIP-5	24500	pg	98000	pg/mL
TRAIL	10450	pg	41800	pg/mL
YKL-40	5125	pg	20500	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 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	Nithya Santhanam	Quality		10 FEB 2023