



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 10
Catalog Number: C0091-2
Lot Number: A00U0333
Expiration Date: 30 SEP 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
Eotaxin-2	6262.5	pg	25050	pg/mL
GRO- α	3137.5	pg	12550	pg/mL
I-309	4725	pg	18900	pg/mL
MCP-2	2325	pg	9300	pg/mL
MCP-3	6400	pg	25600	pg/mL
M-CSF	2412.5	pg	9650	pg/mL
MIF	25750	pg	103000	pg/mL
MIP-5	43375	pg	173500	pg/mL
TRAIL	11112.5	pg	44450	pg/mL
YKL-40	5812.5	pg	23250	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	Karen Hamilla	Quality		21 NOV 2023