



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 10
Catalog Number: C0091-2
Lot Number: A00U0354
Expiration Date: 31 DEC 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	6012.5	pg	24050	pg/mL
GRO- α	2937.5	pg	11750	pg/mL
I-309	4175	pg	16700	pg/mL
MCP-2	2112.5	pg	8450	pg/mL
MCP-3	5400	pg	21600	pg/mL
M-CSF	2237.5	pg	8950	pg/mL
MIF	24875	pg	99500	pg/mL
MIP-5	38750	pg	155000	pg/mL
TRAIL	9762.5	pg	39050	pg/mL
YKL-40	6087.5	pg	24350	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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 JAN 2024