



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 10
Catalog Number: C0091-2
Lot Number: A00U0398
Expiration Date: 30 NOV 2029
Storage Temperature: 2-8°C
Intended Assay Species: Human, NHP

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin-2	6825	pg	27300	pg/mL
GRO-α	3137.5	pg	12550	pg/mL
I-309	4087.5	pg	16350	pg/mL
MCP-2	3050	pg	12200	pg/mL
MCP-3	6750	pg	27000	pg/mL
M-CSF	2437.5	pg	9750	pg/mL
MIF	25750	pg	103000	pg/mL
MIP-5	38875	pg	155500	pg/mL
TRAIL	12150	pg	48600	pg/mL
YKL-40	6725	pg	26900	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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 DEC 2024