



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 10
Catalog Number: C0091-2
Lot Number: A00U0425
Expiration Date: 31 MAR 2030
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	6912.5	pg	27650	pg/mL
GRO- α	3175	pg	12700	pg/mL
I-309	3262.5	pg	13050	pg/mL
MCP-2	2225	pg	8900	pg/mL
MCP-3	5562.5	pg	22250	pg/mL
M-CSF	2475	pg	9900	pg/mL
MIF	33000	pg	132000	pg/mL
MIP-5	35625	pg	142500	pg/mL
TRAIL	11875	pg	47500	pg/mL
YKL-40	6612.5	pg	26450	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		24 APR 2025