



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 10
Catalog Number: C0091-2
Lot Number: A00U0441
Expiration Date: 31 MAY 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	6900	pg	27600	pg/mL
GRO- α	3537.5	pg	14150	pg/mL
I-309	3437.5	pg	13750	pg/mL
MCP-2	2275	pg	9100	pg/mL
MCP-3	5862.5	pg	23450	pg/mL
M-CSF	2537.5	pg	10150	pg/mL
MIF	33625	pg	134500	pg/mL
MIP-5	37125	pg	148500	pg/mL
TRAIL	12300	pg	49200	pg/mL
YKL-40	6662.5	pg	26650	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		18 JUL 2025