



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 10
Catalog Number: C0091-2
Lot Number: A00U0491
Expiration Date: 31 JUL 2031
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	7675	pg	30700	pg/mL
GRO- α	3300	pg	13200	pg/mL
I-309	3100	pg	12400	pg/mL
MCP-2	2600	pg	10400	pg/mL
MCP-3	5600	pg	22400	pg/mL
M-CSF	2600	pg	10400	pg/mL
MIF	33500	pg	134000	pg/mL
MIP-5	35500	pg	142000	pg/mL
TRAIL	12250	pg	49000	pg/mL
YKL-40	6725	pg	26900	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		11 SEP 2026