



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 10

Catalog Number: C0091 -2

Lot Number: A00U0286

Expiration Date: 31 AUG 2027

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	5675	pg	22700	pg/mL
GRO- α	2550	pg	10200	pg/mL
I-309	2250	pg	9000	pg/mL
MCP-2	1950	pg	7800	pg/mL
MCP-3	2775	pg	11100	pg/mL
M-CSF	2095	pg	8380	pg/mL
MIF	26000	pg	104000	pg/mL
MIP-5	26625	pg	106500	pg/mL
TRAIL	10350	pg	41400	pg/mL
YKL-40	5150	pg	20600	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		04 NOV 2022