



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 10

Catalog Number: C0091-2

Lot Number: A00U0242

Expiration Date: 30 NOV 2026

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	6413	pg	25650	pg/mL
GRO- α	2613	pg	10450	pg/mL
I-309	2650	pg	10600	pg/mL
MCP-2	2088	pg	8350	pg/mL
MCP-3	4413	pg	17650	pg/mL
M-CSF	2263	pg	9050	pg/mL
MIF	28000	pg	112000	pg/mL
MIP-5	28250	pg	113000	pg/mL
TRAIL	11088	pg	44350	pg/mL
YKL-40	5963	pg	23850	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		06 APR 2022