



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 10

Catalog Number: C0091-2

Lot Number: A00U0163

Expiration Date: 31 JUL 2025

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	8475	pg	33900	pg/mL
GRO- α	3213	pg	12850	pg/mL
I-309	3588	pg	14350	pg/mL
MCP-2	2275	pg	9100	pg/mL
MCP-3	5838	pg	23350	pg/mL
M-CSF	2163	pg	8650	pg/mL
MIF	27875	pg	111500	pg/mL
MIP-5	36125	pg	144500	pg/mL
TRAIL	12163	pg	48650	pg/mL
YKL-40	5950	pg	23800	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		13 AUG 2020