



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 10

Catalog Number: C0091-2

Lot Number: A00U0125

Expiration Date: 30 SEP 2024

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	8913	pg	35650	pg/mL
GRO- α	3175	pg	12700	pg/mL
I-309	3575	pg	14300	pg/mL
MCP-2	2638	pg	10550	pg/mL
MCP-3	5950	pg	23800	pg/mL
M-CSF	2225	pg	8900	pg/mL
MIF	30125	pg	120500	pg/mL
MIP-5	39625	pg	158500	pg/mL
TRAIL	11450	pg	45800	pg/mL
YKL-40	6188	pg	24750	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	Xiomara Talavera	Quality		07 OCT 2019