



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 10

**Catalog Number:** C0091-2

**Lot Number:** A00U0116

**Expiration Date:** 31 JUL 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	9 100	pg	36 400	pg/mL
GRO- $\alpha$	3 175	pg	12 700	pg/mL
I-309	3 650	pg	14 600	pg/mL
MCP-2	2 575	pg	10 300	pg/mL
MCP-3	5 800	pg	23 200	pg/mL
M-CSF	2 150	pg	8 600	pg/mL
MIF	33 375	pg	133 500	pg/mL
MIP-5	38 250	pg	153 000	pg/mL
TRAIL	12 475	pg	49 900	pg/mL
YKL-40	5 950	pg	23 800	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		30 JUL 2019