



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 10

**Catalog Number:** C0091-2

**Lot Number:** A00U0097

**Expiration Date:** 31 MAR 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	9 513	pg	38 050	pg/mL
GRO- $\alpha$	3 175	pg	12 700	pg/mL
I-309	2 963	pg	11 850	pg/mL
MCP-2	2 638	pg	10 550	pg/mL
MCP-3	7 175	pg	28 700	pg/mL
M-CSF	2 738	pg	10 950	pg/mL
MIF	32 250	pg	129 000	pg/mL
MIP-5	40 000	pg	160 000	pg/mL
TRAIL	14 875	pg	59 500	pg/mL
YKL-40	5 988	pg	23 950	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>	Dawn Stollar	Quality		10 APR 2019