



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 10

**Catalog Number:** C0091-2

**Lot Number:** A00U0150

**Expiration Date:** 31 MAR 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     | 8275                            | pg              | 33100   | pg/mL           |
| GRO- $\alpha$ | 3050                            | pg              | 12200   | pg/mL           |
| I-309         | 3063                            | pg              | 12250   | pg/mL           |
| MCP-2         | 2350                            | pg              | 9400  | pg/mL           |
| MCP-3         | 6250                            | pg              | 25000   | pg/mL           |
| M-CSF         | 2188                            | pg              | 8750  | pg/mL           |
| MIF           | 29625                           | pg              | 118500  | pg/mL           |
| MIP-5         | 35750                           | pg              | 143000  | pg/mL           |
| TRAIL         | 12250                           | pg              | 49000   | pg/mL           |
| YKL-40        | 6588                            | pg              | 26350   | 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  |           | 27 APR 2020 |