



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 10

Catalog Number: C0091-2

Lot Number: A00U0222

Expiration Date: 31 OCT 2026

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 | 6588 | pg | 26350 | pg/mL |
| GRO- α | 2688 | pg | 10750 | pg/mL |
| I-309 | 2713 | pg | 10850 | pg/mL |
| MCP-2 | 2088 | pg | 8350 | pg/mL |
| MCP-3 | 4488 | pg | 17950 | pg/mL |
| M-CSF | 2213 | pg | 8850 | pg/mL |
| MIF | 27500 | pg | 110000 | pg/mL |
| MIP-5 | 27000 | pg | 108000 | pg/mL |
| TRAIL | 11213 | pg | 44850 | pg/mL |
| YKL-40 | 5725 | pg | 22900 | 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 | | 19 NOV 2021 |