



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 10

Catalog Number: C0091-2

Lot Number: A00U0211

Expiration Date: 30 APR 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 | 6975 | pg | 27900 | pg/mL |
| GRO- α | 3000 | pg | 12000 | pg/mL |
| I-309 | 2950 | pg | 11800 | pg/mL |
| MCP-2 | 2263 | pg | 9050 | pg/mL |
| MCP-3 | 4975 | pg | 19900 | pg/mL |
| M-CSF | 2238 | pg | 8950 | pg/mL |
| MIF | 28125 | pg | 112500 | pg/mL |
| MIP-5 | 31250 | pg | 125000 | pg/mL |
| TRAIL | 11613 | pg | 46450 | pg/mL |
| YKL-40 | 6263 | pg | 25050 | 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 | | 28 JUN 2021 |