



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Chemokine Panel 1 (human) Calibrator Blend

Catalog Number: C0047-2

Lot Number: A0080204

Expiration Date: 30 APR 2024

Storage Temperature: 2-8°C

Concentration Assignment:

| Analyte | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|----------------|---|-----------------|
| Eotaxin | 1790 | pg/mL |
| MIP-1 β | 1280 | pg/mL |
| Eotaxin-3 | 5710 | pg/mL |
| TARC | 1870 | pg/mL |
| IP-10 | 2370 | pg/mL |
| MIP-1 α | 1090 | pg/mL |
| IL-8 | 66 700 | pg/mL |
| MCP-1 | 544 | pg/mL |
| MDC | 9 350 | pg/mL |
| MCP-4 | 683 | pg/mL |

Reconstitution Volume: 1000 μ L

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|--------------|----------|-----------|-------------|
| Review/Approval | Dawn Stollar | Quality | | 22 MAY 2019 |