



CERTIFICATE OF ANALYSIS

Product Description: Chemokine Panel 1 (human) Calibrator Blend
Catalog Number: C0047-2
Lot Number: A0080369
Expiration Date: 30 JUN 2028
Storage Temperature: 2-8°C

Concentration Assignment:

| Analyte | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|------------|-------------------------------------------------|-----------------|
| Eotaxin | 1860 | pg/mL |
| MIP-1β | 1190 | pg/mL |
| Eotaxin-3 | 6230 | pg/mL |
| TARC/CCL17 | 1790 | pg/mL |
| IP-10 | 2220 | pg/mL |
| MIP-1α | 1170 | pg/mL |
| IL-8 (HA) | 74600 | pg/mL |
| MCP-1 | 524 | pg/mL |
| MDC | 6060 | pg/mL |
| MCP-4 | 718 | pg/mL |

Reconstitution Volume: 1000 µL

This product was manufactured and tested according to MSD documents. This lot meets MSD’s specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality | | 19 JAN 2024 |