



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (mouse) Control Pack

Catalog Number: C4245-1

Lot Number: CP000055

Expiration Date: 30 JUN 2022

Concentration of Materials Tested:

| Units = pg/mL | Assay | | | | | | | | |
|-------------------------|------------|-------------|-------------|----------------------|-------------|----------------|----------------------|-------------|-------------|
| | ms IL-9 | ms MCP-1 | ms IL-33 | ms IL-27p28/IL-30 | ms IL-15 | ms IL-17A/F | ms MIP-1 α | ms IP-10 | ms MIP-2 |
| Control 1 (A00C0504) | 3430 | 489 | 2250 | 8890 | 30 000 | 2170 | 585 | 883 | 563 |
| Control 2 (A00C0505) | 835 | 149 | 252 | 1310 | 4440 | 228 | 50.1 | 251 | 66.6 |
| Control 3 (A00C0506) | 219 | 64.9 | 33.7 | 193 | 795 | 20.3 | 4.41 | 82.7 | 10.0 |

Reconstitution Volume: 250 μ L

| Control Pack Components | | | |
|------------------------------------|------------|---------------------|-----------------|
| Description | Lot Number | Storage Temperature | Expiration Date |
| Cytokine Panel 1 (mouse) Control 1 | A00C0504 | 2-8°C | 30 JUN 2022 |
| Cytokine Panel 1 (mouse) Control 2 | A00C0505 | 2-8°C | 30 JUN 2022 |
| Cytokine Panel 1 (mouse) Control 3 | A00C0506 | 2-8°C | 30 JUN 2022 |

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

| | Name | Function | Signature | Date |
|------------------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 21 AUG 2017 |