

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse MCP-1 Kit

Kit Catalog Numbers: K152NN-Series

Kit Lot Number: K0081519

Expiration Date: 31 Aug 2022

Acceptance Data/ Percent Recovery Specifications:

| Spot | Name | Calibrator | | Cytokine Panel 1 (mouse) Control 1 | | Cytokine Panel 1 (mouse) Control 2 | | Cytokine Panel 1 (mouse) Control 3 | |
|------|-------|------------|-------|------------------------------------|-------|------------------------------------|-------|------------------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | MCP-1 | 80% - 120% | < 20% | 75% - 125% | < 20% | 75% - 125% | < 20% | 75% - 125% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|-------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | MCP-1 | 4.42 | 75% - 125% | < 20% | 325 | 75% - 125% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|-------|------------------------------------|------|-------|
| 1 | MCP-1 | Calibrator | 447 | pg/mL |
| | | Cytokine Panel 1 (mouse) Control 1 | 443 | pg/mL |
| | | Cytokine Panel 1 (mouse) Control 2 | 135 | pg/mL |
| | | Cytokine Panel 1 (mouse) Control 3 | 56 | pg/mL |

Controls tested may not be the same lot as shipped.

Standard Curve Data:

| MCP-1 - Spot 1 | | |
|----------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 447 | 339,000 | 1.9 |
| 112 | 48,900 | 8.1 |
| 27.9 | 8,950 | 5.1 |
| 6.98 | 2,110 | 4.0 |
| 1.75 | 627 | 3.8 |
| 0.437 | 273 | 4.2 |
| 0.109 | 193 | 3.0 |
| 0.00 | 160 | 3.4 |

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------------|---------------------|-----------------|
| Mouse MCP-1 Plate | Z0047366 | 2-8°C | 31 Jan 2023 |
| Anti-ms MCP-1 Antibody | D0081498 | 2-8°C | 30 Nov 2023 |
| Cytokine Panel 1 (mouse) Calibrator Blend | A0080218 | 2-8°C | 31 Oct 2024 |
| Diluent 41 | M0410033 | ≤ -10°C | 31 Oct 2022 |
| Diluent 45 | M0450033 | ≤ -10°C | 31 Aug 2022 |
| MSD Read Buffer T | Not Kit Specific | Not Kit Specific | N/A |

| Additional Components Included in V-PLEX Plus | |
|---|------------------|
| Plate Seals | Not Kit Specific |
| Wash Buffer (20X),100mL | Not Kit Specific |
| Cytokine Panel 1 (mouse) Control 1 | Not Kit Specific |
| Cytokine Panel 1 (mouse) Control 2 | Not Kit Specific |
| Cytokine Panel 1 (mouse) Control 3 | Not Kit Specific |

Statement:

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

| | Name | Function | Signature | Date |
|------------------------|---------------|----------|--|-------------|
| Review/Approval | Karen Hamilla | Quality |  | 10 SEP 2020 |