

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-Plex Plus Mouse MCP-1 Kit  
**Kit Lot Number:** K0082289  
**Expiration Date:** 30 Jun 2025  
**Kit Catalog Numbers:** K152NN-Series

**Precision and Accuracy 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 Calibrator:**


| Spot | Name  | Conc. | Units |
|------|-------|-------|-------|
| 1    | MCP-1 | 479   | pg/mL |

**Lot-Specific Components:**

| Description                                   | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| Mouse MCP-1 SECTOR Plate                      | Z0047862   | 2-8°C               | 31 Oct 2025     |
| Anti-ms-MCP-1 Antibody                        | D0081800   | 2-8°C               | 31 Jul 2025     |
| Cytokine Panel 1 (mouse) Kit Calibrator Blend | A0080353   | 2-8°C               | 29 Feb 2028     |
| Diluent 41                                    | M0410052   | ≤ -10°C             | 30 Sep 2025     |
| Diluent 45                                    | M0450051   | ≤ -10°C             | 30 Jun 2025     |

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

|                 | Name          | Function | Signature  | Date        |
|-----------------|---------------|----------|--|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 15 JUN 2023 |