

CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Kit Lot Number: K0081876

Expiration Date: 29 Feb 2024

Kit Catalog Numbers: K151NN-Series; K156NN-Series

Precision and Accuracy Specifications:

| Spot | Name | Calibrator | | Chemokine Panel 1(human) Control 1 | | Chemokine Panel 1(human) Control 2 | | Chemokine Panel 1(human) Control 3 | |
|------|-------|------------|-------|------------------------------------|-------|------------------------------------|-------|------------------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | MCP-1 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|-------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | MCP-1 | 1.09 | 80% - 120% | < 20% | 375 | 80% - 120% | < 20% |

Stock Concentration of Calibrator:


| Spot | Name | Conc. | Units |
|------|-------|-------|-------|
| 1 | MCP-1 | 546 | pg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Human MCP-1 Plate | Z0047600 | 2-8°C | 31 Mar 2024 |
| Anti-hu MCP-1 Antibody | D0081541 | 2-8°C | 29 Feb 2024 |
| Chemokine Panel 1 (human) Calibrator Blend | A0080306 | 2-8°C | 30 Sep 2026 |
| Diluent 43 | M0430046 | ≤ -10°C | 29 Feb 2024 |
| Diluent 3 | M0030110 | ≤ -10°C | 31 Mar 2024 |

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 |  | 20 JAN 2022 |