

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human MCP-1 Kit
Kit Catalog Numbers: K151NN-Series, K156NN-Series
Kit Lot Number: K0081407
Expiration Date: 31 May 2022

Acceptance Data/ Percent Recovery 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 Material Tested :

| Spot | Name | Item | Conc | Units |
|------|-------|-------------------------------------|------|-------|
| 1 | MCP-1 | Calibrator | 552 | pg/mL |
| | | Chemokine Panel 1 (human) Control 1 | 468 | pg/mL |
| | | Chemokine Panel 1 (human) Control 2 | 71.4 | pg/mL |
| | | Chemokine Panel 1 (human) Control 3 | 12.1 | 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 |
| 552 | 215,000 | 2.8 |
| 138 | 54,500 | 2.0 |
| 34.5 | 14,300 | 3.0 |
| 8.63 | 3,740 | 3.1 |
| 2.16 | 1,050 | 2.0 |
| 0.539 | 303 | 2.6 |
| 0.135 | 120 | 3.8 |
| 0.00 | 57 | 7.3 |

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------------|---------------------|-----------------|
| Human MCP-1 Plate | Z0047254 | 2-8°C | 30 Jun 2022 |
| Anti-hu MCP-1 Antibody | D0081388 | 2-8°C | 31 Mar 2023 |
| Chemokine Panel 1 (human) Calibrator Blend | A0080228 | 2-8°C | 31 Aug 2024 |
| Diluent 43 | M0430036 | ≤ -10°C | 31 May 2022 |
| Diluent 3 | M0030094 | ≤ -10°C | 31 May 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 |
| Chemokine Panel 1 (human) Control 1 | Not Kit Specific |
| Chemokine Panel 1 (human) Control 2 | Not Kit Specific |
| Chemokine Panel 1 (human) 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 |  | 15 JUN 2020 |