

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: K0081340
Expiration Date: 31 Jan 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 | 544 | 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 |
| 544 | 196,000 | 4.0 |
| 136 | 48,300 | 8.6 |
| 34.0 | 12,500 | 3.4 |
| 8.50 | 3,190 | 5.9 |
| 2.13 | 947 | 1.9 |
| 0.531 | 317 | 3.2 |
| 0.133 | 148 | 3.0 |
| 0.00 | 89 | 3.5 |

CERTIFICATE OF ANALYSIS

Components:

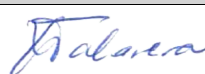
| Description | Lot Number | Storage Temperature | Expiration Date |
|--------------------------------------------|------------------|---------------------|-----------------|
| Human MCP-1 Plate | Z0047191 | 2-8°C | 31 Mar 2022 |
| Anti-hu MCP-1 Antibody | D0081124 | 2-8°C | 31 Jan 2022 |
| Chemokine Panel 1 (human) Calibrator Blend | A0080204 | 2-8°C | 30 Apr 2024 |
| Diluent 43 | M0430034 | ≤ -10°C | 28 Feb 2022 |
| Diluent 3 | M0030092 | ≤ -10°C | 31 Mar 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 | Xiomara Talavera | Quality |  | 12 DEC 2019 |