

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Kit
Kit Lot Number: K0082077
Expiration Date: 31 May 2024
Kit Catalog Numbers: K15047-Series; K15055-Series; K156NO-Series; K151NO-Series; K151NP-Series; K151NS-Series; K156NP-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 | Eotaxin | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |
| 2 | MIP-1β | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |
| 3 | Eotaxin-3 | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |
| 4 | TARC | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |
| 5 | IP-10 | 80% - 120% | < 20% | 75% - 125% | < 20% | 75% - 125% | < 20% | 75% - 125% | < 20% |
| 6 | MIP-1α | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |
| 7 | IL-8 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |
| 8 | MCP-1 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |
| 9 | MDC | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |
| 10 | MCP-4 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|-----------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | Eotaxin | 12.3 | 80% - 120% | < 20% | 1120 | 80% - 120% | < 20% |
| 2 | MIP-1β | 1.88 | 80% - 120% | < 20% | 520 | 80% - 120% | < 20% |
| 3 | Eotaxin-3 | 10.2 | 80% - 120% | < 20% | 3750 | 80% - 120% | < 20% |
| 4 | TARC | 3.32 | 80% - 120% | < 20% | 1120 | 75% - 125% | < 20% |
| 5 | IP-10 | 1.37 | 80% - 120% | < 20% | 500 | 80% - 120% | < 20% |
| 6 | MIP-1α | 13.8 | 80% - 120% | < 20% | 743 | 80% - 120% | < 20% |
| 7 | IL-8 | 713 | 75% - 125% | < 20% | 43400 | 75% - 125% | < 20% |
| 8 | MCP-1 | 1.09 | 80% - 120% | < 20% | 375 | 80% - 120% | < 20% |
| 9 | MDC | 88.3 | 80% - 120% | < 20% | 3700 | 75% - 125% | < 20% |
| 10 | MCP-4 | 1.49 | 80% - 120% | < 20% | 472 | 80% - 120% | < 20% |

Non-Specific Binding Specifications:

| Spot | Name | NSB Spec |
|------|-----------|----------|
| 1 | Eotaxin | < 1.0% |
| 2 | MIP-1β | < 1.0% |
| 3 | Eotaxin-3 | < 1.0% |
| 4 | TARC | < 1.0% |
| 5 | IP-10 | < 1.0% |
| 6 | MIP-1α | < 1.0% |
| 7 | IL-8 | < 1.0% |
| 8 | MCP-1 | < 1.0% |
| 9 | MDC | < 1.0% |
| 10 | MCP-4 | < 1.0% |

CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:


| Spot | Name | Conc. | Units |
|------|----------------|-------|-------|
| 1 | Eotaxin | 1960 | pg/mL |
| 2 | MIP-1 β | 1150 | pg/mL |
| 3 | Eotaxin-3 | 6460 | pg/mL |
| 4 | TARC | 1990 | pg/mL |
| 5 | IP-10 | 2510 | pg/mL |
| 6 | MIP-1 α | 1160 | pg/mL |
| 7 | IL-8 | 70600 | pg/mL |
| 8 | MCP-1 | 558 | pg/mL |
| 9 | MDC | 5880 | pg/mL |
| 10 | MCP-4 | 734 | pg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|----------------------------|-----------------|
| Chemokine Panel 1 (human) Plate | Z0047502 | 2-8°C | 31 Dec 2024 |
| Anti-hu Eotaxin Antibody | D0081765 | 2-8°C | 30 Apr 2025 |
| Anti-hu MIP-1 β Antibody | D0081872 | 2-8°C | 31 Dec 2025 |
| Anti-hu Eotaxin-3 Antibody | D0081677 | 2-8°C | 30 Sep 2024 |
| Anti-hu TARC Antibody | D0081767 | 2-8°C | 30 Sep 2025 |
| Anti-hu IP-10 Antibody | D0081809 | 2-8°C | 31 Jul 2025 |
| Anti-hu MIP-1 α Antibody | D0081835 | 2-8°C | 30 Sep 2025 |
| Anti-hu IL-8 (HA) Antibody | D0081810 | 2-8°C | 31 Jul 2025 |
| Anti-hu MCP-1 Antibody | D0081811 | 2-8°C | 31 Jul 2025 |
| Anti-hu MDC Antibody | D0081836 | 2-8°C | 30 Sep 2025 |
| Anti-hu MCP-4 Antibody | D0081837 | 2-8°C | 30 Nov 2025 |
| Chemokine Panel 1 (human) Calibrator Blend | A0080320 | 2-8°C | 31 Mar 2027 |
| Diluent 3 | M0030114 | $\leq -10^{\circ}\text{C}$ | 30 Sep 2024 |
| Diluent 43 | M0430047 | $\leq -10^{\circ}\text{C}$ | 31 May 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 |  | 02 SEP 2022 |