



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus Chemokine Panel 1 (hu) Gen. B Kit, QuickPlex  
**Lot Number:** K0082638  
**Expiration Date:** 31 MAY 2027  
**Catalog Number:** K156ANJ-Series, K15705-Series, K156ANK-Series, K151ANQ-Series, K156ANR-Series, K156ANL-Series, K156ANM-Series, K156ANP-Series, K15712-Series, K156ANQ-Series, K151ANR-Series, K156ANG-Series, K156ANH-Series, K151ANG-Series, K151ANH-Series, K151ANJ-Series, K151ANK-Series, K151ANL-Series, K151ANM-Series, K151ANP-Series

**Lot-Specific Components:**

| Description                                                  | Lot Number | Storage Temperature        | Expiration Date |
|--------------------------------------------------------------|------------|----------------------------|-----------------|
| Chemokine Panel 1 (hu) 96-Well 10-Spot QuickPlex Ultra Plate | Z0048072   | 2-8°C                      | 31 JUL 2027     |
| Chemokine Panel 1 (human) Gen. B Calibrator Blend            | A0080416   | 2-8°C                      | 31 DEC 2029     |
| Human Eotaxin Antibody                                       | D0082316   | 2-8°C                      | 31 JAN 2028     |
| Human Eotaxin-3 Antibody                                     | D0082397   | 2-8°C                      | 31 JUL 2028     |
| Human IP-10 Antibody                                         | D0082417   | 2-8°C                      | 31 JUL 2028     |
| Human MCP-1 Antibody                                         | D0082322   | 2-8°C                      | 31 JAN 2028     |
| Human MCP-4 Antibody                                         | D0082403   | 2-8°C                      | 31 JUL 2028     |
| Human MDC Antibody                                           | D0082402   | 2-8°C                      | 30 JUN 2028     |
| Human MIP-1 $\alpha$ Antibody                                | D0082321   | 2-8°C                      | 31 JAN 2028     |
| Human MIP-1 $\beta$ Antibody                                 | D0082317   | 2-8°C                      | 29 FEB 2028     |
| Human TARC Antibody                                          | D0082398   | 2-8°C                      | 30 JUN 2028     |
| Diluent 3                                                    | M0030137   | $\leq -10^{\circ}\text{C}$ | 31 JUL 2027     |
| Diluent 57                                                   | M0570018   | $\leq -10^{\circ}\text{C}$ | 31 MAY 2027     |

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

**Precision, Accuracy, and Sensitivity Specifications:**

| Spot | Name           | Calibrator |       | Chemokine Panel 1 (human)<br>Gen. B Control 1 |       | Chemokine Panel 1<br>(human) Gen. B Control<br>2 |       | Chemokine Panel 1<br>(human) Gen. B Control<br>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-1 $\beta$  | 80% - 120% | < 20% | 80% - 120%                                    | < 20% | 80% - 120%                                       | < 20% | 80% - 120%                                       | < 20% |
| 3    | Eotaxin-3      | 80% - 120% | < 20% | 75% - 125%                                    | < 20% | 75% - 125%                                       | < 20% | 75% - 125%                                       | < 20% |
| 4    | TARC/CCL17     | 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-1 $\alpha$ | 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% |



## CERTIFICATE OF ANALYSIS

| Spot | Name           | LLOQ  |            |       | ULOQ  |            |       |
|------|----------------|-------|------------|-------|-------|------------|-------|
|      |                | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | Eotaxin        | 2.14  | 80% - 120% | < 20% | 910   | 80% - 120% | < 20% |
| 2    | MIP-1 $\beta$  | 1.20  | 80% - 120% | < 20% | 390   | 80% - 120% | < 20% |
| 3    | Eotaxin-3      | 11.1  | 80% - 120% | < 20% | 11100 | 80% - 120% | < 20% |
| 4    | TARC/CCL17     | 2.14  | 80% - 120% | < 20% | 481   | 75% - 125% | < 20% |
| 5    | IP-10          | 1.29  | 80% - 120% | < 20% | 1240  | 80% - 120% | < 20% |
| 6    | MIP-1 $\alpha$ | 0.357 | 80% - 120% | < 20% | 403   | 80% - 120% | < 20% |
| 8    | MCP-1          | 1.45  | 80% - 120% | < 20% | 247   | 80% - 120% | < 20% |
| 9    | MDC            | 6.43  | 80% - 120% | < 20% | 2930  | 75% - 125% | < 20% |
| 10   | MCP-4          | 0.518 | 80% - 120% | < 20% | 234   | 80% - 120% | < 20% |

## Non-Specific Binding Specifications:

| Spot | Name           | Spec   |
|------|----------------|--------|
| 1    | Eotaxin        | < 1.0% |
| 2    | MIP-1 $\beta$  | < 1.0% |
| 3    | Eotaxin-3      | < 1.0% |
| 4    | TARC/CCL17     | < 1.0% |
| 5    | IP-10          | < 1.0% |
| 6    | MIP-1 $\alpha$ | < 1.0% |
| 8    | MCP-1          | < 1.0% |
| 9    | MDC            | < 1.0% |
| 10   | MCP-4          | < 1.0% |

## Stock Concentration of Calibrator:

| Spot | Name           | Conc. | Units |
|------|----------------|-------|-------|
| 1    | Eotaxin        | 1450  | pg/mL |
| 2    | MIP-1 $\beta$  | 583   | pg/mL |
| 3    | Eotaxin-3      | 21500 | pg/mL |
| 4    | TARC/CCL17     | 744   | pg/mL |
| 5    | IP-10          | 1860  | pg/mL |
| 6    | MIP-1 $\alpha$ | 618   | pg/mL |
| 8    | MCP-1          | 379   | pg/mL |
| 9    | MDC            | 4620  | pg/mL |
| 10   | MCP-4          | 358   | pg/mL |

Dilution to top of Curve: 1X

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

|                 | Name          | Function | Signature                                                                            | Date        |
|-----------------|---------------|----------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 11 APR 2025 |