

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Proinflammatory Panel 2 (rat) Kit**Lot Number:** K0082590**Expiration Date:** 31 AUG 2026**Catalog Number:** K15059-Series, K153QP-Series, K153QU-Series, K153QO-Series, K153QX-Series, K153QR-Series, K153QS-Series, K153QT-Series, K153OD-Series**Lot-Specific Components:**

| Description                                    | Lot Number | Storage Temperature        | Expiration Date |
|------------------------------------------------|------------|----------------------------|-----------------|
| Anti-rat IFN- $\gamma$ Antibody                | D0082368   | 2-8°C                      | 30 APR 2028     |
| Anti-rat IL-10 Antibody                        | D0082155   | 2-8°C                      | 31 MAR 2027     |
| Anti-rat IL-13 Antibody                        | D0082156   | 2-8°C                      | 31 MAY 2027     |
| Anti-rat IL-1 $\beta$ Antibody                 | D0082203   | 2-8°C                      | 31 AUG 2027     |
| Anti-rat IL-4 Antibody                         | D0082240   | 2-8°C                      | 31 DEC 2027     |
| Anti-rat IL-5 Antibody                         | D0082152   | 2-8°C                      | 31 MAR 2027     |
| Anti-rat IL-6 Antibody                         | D0082047   | 2-8°C                      | 30 SEP 2026     |
| Anti-rat KC/GRO Antibody                       | D0082154   | 2-8°C                      | 31 MAR 2027     |
| Anti-rat TNF- $\alpha$ Antibody                | D0082157   | 2-8°C                      | 31 MAR 2027     |
| Proinflammatory Panel 1 (rat) Calibrator Blend | A0080413   | 2-8°C                      | 30 APR 2030     |
| Proinflammatory Panel 2 (rat) Plate            | Z0048040   | 2-8°C                      | 30 APR 2027     |
| Diluent 40                                     | M0400021   | $\leq -10^{\circ}\text{C}$ | 31 OCT 2026     |
| Diluent 42                                     | M0420020   | $\leq -10^{\circ}\text{C}$ | 30 SEP 2026     |
| Blocker H                                      | R0100025   | Room Temperature           | 31 AUG 2026     |

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

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

| Spot | Name          | Calibrator |       | Proinflammatory Panel 1 (rat) Control 1 |       | Proinflammatory Panel 1 (rat) Control 2 |       | Proinflammatory Panel 1 (rat) Control 3 |       |
|------|---------------|------------|-------|-----------------------------------------|-------|-----------------------------------------|-------|-----------------------------------------|-------|
|      |               | Recovery   | CV    | Recovery                                | CV    | Recovery                                | CV    | Recovery                                | CV    |
| 1    | IFN- $\gamma$ | 80% - 120% | < 20% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% |
| 2    | IL-1 $\beta$  | 80% - 120% | < 20% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% |
| 4    | IL-4          | 80% - 120% | < 20% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% |
| 5    | IL-5          | 80% - 120% | < 20% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% |
| 6    | IL-6          | 80% - 120% | < 20% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% |
| 7    | KC/GRO        | 80% - 120% | < 20% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% |
| 8    | IL-10         | 80% - 120% | < 20% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% |
| 9    | IL-13         | 80% - 120% | < 20% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% |
| 10   | TNF- $\alpha$ | 80% - 120% | < 20% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% | 75% - 125%                              | < 25% |



## CERTIFICATE OF ANALYSIS

| Spot | Name          | LLOQ  |            |       | ULOQ  |            |       |
|------|---------------|-------|------------|-------|-------|------------|-------|
|      |               | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | IFN- $\gamma$ | 39.7  | 75% - 125% | < 25% | 3750  | 75% - 125% | < 25% |
| 2    | IL-1 $\beta$  | 102   | 75% - 125% | < 25% | 8100  | 75% - 125% | < 25% |
| 4    | IL-4          | 8.00  | 75% - 125% | < 25% | 723   | 75% - 125% | < 25% |
| 5    | IL-5          | 82.0  | 75% - 125% | < 25% | 3000  | 75% - 125% | < 25% |
| 6    | IL-6          | 96.9  | 75% - 125% | < 25% | 8550  | 75% - 125% | < 25% |
| 7    | KC/GRO        | 21.7  | 75% - 125% | < 25% | 728   | 75% - 125% | < 25% |
| 8    | IL-10         | 163   | 75% - 125% | < 25% | 15700 | 75% - 125% | < 25% |
| 9    | IL-13         | 12.5  | 75% - 125% | < 25% | 1080  | 75% - 125% | < 25% |
| 10   | TNF- $\alpha$ | 9.10  | 75% - 125% | < 25% | 793   | 75% - 125% | < 25% |

## Non-Specific Binding Specifications:

| Spot | Name          | Spec   |
|------|---------------|--------|
| 1    | IFN- $\gamma$ | < 1.0% |
| 2    | IL-1 $\beta$  | < 1.0% |
| 4    | IL-4          | < 1.0% |
| 5    | IL-5          | < 1.0% |
| 6    | IL-6          | < 1.0% |
| 7    | KC/GRO        | < 1.0% |
| 8    | IL-10         | < 1.0% |
| 9    | IL-13         | < 1.0% |
| 10   | TNF- $\alpha$ | < 1.0% |

## Stock Concentration of Calibrator:

| Spot | Name          | Conc. | Units |
|------|---------------|-------|-------|
| 1    | IFN- $\gamma$ | 6300  | pg/mL |
| 2    | IL-1 $\beta$  | 11900 | pg/mL |
| 4    | IL-4          | 1050  | pg/mL |
| 5    | IL-5          | 9540  | pg/mL |
| 6    | IL-6          | 13400 | pg/mL |
| 7    | KC/GRO        | 2950  | pg/mL |
| 8    | IL-10         | 23900 | pg/mL |
| 9    | IL-13         | 1670  | pg/mL |
| 10   | TNF- $\alpha$ | 1330  | 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  |  | 30 JUN 2025 |