

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX and V-PLEX Plus Proinflammatory Panel 1 (human) Kit  
**Lot Number:** K0082745  
**Expiration Date:** 30 SEP 2027  
**Catalog Number:** K15049-Series; K15052-Series; K15053-Series; K15056-Series; K151QR-Series;  
 K151OD-Series; K15345-Series; K15348-Series; K15346-Series; K15349-Series

**Lot-Specific Components:**

| Description                                               | Lot Number | Storage Temperature        | Expiration Date |
|-----------------------------------------------------------|------------|----------------------------|-----------------|
| Anti-hu IFN- $\gamma$ Antibody                            | D0082440   | 2-8°C                      | 31 AUG 2028     |
| Anti-hu IL-10 Antibody                                    | D0082432   | 2-8°C                      | 31 AUG 2028     |
| Anti-hu IL-12p70 Antibody                                 | D0082342   | 2-8°C                      | 29 FEB 2028     |
| Anti-hu IL-13 Antibody                                    | D0082471   | 2-8°C                      | 31 OCT 2028     |
| Anti-hu IL-1 $\beta$ Antibody                             | D0082366   | 2-8°C                      | 31 MAR 2028     |
| Anti-hu IL-2 Antibody                                     | D0082412   | 2-8°C                      | 31 JUL 2028     |
| Anti-hu IL-4 Antibody                                     | D0082441   | 2-8°C                      | 31 AUG 2028     |
| Anti-hu IL-6 Antibody                                     | D0082442   | 2-8°C                      | 31 AUG 2028     |
| Anti-hu IL-8 Antibody                                     | D0082431   | 2-8°C                      | 31 AUG 2028     |
| Anti-hu TNF- $\alpha$ Antibody                            | D0082434   | 2-8°C                      | 31 AUG 2028     |
| Proinflammatory Panel 1 (hu) 96-Well 10-Spot SECTOR Plate | Z0048144   | 2-8°C                      | 31 DEC 2027     |
| Proinflammatory Panel 1 (human) Calibrator Blend          | A0080428   | 2-8°C                      | 30 APR 2030     |
| Diluent 2                                                 | M0020134   | $\leq -10^{\circ}\text{C}$ | 30 SEP 2027     |
| Diluent 3                                                 | M0030140   | $\leq -10^{\circ}\text{C}$ | 30 SEP 2027     |

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<br>(human) Control 1 |       | Proinflammatory Panel 1<br>(human) Control 2 |       | Proinflammatory Panel 1<br>(human) Control 3 |       |
|------|---------------|------------|-------|----------------------------------------------|-------|----------------------------------------------|-------|----------------------------------------------|-------|
|      |               | Recovery   | CV    | Recovery                                     | CV    | Recovery                                     | CV    | Recovery                                     | CV    |
| 1    | IFN- $\gamma$ | 80% - 120% | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% |
| 2    | IL-1 $\beta$  | 80% - 120% | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% |
| 3    | IL-2          | 80% - 120% | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% |
| 4    | IL-4          | 80% - 120% | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% |
| 5    | IL-6          | 80% - 120% | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% |
| 6    | IL-8          | 80% - 120% | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% |
| 7    | IL-10         | 80% - 120% | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% |
| 8    | IL-12p70      | 80% - 120% | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% |
| 9    | IL-13         | 80% - 120% | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% | 80% - 120%                                   | < 20% |
| 10   | TNF- $\alpha$ | 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    | IFN- $\gamma$ | 1.76  | 80% - 120% | < 20% | 938   | 80% - 120% | < 20% |
| 2    | IL-1 $\beta$  | 0.646 | 80% - 120% | < 20% | 375   | 80% - 120% | < 20% |
| 3    | IL-2          | 0.89  | 80% - 120% | < 20% | 938   | 80% - 120% | < 20% |
| 4    | IL-4          | 0.218 | 80% - 120% | < 20% | 158   | 80% - 120% | < 20% |
| 5    | IL-6          | 0.633 | 80% - 120% | < 20% | 488   | 80% - 120% | < 20% |
| 6    | IL-8          | 0.591 | 80% - 120% | < 20% | 375   | 80% - 120% | < 20% |
| 7    | IL-10         | 0.298 | 80% - 120% | < 20% | 233   | 80% - 120% | < 20% |
| 8    | IL-12p70      | 1.22  | 80% - 120% | < 20% | 315   | 80% - 120% | < 20% |
| 9    | IL-13         | 4.21  | 80% - 120% | < 20% | 353   | 80% - 120% | < 20% |
| 10   | TNF- $\alpha$ | 0.69  | 80% - 120% | < 20% | 248   | 80% - 120% | < 20% |

## Non-Specific Binding Specifications:

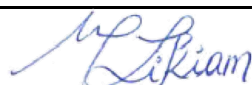
| Spot | Name          | Spec   |
|------|---------------|--------|
| 1    | IFN- $\gamma$ | < 1.0% |
| 2    | IL-1 $\beta$  | < 1.0% |
| 3    | IL-2          | < 1.0% |
| 4    | IL-4          | < 1.0% |
| 5    | IL-6          | < 1.0% |
| 6    | IL-8          | < 1.0% |
| 7    | IL-10         | < 1.0% |
| 8    | IL-12p70      | < 1.0% |
| 9    | IL-13         | < 1.0% |
| 10   | TNF- $\alpha$ | < 1.0% |

## Stock Concentration of Calibrator:

| Spot | Name          | Conc. | Units |
|------|---------------|-------|-------|
| 1    | IFN- $\gamma$ | 1490  | pg/mL |
| 2    | IL-1 $\beta$  | 605   | pg/mL |
| 3    | IL-2          | 1600  | pg/mL |
| 4    | IL-4          | 258   | pg/mL |
| 5    | IL-6          | 809   | pg/mL |
| 6    | IL-8          | 571   | pg/mL |
| 7    | IL-10         | 404   | pg/mL |
| 8    | IL-12p70      | 507   | pg/mL |
| 9    | IL-13         | 583   | pg/mL |
| 10   | TNF- $\alpha$ | 356   | 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/<br>Approval | Michelle Likiam | Quality  |  | 30 JUL 2025 |