

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX & V-PLEX Plus Proinflammatory Panel 1 (human) Kit**Lot Number:** K0082871**Expiration Date:** 30 NOV 2027**Catalog Number:** K15049-Series, K15052-Series, K15053-Series, K15056-Series, K151QR-Series, K151OD-Series, K15345-Series, K15348-Series, K15346-Series, K15349-Series, K156AR-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------|----------------------------|-----------------|
| Anti-hu IFN- γ Antibody | D0082502 | 2-8°C | 31 JAN 2029 |
| Anti-hu IL-10 Antibody | D0082432 | 2-8°C | 31 AUG 2028 |
| Anti-hu IL-12p70 Antibody | D0082433 | 2-8°C | 31 AUG 2028 |
| Anti-hu IL-13 Antibody | D0082471 | 2-8°C | 31 OCT 2028 |
| Anti-hu IL-1 β Antibody | D0082512 | 2-8°C | 30 JUN 2029 |
| Anti-hu IL-2 Antibody | D0082528 | 2-8°C | 31 MAR 2029 |
| Anti-hu IL-4 Antibody | D0082441 | 2-8°C | 31 AUG 2028 |
| Anti-hu IL-6 Antibody | D0082503 | 2-8°C | 31 JAN 2029 |
| Anti-hu IL-8 Antibody | D0082431 | 2-8°C | 31 AUG 2028 |
| Anti-hu TNF- α Antibody | D0082515 | 2-8°C | 28 FEB 2029 |
| Proinflammatory Panel 1 (hu) 96-Well 10-Spot SECTOR Plate | Z0048217 | 2-8°C | 30 JUN 2028 |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080434 | 2-8°C | 31 MAY 2030 |
| Diluent 2 | M0020135 | $\leq -10^{\circ}\text{C}$ | 30 NOV 2027 |
| Diluent 3 | M0030141 | $\leq -10^{\circ}\text{C}$ | 30 NOV 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 (human) Control 1 | | Proinflammatory Panel 1 (human) Control 2 | | Proinflammatory Panel 1 (human) Control 3 | |
|------|---------------|------------|-------|---|-------|---|-------|---|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | IFN- γ | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |
| 2 | IL-1 β | 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- α | 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- γ | 1.76 | 80% - 120% | < 20% | 938 | 80% - 120% | < 20% |
| 2 | IL-1 β | 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- α | 0.69 | 80% - 120% | < 20% | 248 | 80% - 120% | < 20% |

Non-Specific Binding Specifications:

| Spot | Name | Spec |
|------|---------------|--------|
| 1 | IFN- γ | < 1.0% |
| 2 | IL-1 β | < 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- α | < 1.0% |

Stock Concentration of Calibrator:

| Spot | Name | Conc. | Units |
|------|---------------|-------|-------|
| 1 | IFN- γ | 1170 | pg/mL |
| 2 | IL-1 β | 622 | pg/mL |
| 3 | IL-2 | 1560 | pg/mL |
| 4 | IL-4 | 276 | pg/mL |
| 5 | IL-6 | 780 | pg/mL |
| 6 | IL-8 | 555 | pg/mL |
| 7 | IL-10 | 381 | pg/mL |
| 8 | IL-12p70 | 494 | pg/mL |
| 9 | IL-13 | 596 | pg/mL |
| 10 | TNF- α | 346 | 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 | | 20 FEB 2026 |