

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 1 (human) Kit
Kit Lot Number: K0081913
Expiration Date: 31 Mar 2024
Kit Catalog Numbers: K151QR-Series; K151OD-Series; K15049-Series; K15052-Series; K15053-Series; K15346-Series; K15349-Series; K15345-Series; K15348-Series; K15056-Series

Precision and Accuracy 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% |

| 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 | 0.890 | 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 | 0.690 | 80% - 120% | < 20% | 248 | 80% - 120% | < 20% |

Non-Specific Binding Specifications:

| Spot | Name | NSB 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% |

CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:


| Spot | Name | Conc. | Units |
|------|---------------|-------|-------|
| 1 | IFN- γ | 1360 | pg/mL |
| 2 | IL-1 β | 649 | pg/mL |
| 3 | IL-2 | 1390 | pg/mL |
| 4 | IL-4 | 201 | pg/mL |
| 5 | IL-6 | 723 | pg/mL |
| 6 | IL-8 | 605 | pg/mL |
| 7 | IL-10 | 405 | pg/mL |
| 8 | IL-12p70 | 498 | pg/mL |
| 9 | IL-13 | 496 | pg/mL |
| 10 | TNF- α | 348 | pg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|----------------------------|-----------------|
| Proinflammatory Panel 1 (human) Plate | Z0047628 | 2-8°C | 30 Apr 2024 |
| Anti-hu IFN- γ Antibody | D0081745 | 2-8°C | 31 May 2025 |
| Anti-hu IL-1 β Antibody | D0081732 | 2-8°C | 31 Jan 2025 |
| Anti-hu IL-2 Antibody | D0081761 | 2-8°C | 31 Mar 2025 |
| Anti-hu IL-4 Antibody | D0081752 | 2-8°C | 31 Mar 2025 |
| Anti-hu IL-6 Antibody | D0081699 | 2-8°C | 28 Feb 2025 |
| Anti-hu IL-8 Antibody | D0081700 | 2-8°C | 30 Nov 2024 |
| Anti-hu IL-10 Antibody | D0081751 | 2-8°C | 30 Jun 2025 |
| Anti-hu IL-12p70 Antibody | D0081701 | 2-8°C | 28 Feb 2025 |
| Anti-hu IL-13 Antibody | D0081651 | 2-8°C | 30 Sep 2024 |
| Anti-hu TNF- α Antibody | D0081733 | 2-8°C | 31 Mar 2025 |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080310 | 2-8°C | 31 Oct 2026 |
| Diluent 2 | M0020119 | $\leq -10^{\circ}\text{C}$ | 31 Mar 2024 |
| Diluent 3 | M0030111 | $\leq -10^{\circ}\text{C}$ | 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 |  | 28 JAN 2022 |