

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Human IL-10 Kit**Lot Number:** K0082660**Expiration Date:** 31 DEC 2026**Catalog Number:** K151QU-Series; K156QU-Series**Lot-Specific Components:**

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Anti-hu IL-10 Antibody | D0082118 | 2-8°C | 31 JAN 2027 |
| Human IL-10 SECTOR Plate | Z0048089 | 2-8°C | 31 AUG 2027 |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080334 | 2-8°C | 31 JUL 2027 |
| Diluent 2 | M0020131 | ≤ -10°C | 31 DEC 2026 |
| Diluent 3 | M0030136 | ≤ -10°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 | | 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 | IL-10 | 80% - 120% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|-------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IL-10 | 0.298 | 80% - 120% | < 20% | 233 | 80% - 120% | < 20% |

Stock Concentration of Calibrator:

| Spot | Name | Conc. | Units |
|------|-------|-------|-------|
| 1 | IL-10 | 386 | 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 | | 03 APR 2025 |