



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IL-2 Kit
Lot Number: K0082499
Expiration Date: 30 JUN 2026
Catalog Number: K151QQ-Series, K156QQ-Series

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Anti-hu IL-2 Antibody | D0082163 | 2-8°C | 31 MAY 2027 |
| Human IL-2 Plate | Z0047970 | 2-8°C | 30 SEP 2026 |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080397 | 2-8°C | 31 MAY 2029 |
| Diluent 2 | M0020129 | ≤ -10°C | 30 JUN 2026 |
| Diluent 3 | M0030127 | ≤ -10°C | 30 SEP 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 (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-2 | 80% - 120% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IL-2 | 0.89 | 80% - 120% | < 20% | 938 | 80% - 120% | < 20% |

Stock Concentration of Calibrator:

| Spot | Name | Conc. | Units |
|------|------|-------|-------|
| 1 | IL-2 | 1470 | 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 | | 08 JUL 2024 |