

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Human VEGF Kit
Lot Number: K0082446
Expiration Date: 30 SEP 2026
Catalog Number: K151RH-Series, K156RH-Series

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| Anti-hu VEGF Antibody | D0082185 | 2-8°C | 30 JUN 2027 |
| Cytokine Panel 1 (human) Calibrator Blend | A0080331 | 2-8°C | 31 JUL 2027 |
| Human VEGF Plate | Z0057452 | 2-8°C | 30 SEP 2026 |
| Diluent 3 | M0030127 | ≤ -10°C | 30 SEP 2026 |
| Diluent 43 | M0430056 | ≤ -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 | | Cytokine Panel 1 (human) Control 1 | | Cytokine Panel 1 (human) Control 2 | | Cytokine Panel 1 (human) Control 3 | |
|------|------|------------|-------|------------------------------------|-------|------------------------------------|-------|------------------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | VEGF | 80% - 120% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | VEGF | 7.7 | 80% - 120% | < 20% | 562 | 80% - 120% | < 20% |

Stock Concentration of Calibrator:

| Spot | Name | Conc. | Units |
|------|------|-------|-------|
| 1 | VEGF | 1280 | 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 | | 17 MAY 2024 |