

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human VEGF Kit  
**Kit Lot Number:** K0081615  
**Expiration Date:** 31 Dec 2021  
**Kit Catalog Numbers:** K151RH-Series; K156RH-Series

**Precision and Accuracy 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% | 80% - 120%                         | < 20% | 80% - 120%                         | < 20% | 80% - 120%                         | < 20% |

| Spot | Name | LLOQ  |            |       | ULOQ  |            |       |
|------|------|-------|------------|-------|-------|------------|-------|
|      |      | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | VEGF | 7.70  | 80% - 120% | < 20% | 562   | 80% - 120% | < 20% |

**Stock Concentration of Calibrator:**


| Spot | Name | Conc. | Units |
|------|------|-------|-------|
| 1    | VEGF | 1080  | pg/mL |

**Lot-Specific Components:**

| Description                               | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| Human VEGF Plate                          | Z0056556   | 2-8°C               | 31 Dec 2021     |
| Anti-hu VEGF Antibody                     | D0081393   | 2-8°C               | 30 Apr 2023     |
| Cytokine Panel 1 (human) Calibrator Blend | A0080188   | 2-8°C               | 30 Nov 2023     |
| Diluent 43                                | M0430043   | ≤ -10°C             | 31 Jul 2023     |
| Diluent 3                                 | M0030104   | ≤ -10°C             | 30 Jun 2023     |

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  |  | 20 APR 2021 |