

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Angiogenesis Panel 1 (human) Kit

**Kit Lot Number:** K0081655

**Expiration Date:** 31 Dec 2022

**Kit Catalog Numbers:** K15190-Series; K151KW-Series; K151ME-Series; K151LT-Series; K151LU-Series; K151RZ-Series

**Precision and Accuracy Specifications:**

|      |        | Calibrator |       | Angiogenesis Control 1 |       | Angiogenesis Control 2 |       | Angiogenesis Control 3 |       |
|------|--------|------------|-------|------------------------|-------|------------------------|-------|------------------------|-------|
| Spot | Name   | Recovery   | CV    | Recovery               | CV    | Recovery               | CV    | Recovery               | CV    |
| 1    | VEGF-A | 80% - 120% | < 20% | 75% - 125%             | < 25% | 75% - 125%             | < 25% | 75% - 125%             | < 25% |
| 2    | VEGF-C | 80% - 120% | < 20% | 75% - 125%             | < 25% | 75% - 125%             | < 25% | 75% - 125%             | < 25% |
| 3    | VEGF-D | 80% - 120% | < 20% | 75% - 125%             | < 25% | 75% - 125%             | < 25% | 75% - 125%             | < 25% |
| 4    | Tie-2  | 80% - 120% | < 20% | 75% - 125%             | < 25% | 75% - 125%             | < 25% | 75% - 125%             | < 25% |
| 5    | Flt-1  | 80% - 120% | < 20% | 75% - 125%             | < 25% | 75% - 125%             | < 25% | 75% - 125%             | < 25% |
| 6    | PlGF   | 80% - 120% | < 20% | 75% - 125%             | < 25% | 75% - 125%             | < 25% | 75% - 125%             | < 25% |
| 7    | bFGF   | 80% - 120% | < 20% | 75% - 125%             | < 25% | 75% - 125%             | < 25% | 75% - 125%             | < 25% |

|      |        | LLOQ  |            |       | ULOQ  |            |       |
|------|--------|-------|------------|-------|-------|------------|-------|
| Spot | Name   | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | VEGF-A | 5.00  | 80% - 120% | < 20% | 1510  | 80% - 120% | < 20% |
| 2    | VEGF-C | 146   | 80% - 120% | < 20% | 17500 | 80% - 120% | < 20% |
| 3    | VEGF-D | 67.1  | 80% - 120% | < 20% | 18800 | 80% - 120% | < 20% |
| 4    | Tie-2  | 396   | 80% - 120% | < 20% | 63400 | 80% - 120% | < 20% |
| 5    | Flt-1  | 10.0  | 80% - 120% | < 25% | 6410  | 80% - 120% | < 20% |
| 6    | PlGF   | 1.50  | 80% - 120% | < 25% | 800   | 80% - 120% | < 25% |
| 7    | bFGF   | 2.60  | 80% - 120% | < 25% | 1780  | 80% - 120% | < 20% |

**Non-Specific Binding Specifications:**

| Spot | Name   | VEGF-A  | VEGF-C | VEGF-D | Tie-2  | VEGFR-1/Flt-1 | PlGF   | bFGF   |
|------|--------|---------|--------|--------|--------|---------------|--------|--------|
| 1    | VEGF-A | NA      | < 0.5% | < 0.5% | < 0.5% | < 10.0%       | < 0.5% | < 0.5% |
| 2    | VEGF-C | < 0.5%  | NA     | < 5.0% | < 0.5% | < 5.0%        | < 0.5% | < 0.5% |
| 3    | VEGF-D | < 0.5%  | < 0.5% | NA     | < 0.5% | < 0.5%        | < 0.5% | < 0.5% |
| 4    | Tie-2  | < 0.5%  | < 0.5% | < 0.5% | NA     | < 0.5%        | < 0.5% | < 0.5% |
| 5    | Flt-1  | < 10.0% | < 0.5% | < 0.5% | < 0.5% | NA            | < 0.5% | < 0.5% |
| 6    | PlGF   | < 0.5%  | < 0.5% | < 0.5% | < 0.5% | < 0.5%        | NA     | < 0.5% |
| 7    | bFGF   | < 0.5%  | < 0.5% | < 0.5% | < 0.5% | < 0.5%        | < 0.5% | NA     |

**CERTIFICATE OF ANALYSIS**

**Stock Concentration of Calibrator:**


| Spot | Name   | Conc.   | Units |
|------|--------|---------|-------|
| 1    | VEGF-A | 40200   | pg/mL |
| 2    | VEGF-C | 536000  | pg/mL |
| 3    | VEGF-D | 612000  | pg/mL |
| 4    | Tie-2  | 2020000 | pg/mL |
| 5    | Flt-1  | 204000  | pg/mL |
| 6    | PIGF   | 24000   | pg/mL |
| 7    | bFGF   | 54400   | pg/mL |

**Lot-Specific Components:**

| Description                                   | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| Angiogenesis Panel 1 (human) Plate            | Z0056875   | 2-8°C               | 31 Jul 2023     |
| Anti-hu VEGF Antibody                         | D0081494   | 2-8°C               | 30 Sep 2023     |
| Anti-hu VEGF-C Antibody                       | D0081517   | 2-8°C               | 31 Oct 2023     |
| Anti-hu VEGF-D Antibody                       | D0081576   | 2-8°C               | 31 Mar 2024     |
| Anti-hu Tie-2 Antibody                        | D0081577   | 2-8°C               | 30 Jun 2024     |
| Anti-hu Flt-1 Antibody                        | D0081578   | 2-8°C               | 31 Mar 2024     |
| Anti-hu PIGF Antibody                         | D0081492   | 2-8°C               | 31 Oct 2023     |
| Anti-hu bFGF Antibody                         | D0081493   | 2-8°C               | 30 Sep 2023     |
| Angiogenesis Panel 1 (human) Calibrator Blend | A0080260   | ≤ -70°C             | 31 Dec 2022     |
| Diluent 7                                     | M0070066   | ≤ -10°C             | 30 Jun 2023     |
| Diluent 11                                    | M0110048   | ≤ -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  |  | 21 JUL 2021 |