

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP VEGF Kit
Kit Catalog Numbers: K156RH-Series
Kit Lot Number: K0081032
Expiration Date: 31 May 2020

Acceptance Data/ Percent Recovery 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 | NHP VEGF | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

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

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|----------|------------------------------------|------|-------|
| 1 | NHP VEGF | Calibrator | 1490 | pg/mL |
| | | Cytokine Panel 1 (human) Control 1 | 410 | pg/mL |
| | | Cytokine Panel 1 (human) Control 2 | 140 | pg/mL |
| | | Cytokine Panel 1 (human) Control 3 | 50 | pg/mL |

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

Standard Curve Data:

| NHP VEGF - Spot 1 | | |
|-------------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 1490 | 331,000 | 1.5 |
| 373 | 62,800 | 1.5 |
| 93.1 | 13,000 | 3.1 |
| 23.3 | 3,100 | 3.7 |
| 5.82 | 819 | 2.0 |
| 1.46 | 244 | 3.3 |
| 0.364 | 106 | 3.2 |
| 0.00 | 59 | 6.8 |

CERTIFICATE OF ANALYSIS

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------------|---------------------|-----------------|
| Human VEGF Plate | Z0056286 | 2-8°C | 30 Nov 2020 |
| Anti-hu VEGF Antibody | D0081058 | 2-8°C | 31 May 2020 |
| Cytokine Panel 1 (human) Calibrator Blend | A0080144 | 2-8°C | 31 Aug 2022 |
| Diluent 43 | M0430026 | ≤ -10°C | 31 Aug 2020 |
| Diluent 3 | M0030082 | ≤ -10°C | 31 Oct 2020 |
| MSD Read Buffer T | Not Kit Specific | Not Kit Specific | N/A |

| Additional Components Included in V-PLEX Plus | |
|---|------------------|
| Plate Seals | Not Kit Specific |
| Wash Buffer (20X), 100mL | Not Kit Specific |
| Cytokine Panel 1 (human) Control 1 | Not Kit Specific |
| Cytokine Panel 1 (human) Control 2 | Not Kit Specific |
| Cytokine Panel 1 (human) Control 3 | Not Kit Specific |

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|--------------|-------------------|--|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance |  | 12 JUL 2018 |