

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP IL-15 Kit
Kit Catalog Numbers: K156RD-Series
Kit Lot Number: K0081036
Expiration Date: 31 Oct 2019

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 IL-15 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|-----------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | NHP IL-15 | 0.774 | 80% - 120% | < 20% | 525 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|-----------|------------------------------------|------|-------|
| 1 | NHP IL-15 | Calibrator | 796 | pg/mL |
| | | Cytokine Panel 1 (human) Control 1 | 288 | pg/mL |
| | | Cytokine Panel 1 (human) Control 2 | 46.3 | pg/mL |
| | | Cytokine Panel 1 (human) Control 3 | 8.87 | 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 IL-15 - Spot 1 | | |
|--------------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 796 | 234,000 | 1.8 |
| 199 | 39,200 | 2.1 |
| 49.8 | 9,350 | 1.6 |
| 12.4 | 2,350 | 1.3 |
| 3.11 | 664 | 1.8 |
| 0.777 | 215 | 2.9 |
| 0.194 | 99 | 3.9 |
| 0.00 | 61 | 5.3 |

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|-------------------------------------------|------------------|---------------------|-----------------|
| Human IL-15 Plate | Z0046856 | 2-8°C | 31 Dec 2020 |
| Anti-hu IL15 Antibody | D0080983 | 2-8°C | 31 Oct 2019 |
| Cytokine Panel 1 (human) Calibrator Blend | A0080073 | 2-8°C | 31 Jan 2021 |
| 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 | <i>Dawn Stollar</i> | 22 AUG 2018 |