

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® AND V-PLEX Plus Human IL-15 Kit
Kit Catalog Numbers: K156RD-Series, K151RD-Series
Kit Lot Number: K0081468
Expiration Date: 30 Sep 2021

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 | 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 | IL-15 | 0.774 | 80% - 120% | < 20% | 525 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|-------|------------------------------------|------|-------|
| 1 | IL-15 | Calibrator | 877 | pg/mL |
| | | Cytokine Panel 1 (human) Control 1 | 311 | pg/mL |
| | | Cytokine Panel 1 (human) Control 2 | 50.5 | pg/mL |
| | | Cytokine Panel 1 (human) Control 3 | 9.58 | pg/mL |

Controls tested may not be the same lot as shipped.

Standard Curve Data:

| IL-15 - Spot 1 | | |
|----------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 877 | 285,000 | 2.4 |
| 219 | 54,600 | 4.9 |
| 54.8 | 13,500 | 6.5 |
| 13.7 | 3,250 | 4.0 |
| 3.43 | 861 | 3.7 |
| 0.856 | 258 | 2.9 |
| 0.214 | 116 | 1.5 |
| 0.00 | 68 | 4.9 |

CERTIFICATE OF ANALYSIS


Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------------|---------------------|-----------------|
| Human IL-15 Plate | Z0047112 | 2-8°C | 30 Nov 2021 |
| Anti-hu IL-15 Antibody | D0081081 | 2-8°C | 30 Sep 2021 |
| Cytokine Panel 1 (human) Calibrator Blend | A0080188 | 2-8°C | 30 Nov 2023 |
| Dilent 43 | M0430035 | ≤ -10°C | 31 Mar 2022 |
| Dileunt 3 | M0030092 | ≤ -10°C | 31 Mar 2022 |
| MSD Read Buffer T | Not Kit Specific | Not Kit Specific | N/A |

| Additional Components Included in V-PLEX Plus | |
|---|------------------|
| MSD Wash Buffer (20X), 100mL | Not Kit Specific |
| Plate Seals | 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 | Karen Hamilla | Quality |  | 20 APR 2020 |