

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-Plex® and V-Plex Plus Human IL-1 α Kit
Kit Catalog Numbers: K151RB-Series
Kit Lot Number: K0081262
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-1 α | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|---------------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IL-1 α | 2.85 | 80% - 120% | < 20% | 278 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|---------------|------------------------------------|------|-------|
| 1 | IL-1 α | Calibrator | 454 | pg/mL |
| | | Cytokine Panel 1 (human) Control 1 | 183 | pg/mL |
| | | Cytokine Panel 1 (human) Control 2 | 58.1 | pg/mL |
| | | Cytokine Panel 1 (human) Control 3 | 16.3 | pg/mL |

Controls tested may not be the same lot as shipped.

Standard Curve Data:

| IL-1 α - Spot 1 | | |
|------------------------|------------|------|
| Conc. pg/mL | ECL Signal | %CV |
| 454 | 74,900 | 4.2 |
| 114 | 21,700 | 2.5 |
| 28.4 | 5,500 | 3.9 |
| 7.09 | 1,510 | 3.1 |
| 1.77 | 478 | 3.2 |
| 0.443 | 233 | 15.0 |
| 0.111 | 160 | 13.0 |
| 0.00 | 133 | 8.6 |

CERTIFICATE OF ANALYSIS


Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------------|---------------------|-----------------|
| Human IL-1 α Plate | Z0047110 | 2-8°C | 30 Nov 2021 |
| Anti-hu IL-1 α Antibody | D0081144 | 2-8°C | 28 Feb 2022 |
| Cytokine Panel 1 (human) Calibrator Blend | A0080188 | 2-8°C | 30 Nov 2023 |
| Diluent 43 | M0430032 | ≤ -10°C | 30 Sep 2021 |
| Diluent 3 | M0030089 | ≤ -10°C | 30 Sep 2021 |
| 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), 100 mL | 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 |  | 23 AUG 2019 |