

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-1 β Kit
Kit Catalog Numbers: K151QP-Series, K156QP-Series
Kit Lot Number: K0081373
Expiration Date: 31 Dec 2021

Acceptance Data/ Percent Recovery Specifications:

| Spot | Name | Calibrator | | Proinflammatory Panel 1 (human) Control 1 | | Proinflammatory Panel 1 (human) Control 2 | | Proinflammatory Panel 1 (human) Control 3 | |
|------|--------------|------------|-------|---|-------|---|-------|---|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | IL-1 β | 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 β | 2.14 | 80% - 120% | < 20% | 375 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|--------------|---|------|-------|
| 1 | IL-1 β | Calibrator | 543 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 1 | 218 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 2 | 51.8 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 3 | 11.9 | pg/mL |

Controls tested may not be the same lot as shipped.

Standard Curve Data:

| IL-1 β - Spot 1 | | |
|-----------------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 543 | 531,000 | 1.7 |
| 136 | 129,000 | 1.6 |
| 33.9 | 31,400 | 1.2 |
| 8.48 | 7,810 | 1.4 |
| 2.12 | 2,040 | 1.4 |
| 0.530 | 591 | 1.7 |
| 0.133 | 237 | 9.9 |
| 0.00 | 106 | 4.6 |

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------------|---------------------|-----------------|
| Human IL-1 β Plate | Z0047222 | 2-8°C | 30 Apr 2022 |
| Anti-hu IL-1 β Antibody | D0081128 | 2-8°C | 31 Jan 2022 |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080229 | 2-8°C | 30 Sep 2024 |
| Diluent 2 | M0020108 | \leq -10°C | 28 Feb 2022 |
| Diluent 3 | M0030090 | \leq -10°C | 31 Dec 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), 100mL | Not Kit Specific |
| Proinflammatory Panel 1 (human) Control 1 | Not Kit Specific |
| Proinflammatory Panel 1 (human) Control 2 | Not Kit Specific |
| Proinflammatory 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 | Carissa Hubbard | Quality |  | 23 DEC 2019 |