

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® AND V-PLEX Plus Human IL-6 Kit
Kit Catalog Numbers: K151QX-Series, K156QX-Series
Kit Lot Number: K0081317
Expiration Date: 28 Feb 2022

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-6 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IL-6 | 0.633 | 80% - 120% | < 20% | 488 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|------|---|------|-------|
| 1 | IL-6 | Calibrator | 755 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 1 | 267 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 2 | 47.7 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 3 | 7.49 | pg/mL |

Controls tested may not be the same lot as shipped.

Standard Curve Data:

| IL-6 - Spot 1 | | |
|---------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 755 | 388,000 | 1.9 |
| 189 | 88,000 | 3.6 |
| 47.2 | 21,100 | 3.1 |
| 11.8 | 5,210 | 3.0 |
| 2.95 | 1,340 | 5.3 |
| 0.737 | 381 | 2.1 |
| 0.184 | 136 | 3.5 |
| 0.00 | 56 | 5.9 |

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------------|---------------------|-----------------|
| Human IL-6 Plate | Z0047164 | 2-8°C | 28 Feb 2022 |
| Anti-hu IL-6 Antibody | D0081210 | 2-8°C | 30 Apr 2022 |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080216 | 2-8°C | 30 Jun 2024 |
| Diluent 2 | M0020109 | ≤ -10°C | 28 Feb 2022 |
| Diluent 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 | |
|---|------------------|
| 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 | Karen Hamilla | Quality |  | 05 DEC 2019 |