



CERTIFICATE OF ANALYSIS

Page 1 of 3

Product Description: V-PLEX® and V-PLEX Plus Human IL-12p70 Kit
Kit Catalog Numbers: K151QV-Series
Kit Lot Number: K0080842
Expiration Date: 31 Jul 2019

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 | hu IL-12p70 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|-------------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | hu IL-12p70 | 1.22 | 80% - 120% | < 20% | 315 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|-------------|-------------------------------------------|------|-------|
| 1 | hu IL-12p70 | Calibrator | 431 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 1 | 168 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 2 | 33.2 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 3 | 6.12 | pg/mL |

Controls tested may not be the same lot as shipped.



CERTIFICATE OF ANALYSIS

Page 2 of 3

Standard Curve Data:

| hu IL-12p70 - Spot 1 | | |
|----------------------|---------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 431 | 89,100 | 3.3 |
| 108 | 23,000 | 6.3 |
| 26.9 | 6,370 | 3.4 |
| 6.73 | 1,620 | 4.2 |
| 1.68 | 463 | 6.0 |
| 0.421 | 162 | 4.7 |
| 0.105 | 89 | 2.6 |
| 0.00 | 64 | 5.7 |

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--------------------------------------------------|------------------|---------------------|-----------------|
| Human IL-12p70 Plate | Z0046564 | 2-8°C | 31 Dec 2019 |
| Anti-hu IL-12p70 Antibody | D0080921 | 2-8°C | 31 Jul 2019 |
| Proinflammatory Panel 1 (Human) Calibrator Blend | A0080061 | 2-8°C | 30 Jun 2020 |
| Diluent 2 | M0020095 | ≤ -10°C | 30 Sep 2019 |
| Diluent 3 | M0030076 | ≤ -10°C | 30 Sep 2019 |
| 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.



CERTIFICATE OF ANALYSIS

Page 3 of 3

| | Name | Function | Signature | Date |
|-----------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 18 JUL 2017 |