

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP IP-10 Kit
Kit Catalog Numbers: K156NV-Series
Kit Lot Number: K0080999
Expiration Date: 30 Jun 2019

Acceptance Data/ Percent Recovery Specifications:

| Spot | Name | Calibrator | | Chemokine Panel 1 (human) Control 1 | | Chemokine Panel 1 (human) Control 2 | | Chemokine Panel 1 (human) Control 3 | |
|------|-----------|------------|-------|-------------------------------------|-------|-------------------------------------|-------|-------------------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | NHP IP-10 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|-----------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | NHP IP-10 | 1.37 | 80% - 120% | < 20% | 500 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|-----------|-------------------------------------|------|-------|
| 1 | NHP IP-10 | Calibrator | 2420 | pg/mL |
| | | Chemokine Panel 1 (human) Control 1 | 606 | pg/mL |
| | | Chemokine Panel 1 (human) Control 2 | 89.2 | pg/mL |
| | | Chemokine Panel 1 (human) Control 3 | 15.8 | pg/mL |

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

Standard Curve Data:

| NHP IP-10 - Spot 1 | | |
|--------------------|------------|------|
| Conc. pg/mL | ECL Signal | %CV |
| 2420 | 323,000 | 5.0 |
| 605 | 118,000 | 6.1 |
| 151 | 31,300 | 5.0 |
| 37.8 | 8,260 | 5.8 |
| 9.45 | 2,210 | 3.3 |
| 2.36 | 615 | 4.4 |
| 0.591 | 208 | 7.6 |
| 0.00 | 76 | 16.0 |

CERTIFICATE OF ANALYSIS

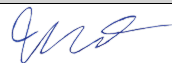
Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------------|---------------------|-----------------|
| Human IP-10 Plate | Z0046809 | 2-8°C | 30 Sep 2020 |
| Anti-hu IP-10 Antibody | D0080863 | 2-8°C | 30 Jun 2019 |
| Chemokine Panel 1 (human) Calibrator Blend | A0080131 | 2-8°C | 31 Jul 2022 |
| Diluent 43 | M0430024 | ≤ -10°C | 31 Jan 2020 |
| Diluent 3 | M0030081 | ≤ -10°C | 30 Sep 2020 |
| 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 |
| Chemokine Panel 1 (human) Control 1 | Not Kit Specific |
| Chemokine Panel 1 (human) Control 2 | Not Kit Specific |
| Chemokine 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 | Jessica Drolet | Quality Assurance |  | 14 May 2018 |