



CERTIFICATE OF ANALYSIS

COA-6069

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human MDC Kit
Kit Catalog Numbers: K151NPD-1, K151NPD-2, K151NPD-4, K151NPG-1, K151NPG-2, K151NPG-4
Kit Lot Number: K0080184
Expiration Date: 31 Dec 2016

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 | hu MDC | 80% - 120% | < 20% | 75% - 125% | < 20% | 75% - 125% | < 20% | 75% - 125% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|--------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | hu MDC | 88.3 | 80% - 120% | < 20% | 7500 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|--------|-------------------------------------|------|-------|
| 1 | hu MDC | Calibrator | 9220 | pg/mL |
| | | Chemokine panel 1 (human) Control 1 | 8670 | pg/mL |
| | | Chemokine panel 1 (human) Control 2 | 2660 | pg/mL |
| | | Chemokine panel 1 (human) Control 3 | 703 | pg/mL |

Controls tested may not be the same lot as shipped.



CERTIFICATE OF ANALYSIS

COA-6069

Page 2 of 2

Standard Curve Data:

| hu MDC - Spot 1 | | |
|-----------------|---------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 9220 | 454,000 | 3.7 |
| 2310 | 120,000 | 4.6 |
| 576 | 21,600 | 6.0 |
| 144 | 4,550 | 3.2 |
| 36.0 | 1,140 | 3.3 |
| 9.00 | 330 | 2.7 |
| 2.25 | 126 | 2.3 |
| 0.00 | 61 | 7.3 |

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------------|---------------------|-----------------|
| Human MDC Plate | Z0045319 | 2-8°C | 30 Nov 2016 |
| Anti-hu MDC Antibody | D0080119 | 2-8°C | 31 Dec 2015 |
| Chemokine panel 1 (human) Calibrator Blend | A0080038 | 2-8°C | 30 Sep 2016 |
| Diluent 43 | M0430009 | ≤ -10°C | 30 Nov 2016 |
| Diluent 3 | M0030061 | ≤ -10°C | 31 Aug 2016 |
| MSD Read Buffer T (4X) | 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 kit was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|-----------------|-------------|-------------------|-----------|-------------|
| Review/Approval | Monika Jain | Quality Assurance | | 15 OCT 2015 |




Meso Scale Diagnostics, L.L.C.

Description: V-PLEX® and V-PLEX Plus Human MDC Kit

Catalog Number(s): K151NP D/G-1/2/4

Lot #: K0080184

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of Lot # K0080184 has been extended to December 31, 2016.

| Function | Name | Signature | Date |
|-------------------|------------------|--|-------------|
| Quality Assurance | Michael Grossman |  | 14 OCT 2015 |

This memo supersedes all other statements of date of expiration for this kit, Lot # K0080184.