



CERTIFICATE OF ANALYSIS

COA-5857

Page 1 of 3

Product Description: V-PLEX® and V-PLEX Plus Human IL-8 Kit
Kit Catalog Numbers: K151RAD-1, K151RAD-2, K151RAD-4, K151RAG-1, K151RAG-2, K151RAG-4
Kit Lot Number: K0080391
Expiration Date: 30 Nov 2016

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-8 | 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-8 | 1.13 | 80% - 120% | < 20% | 375 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|---------|---|------|-------|
| 1 | hu IL-8 | Calibrator | 529 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 1 | 129 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 2 | 37.6 | pg/mL |
| | | Proinflammatory Panel 1 (human) Control 3 | 5.19 | pg/mL |

Controls tested may not be the same lot as shipped.



CERTIFICATE OF ANALYSIS

COA-5857

Page 2 of 3

Standard Curve Data:

| hu IL-8 - Spot 1 | | |
|------------------|---------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 529 | 238,501 | 4.7 |
| 132 | 61,893 | 3.2 |
| 33.1 | 14,784 | 5.2 |
| 8.27 | 3,703 | 7.0 |
| 2.07 | 1,019 | 2.7 |
| 0.517 | 291 | 3.8 |
| 0.129 | 116 | 9.4 |
| 0 | 54 | 6.4 |

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------------|---------------------|-----------------|
| Human IL-8 Plate | Z0044973 | 2-8°C | 29 Feb 2016 |
| Anti-hu IL-8 Antibody | D0080238 | 2-8°C | 30 Apr 2016 |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080031 | 2-8°C | 31 Mar 2016 |
| Diluent 2 | M0020079 | ≤ -10°C | 30 Nov 2015 |
| Diluent 3 | M0030059 | ≤ -10°C | 31 Mar 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 |
| 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 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.



CERTIFICATE OF ANALYSIS

COA-5857

Page 3 of 3

| | Name | Function | Signature | Date |
|------------------------|-------------|-------------------|--------------------|-------------|
| Review/Approval | Monika Jain | Quality Assurance | <i>Monika Jain</i> | 13 Aug 2015 |




Meso Scale Diagnostics, L.L.C.

Description: V-PLEX® and V-PLEX Plus Human IL-8 Kit

Catalog Number(s): K151RAD/G – 1/2/4

Lot #: K0080391

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of Lot # K0080391 has been extended to November 30, 2016.

| Function | Name | Signature | Date |
|-------------------|------------------|--|-------------|
| Quality Assurance | Michael Grossman |  | 13 AUG 2015 |

This memo supersedes all other statements of date of expiration for this kit, Lot # K0080391.