



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® Human MMP-2 Ultra-Sensitive Kit

**Catalog Number:** K151FYC-Series

**Lot Number:** K0035727

**Expiration Date:** N/A (See expiration date on specific components below)

**Statement:** This product was manufactured and tested according to approved MSD documents. This lot meets MSD’s specifications.

Kit Contents:

| Plate        |          |                  |         |                         |             |
|--------------|----------|------------------|---------|-------------------------|-------------|
| <b>Lot #</b> | Z0046822 | <b>Catalog #</b> | N45033A | <b>Expiration Date:</b> | 31 OCT 2020 |

| Antibody Blend              |                    |
|-----------------------------|--------------------|
| <b>Lot Number:</b>          | D0034040           |
| <b>Catalog Number:</b>      | D21FY              |
| <b>Expiration Date:</b>     | 30 NOV 2021        |
| <b>Storage Temperature:</b> | 2-8°C              |
| <b>Spot #</b>               | <b>Description</b> |
| 1                           | MMP-2              |

| Calibrator Blend            |             |
|-----------------------------|-------------|
| <b>Lot Number:</b>          | A0032396    |
| <b>Catalog Number:</b>      | C0033       |
| <b>Expiration Date:</b>     | 30 NOV 2020 |
| <b>Storage Temperature:</b> | ≤ -70°C     |

Specifications:

Release method:

| Coating Uniformity             | Specifications |
|--------------------------------|----------------|
| Mean Intraplate CV             | ≤ 15%          |
| Mean Interplate CV             | ≤ 18%          |
| No more than one plate with CV | > 18%          |
| No plate with CV               | > 20%          |

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

|                        | Name         | Function          | Signature | Date        |
|------------------------|--------------|-------------------|-----------|-------------|
| <b>Review/Approval</b> | Dawn Stollar | Quality Assurance |           | 03 MAR 2019 |