



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD<sup>®</sup> Human Angiopoietin-1 Kit

**Catalog Number:** K151LPD-Series

**Lot Number:** K00E0307

**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> | Z00E0246 | <b>Catalog #</b> | N451LPA | <b>Expiration Date:</b> | 31 MAR 2019 |

| Antibody Blend              |             |
|-----------------------------|-------------|
| <b>LotNumber:</b>           | D00E0286    |
| <b>Catalog Number:</b>      | D21LP       |
| <b>Expiration Date:</b>     | 30 APR 2019 |
| <b>Storage Temperature:</b> | 2-8°C       |

| Calibrator Blend            |             |
|-----------------------------|-------------|
| <b>Lot Number:</b>          | A00E0152    |
| <b>Catalog Number:</b>      | C01LP       |
| <b>Expiration Date:</b>     | 30 APR 2019 |
| <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 | <i>Dawn Stollar</i> | 22 MAY 2017 |