



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD[®] Human Adiponectin Kit

Catalog Number: K151BXC - Series

Lot Number: K0013754

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 | | | | | |
|-------|----------|-----------|---------|------------------|-------------|
| Lot # | Z0055252 | Catalog # | N451BXA | Expiration Date: | 30 JUN 2017 |

| Antibody Blend | |
|----------------------|-------------|
| Lot Number: | D0013343 |
| Catalog Number: | D21BX |
| Expiration Date: | 30 JUN 2017 |
| Storage Temperature: | 2-8°C |

| Calibrator Blend | |
|----------------------|-------------|
| Lot Number: | A0010376 |
| Catalog Number: | C01BX |
| Expiration Date: | 30 JUN 2017 |
| Storage Temperature: | ≤ -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 |
|-----------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 30 MAY 2017 |