



## CERTIFICATE OF ANALYSIS

COA - 07893

**Product Description:** MSD<sup>®</sup> Mouse Adiponectin Kit

**Catalog #/Size:** K152BXC - Series

**Lot Number:** K0014019

**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 Number: Z0055784

Plate Lot Catalog Number: L452BXB

Plate Lot Expiration Date: 31 MAY 2019

Antibody blend Lot Number: D0013626

Antibody blend Catalog Number: D22BX

Antibody blend Expiration Date: 31 JUL 2019

Calibrator blend Lot Number: A0010410

Calibrator blend Catalog Number: C02BX

Calibrator blend Expiration Date: 31 MAY 2019

### Specifications:

Release method: Coating uniformity

Mean Intraplate CV:  $\leq 15\%$

Mean Interplate CV:  $\leq 18\%$

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>Preparer</b>	Xiomara Talavera	Quality Control	<i>X Talavera</i>	21 FEB 2017
<b>Review/Approval</b>	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	22 FEB 2017