



## CERTIFICATE OF ANALYSIS

COA - 07201

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

**Catalog #/Size:** K152BZC-1 / 1 Plate; K152BZC-2 / 5 Plates; K152BZC-3 / 20 Plates

**Lot Number:** K0013918

**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: Z0055609

Plate Lot Catalog Number: L452BZA

Plate Lot Expiration Date: 31 OCT 2018

Antibody blend Lot Number: D0013521

Antibody blend Catalog Number: D22BZ

Antibody blend Expiration Date: 31 JUL 2018

Calibrator blend Lot Number: A0010387

Calibrator blend Catalog Number: C00BZ

Calibrator blend Expiration Date: 30 APR 2018

### 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>	10 AUG 2016
<b>Review/Approval</b>	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	11 AUG 2016