



## CERTIFICATE OF ANALYSIS

COA - 07093

**Product Description:** MSD<sup>®</sup> Hu Active GLP-1(7-36)amide, Insulin, Glucagon, Leptin Kit

**Catalog #/Size:** K15173C-1 / 1 Plate; K15173C-2 / 5 Plates; K15173C-3 / 20 Plates

**Lot Number:** K0013763

**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: Z0055265

Plate Lot Catalog Number: N45173A

Plate Lot Expiration Date: 31 JUL 2017

Antibody blend Lot Number: D0013350

Antibody blend Catalog Number: D20FA

Antibody blend Expiration Date: 31 JUL 2017

Antibody blend Lot Number: D0013352

Antibody blend Catalog Number: D21BZ

Antibody blend Expiration Date: 31 JUL 2017

Antibody blend Lot Number: D0013353

Antibody blend Catalog Number: D20HC

Antibody blend Expiration Date: 31 JUL 2017

Antibody blend Lot Number: D0013356

Antibody blend Catalog Number: D21BY

Antibody blend Expiration Date: 31 JUL 2017

Calibrator blend Lot Number: A0010368

Calibrator blend Catalog Number: C00FA

Calibrator blend Expiration Date: 31 MAR 2017

Calibrator blend Lot Number: A0010377

Calibrator blend Catalog Number: C00BZ

Calibrator blend Expiration Date: 30 JUN 2017

Calibrator blend Lot Number: A0010366

Calibrator blend Catalog Number: C00HC

Calibrator blend Expiration Date: 28 FEB 2017

Calibrator blend Lot Number: A0010367

Calibrator blend Catalog Number: C01BY

Calibrator blend Expiration Date: 31 MAR 2017



## CERTIFICATE OF ANALYSIS

### 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.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		08 JUL 2016
<b>Review/Approval</b>	Vanessa Castleberry	Quality Assurance		11 JUL 2016