



## CERTIFICATE OF ANALYSIS

COA - 06223

**Product Description:** MSD<sup>®</sup> Active GLP-1 (ver. 2) Kit

**Catalog #/Size:** K150JWC-1 / 1 Plate; K150JWC-2 / 5 Plates; K150JWC-4 / 25 Plates

**Lot Number:** K0040301

**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: Z0055346

Plate Lot Catalog Number: L450JWA

Plate Lot Expiration Date: 31 OCT 2017

Antibody blend Lot Number: D0040528

Antibody blend Catalog Number: D20JX

Antibody blend Expiration Date: 31 OCT 2017

Calibrator blend Lot Number: A0040114

Calibrator blend Catalog Number: C00JW

Calibrator blend Expiration Date: 30 APR 2017

### 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>	Rolanda Spratley	Biologics/Chemistry		23 NOV 2015
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		25 NOV 2015