



CERTIFICATE OF ANALYSIS

COA - 07765

Product Description: MSD[®] Active GLP-1 (ver. 2) Kit

Catalog #/Size: K150JWC - Series

Lot Number: K0040371

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: Z0055714

Plate Lot Catalog Number: L450JWA

Plate Lot Expiration Date: 28 FEB 2019

Antibody blend Lot Number: D0040650

Antibody blend Catalog Number: D20JX

Antibody blend Expiration Date: 28 FEB 2019

Calibrator blend Lot Number: A0040170

Calibrator blend Catalog Number: C00JW

Calibrator blend Expiration Date: 30 SEP 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
Preparer	Xiomara Talavera	Quality Control	<i>X Talavera</i>	17 JAN 2017
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	17 JAN 2017