



CERTIFICATE OF ANALYSIS

COA - 06347

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

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

Lot Number: K0040281

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: Z0055214
- Plate Lot Catalog Number: L450JWA
- Plate Lot Expiration Date: 31 MAY 2017
- Antibody blend Lot Number: D0040503
- Antibody blend Catalog Number: D20JX
- Antibody blend Expiration Date: 30 APR 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: ≤ 15%
- Mean Interplate CV: ≤ 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	Rolanda Spratley	Biologics/Chemistry		11 JAN 2016
Review/Approval	Dawn Stollar	Quality Assurance		11 JAN 2016