



CERTIFICATE OF ANALYSIS

COA - 06224

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

Catalog #/Size: K150JVC-1 / 1 Plate; K150JVC-2 / 5 Plates; K150JVC-4 / 25 Plates

Lot Number: K0040294

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: Z0055298

Plate Lot Catalog Number: L450JVA

Plate Lot Expiration Date: 31 AUG 2017

Antibody blend Lot Number: D0040516

Antibody blend Catalog Number: D20JY

Antibody blend Expiration Date: 31 AUG 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
Preparer	Rolanda Spratley	Biologics/Chemistry		23 NOV 2015
Review/Approval	Dawn Stollar	Quality Assurance		25 NOV 2015