



CERTIFICATE OF ANALYSIS

COA - 06346

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: K0040306

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: Z0055366

Plate Lot Catalog Number: L450JVA

Plate Lot Expiration Date: 30 NOV 2017

Antibody blend Lot Number: D0040538

Antibody blend Catalog Number: D20JY

Antibody blend Expiration Date: 31 JAN 2018

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 | | 17 DEC 2015 |
| Review/Approval | Dawn Stollar | Quality Assurance | | 18 DEC 2015 |