



RFE MEMO

Page 1 of 1

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

Catalog Number(s) / Size(s): K150JWC-Series

Lot Number: K0040400

Based on additional testing, the original date of expiration for this Kit has been extended by 12 months. The expiration date of K0040400 has been extended to 30 SEP 2019.

The expiration date for the following kit component has also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
GLP-1 (7-36)amide Calibrator	A0040170	30 SEP 2018	30 SEP 2019

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	04 APR 2018

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K150JWC - Series

Lot Number: K0040400

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 #	Z0055991	Catalog #	L450JWA	Expiration Date:	31 DEC 2019

Antibody Blend	
Lot Number:	D0040716
Catalog Number:	D20JX
Expiration Date:	31 DEC 2019
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0040170
Catalog Number:	C00JW
Expiration Date:	30 SEP 2018
Storage Temperature:	≤ -70°C

Specifications:

Release method:

Coating Uniformity	Specifications
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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	31 JAN 2018