



CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: MSD[®] Human Active GLP-1 (7-36)amide/Insulin/Glucagon/Leptin Kit

Catalog Number: K15173C - Series

Lot Number: K0013865

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 #	Z0055455	Catalog #	N45173A	Expiration Date:	28 FEB 2018

Antibody Blend	
Lot Number:	D0013454
Catalog Number:	D20FA
Expiration Date:	28 FEB 2018
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0010366
Catalog Number:	C00HC
Expiration Date:	28 FEB 2017
Storage Temperature:	≤ -70°C

Antibody Blend	
Lot Number:	D0013456
Catalog Number:	D21BZ
Expiration Date:	28 FEB 2018
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0010367
Catalog Number:	C01BY
Expiration Date:	31 MAR 2017
Storage Temperature:	≤ -70°C

Antibody Blend	
Lot Number:	D0013457
Catalog Number:	D20HC
Expiration Date:	28 FEB 2018
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0010368
Catalog Number:	C00FA
Expiration Date:	31 MAR 2017
Storage Temperature:	≤ -70°C

Antibody Blend	
Lot Number:	D0013458
Catalog Number:	D21BY
Expiration Date:	28 FEB 2018
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0010369
Catalog Number:	C00BZ
Expiration Date:	31 MAR 2017
Storage Temperature:	≤ -70°C



CERTIFICATE OF ANALYSIS

Page 2 of 2

Specifications:

Release method:

Coating Uniformity	Specifications
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
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	01JUN2017