



CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: MSD® Human Active GLP-1/Insulin/Glucagon/Leptin Kit

Catalog Number: K15174C - Series

Lot Number: K0014303

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 #	Z0056249	Catalog #	N45173A	Expiration Date:	31 OCT 2020

Antibody Blend	
Lot Number:	D0013931
Catalog Number:	D20FA
Expiration Date:	31 OCT 2021
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0010446
Catalog Number:	C00FA
Expiration Date:	30 SEP 2020
Storage Temperature:	≤ -70°C

Antibody Blend	
Lot Number:	D0013932
Catalog Number:	D20FB
Expiration Date:	31 DEC 2021
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0010433
Catalog Number:	C00BZ
Expiration Date:	30 APR 2020
Storage Temperature:	≤ -70°C

Antibody Blend	
Lot Number:	D0013933
Catalog Number:	D21BZ
Expiration Date:	31 OCT 2021
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0010451
Catalog Number:	C00HC
Expiration Date:	31 OCT 2020
Storage Temperature:	≤ -70°C

Antibody Blend	
Lot Number:	D0013934
Catalog Number:	D20HC
Expiration Date:	31 OCT 2021
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0010428
Catalog Number:	C01BY
Expiration Date:	30 NOV 2019
Storage Temperature:	≤ -70°C

Antibody Blend	
Lot Number:	D0013935
Catalog Number:	D21BY
Expiration Date:	31 OCT 2021
Storage Temperature:	2-8°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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	21 DEC 2018