



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® Human Leptin Kit

**Catalog Number:** K151BYC - Series

**Lot Number:** K0014274

**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					
<b>Lot #</b>	Z0056212	<b>Catalog #</b>	N45164A	<b>Expiration Date:</b>	30 SEP 2020

Antibody Blend	
<b>Lot Number:</b>	D0013897
<b>Catalog Number:</b>	D21BY
<b>Expiration Date:</b>	31 AUG 2021
<b>Storage Temperature:</b>	2-8°C

Calibrator Blend	
<b>Lot Number:</b>	A0010447
<b>Catalog Number:</b>	C01BY
<b>Expiration Date:</b>	31 OCT 2020
<b>Storage Temperature:</b>	≤ -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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	26 JUL 2018