

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-Plex® and V-Plex Plus GLP-1 Total Kit

**Kit Catalog Numbers:** K1503P-Series

**Kit Lot Number:** K0081260

**Expiration Date:** 31 Jul 2021

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		GLP-1 (total) Control 1		GLP-1 (total) Control 2		GLP-1 (total) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	GLP-1 (total)	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	GLP-1 (total)	0.180	80% - 120%	< 20%	120	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	GLP-1 (total)	Calibrator	192	pM
		GLP-1 (total) Control 1	91.9	pM
		GLP-1 (total) Control 2	16.7	pM
		GLP-1 (total) Control 3	3	pM

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

GLP-1 (total) - Spot 1		
Conc. pM	ECL Signal	%CV
192	657,000	2.5
48.0	160,000	1.4
12.0	35,400	3.5
3.00	8,310	1.9
0.750	2,290	2.5
0.188	787	1.9
0.0469	413	3.1
0.00	297	4.5

**CERTIFICATE OF ANALYSIS**

**Components:**


Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) Plate	Z0056532	2-8°C	30 Nov 2021
GLP-1 (total) Antibody	D0081310	2-8°C	30 Nov 2022
GLP-1 (total) Calibrator	A0080223	2-8°C	31 Jul 2024
Diluent 12	M0120065	≤ -10°C	31 Jul 2021
Diluent 11	M0110040	≤ -10°C	30 Sep 2021
MSD GOLD™ Read Buffer	Not Kit Specific	Not Kit Specific	N/A

**Additional Components Included in V-PLEX Plus**

Plate Seals	Not Kit Specific
Wash Buffer (20x), 100 mL	Not Kit Specific
GLP-1 (total) Control 1	Not Kit Specific
GLP-1 (total) Control 2	Not Kit Specific
GLP-1 (total) Control 3	Not Kit Specific

**Statement:**

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		11 OCT 2019