

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus GLP-1 Active Kit  
**Kit Catalog Numbers:** K1503O-Series  
**Kit Lot Number:** K0080901  
**Expiration Date:** 31 Mar 2020

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		GLP-1 (active) Control 1		GLP-1 (active) Control 2		GLP-1 (active) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	GLP-1 (active)	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 (active)	0.300	80% - 120%	< 20%	120	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	GLP-1 (active)	Calibrator	276	pM
		GLP-1 (active) Control 1	122	pM
		GLP-1 (active) Control 2	44.3	pM
		GLP-1 (active) Control 3	14.6	pM

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

GLP-1 (active) - Spot 1		
Conc. pM	ECL Signal	%CV
276	1,130,000	2.2
69.0	228,000	3.0
17.3	45,300	3.5
4.31	10,100	2.3
1.08	2,660	2.5
0.270	831	2.4
0.0674	405	3.0
0.00	264	4.1

**CERTIFICATE OF ANALYSIS**

Page 2 of 2

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (active) SECTOR Plate	Z0056083	2-8°C	31 Mar 2020
GLP-1 (active) Antibody	D0081039	2-8°C	31 Mar 2021
GLP-1 (active) Calibrator Blend	A0080155	2-8°C	30 Nov 2022
Diluent 13	M0130042	≤ -10°C	31 Mar 2020
Diluent 12	M0120061	≤ -10°C	31 Mar 2020
Aprotinin	Not Kit Specific	Not Kit Specific	N/A
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), 100mL	Not Kit Specific
GLP-1 (active) Control 1	Not Kit Specific
GLP-1 (active) Control 2	Not Kit Specific
GLP-1 (active) 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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	03 JAN 2018