

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: K0081176

Expiration Date: 30 Jun 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	195	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
195	643,000	1.6
48.8	160,000	1.8
12.2	36,200	2.3
3.05	8,530	3.1
0.762	2,390	2.4
0.190	819	2.8
0.0476	421	3.0
0.00	303	3.9

CERTIFICATE OF ANALYSIS

Components:


Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) Plate	Z0056445	2-8°C	30 Jun 2021
GLP-1 (total) Antibody	D0081231	2-8°C	30 Jun 2022
GLP-1 (total) Calibrator	A0080194	2-8°C	31 Dec 2023
Diluent 12	M0120065	≤ -10°C	31 Jul 2021
Diluent 11	M0110038	≤ -10°C	31 Jul 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), 100mL	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
Review/Approval	Dawn Stollar	Quality Assurance		28 MAR 2019