

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: K0081199
Expiration Date: 29 Feb 2020

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	96.6	pM
		GLP-1 (total) Control 2	18.2	pM
		GLP-1 (total) Control 3	2.91	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	675,000	3.4
48.8	161,000	2.4
12.2	35,100	3.5
3.05	7,800	2.1
0.762	2,070	2.7
0.190	655	3.3
0.0476	310	2.3
0.00	206	2.5

CERTIFICATE OF ANALYSIS

Components:


Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) Plate	Z0056117	2-8°C	30 Jun 2020
GLP-1 (total) Antibody	D0081063	2-8°C	30 Jun 2020
GLP-1 (total) Calibrator	A0080194	2-8°C	31 Dec 2023
Diluent 12	M0120061	≤ -10°C	31 Mar 2020
Diluent 11	M0110034	≤ -10°C	29 Feb 2020
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		13 MAR 2019