

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® AND V-PLEX Plus GLP-1 Total Kit
Kit Lot Number: K0081624
Expiration Date: 31 Mar 2023
Kit Catalog Numbers: K1503P Series

Precision and Accuracy Specifications:

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

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

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	GLP-1 (total)	204	pM

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) Plate	Z0056835	2-8°C	31 May 2023
GLP-1 (total) Antibody	D0081606	2-8°C	30 Jun 2024
GLP-1 (total) Calibrator	A0080246	2-8°C	31 Dec 2024
Diluent 12	M0120069	≤ -10°C	31 Mar 2023
Diluent 11	M0110048	≤ -10°C	30 Jun 2023
MSD GOLD Read Buffer A	Not Kit Specific	Not Kit Specific	N/A

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 APR 2021