

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus GLP-1 Active Kit
Kit Lot Number: K0081835
Expiration Date: 31 Dec 2023
Kit Catalog Numbers: K1503O-Series

Precision and Accuracy Specifications:

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

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	GLP-1 (active)	267	pM

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (active) Plate	Z0057040	2-8°C	29 Feb 2024
GLP-1 (active) Antibody	D0081744	2-8°C	28 Feb 2025
GLP-1 (active) Calibrator	A0080195	2-8°C	31 Jan 2024
Diluent 13	M0130052	≤ -10°C	31 Dec 2023
Diluent 12	M0120072	≤ -10°C	31 Dec 2023

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		17 NOV 2021