



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus GLP-1 Total Kit
Kit Lot Number: K0082265
Expiration Date: 31 Aug 2025
Kit Catalog Numbers: K1503P-Series

Precision and Accuracy Specifications:

		Calibrator		GLP-1 (total) Control 1		GLP-1 (total) Control 2		GLP-1 (total) 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)	189	pM

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) SECTOR Plate	Z0057353	2-8°C	30 Sep 2025
GLP-1 (total) Antibody	D0082037	2-8°C	31 Oct 2026
GLP-1 (total) Calibrator	A0080330	2-8°C	31 May 2027
Diluent 11	M0110061	≤ -10°C	31 Aug 2025
Diluent 12	M0120080	≤ -10°C	31 Aug 2025

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		23 JUN 2023