



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus GLP-1 Total Kit
Lot Number: K0082540
Expiration Date: 31 AUG 2026
Catalog Number: K1503P-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) Antibody	D0082225	2-8°C	31 AUG 2027
GLP-1 (total) Calibrator	A0080394	2-8°C	30 APR 2029
GLP-1 (total) SECTOR Plate	Z0057451	2-8°C	31 AUG 2026
Diluent 11	M0110068	≤ -10°C	31 DEC 2026
Diluent 12	M0120086	≤ -10°C	31 DEC 2026

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

Precision, Accuracy, and Sensitivity 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)	0.18	80% - 120%	< 20%	120	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	GLP-1 (total)	223	pM

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 SEP 2024