



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus GLP-1 Active Kit
Lot Number: K0082472
Expiration Date: 31 JUL 2026
Catalog Number: K1503O-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (active) Antibody	D0082071	2-8°C	30 NOV 2026
GLP-1 (active) Calibrator	A0080398	2-8°C	31 MAY 2029
GLP-1 (active) Plate	Z0057466	2-8°C	31 OCT 2026
Diluent 12	M0120084	≤ -10°C	30 SEP 2026
Diluent 13	M0130059	≤ -10°C	31 JUL 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 (active) Control 1		GLP-1 (active) Control 2		GLP-1 (active) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	GLP-1 (Active)	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 (Active)	0.3	80% - 120%	< 20%	120	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	GLP-1 (Active)	247	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		13 AUG 2024