



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX GLP-1 Total Kit, QuickPlex**Lot Number:** K0082725**Expiration Date:** 30 SEP 2026**Catalog Number:** K1503P-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) 96-Well Sm Spot QuickPlex Ultra Plate	Z0057574	2-8°C	30 SEP 2027
GLP-1 (total) Antibody	D0082485	2-8°C	31 OCT 2028
GLP-1 (total) Calibrator	A0080404	2-8°C	31 AUG 2029
Diluent 11	M0110071	≤ -10°C	31 JUL 2027
Diluent 12	M0120084	≤ -10°C	30 SEP 2026

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

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

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	GLP-1 (total)	213	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		30 JUN 2025