



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus GLP-1 (total) Kit, SECTOR
Lot Number: K0082809
Expiration Date: 30 NOV 2027
Catalog Number: K1503P-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) 96-Well Small Spot SECTOR Plate	Z0057604	2-8°C	31 JAN 2028
GLP-1 (total) Antibody	D0082501	2-8°C	31 JAN 2029
GLP-1 (total) Calibrator	A0080404	2-8°C	31 AUG 2029
Diluent 11	M0110073	≤ -10°C	29 FEB 2028
Diluent 12	M0120091	≤ -10°C	30 NOV 2027

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		21 NOV 2025