



Extension Memo

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus GLP-1 Total Kit
Catalog Number(s): K1503P-Series
Lot Number: K0082604
New Expiration Date: 31 Mar 2026

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
GLP-1 (total) SECTOR Plate	Z0057274	31 Mar 2025	31 Mar 2026
Diluent 11	M0110059	30 Apr 2025	31 Mar 2026
Diluent 12	M0120078	30 Apr 2025	31 Mar 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		18 Dec 2024

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus GLP-1 Total Kit
Lot Number: K0082604
Expiration Date: 31 MAR 2025
Catalog Number: K1503P-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) Antibody	D0081938	2-8°C	31 MAR 2026
GLP-1 (total) Calibrator	A0080404	2-8°C	31 AUG 2029
GLP-1 (total) SECTOR Plate	Z0057274	2-8°C	31 MAR 2025
Diluent 11	M0110059	≤ -10°C	30 APR 2025
Diluent 12	M0120078	≤ -10°C	30 APR 2025

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)	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		18 DEC 2024