



EXTENSION MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus GLP-1 Active Kit

Catalog Number(s) / Size(s): K1503O-Series

Lot Number: K0082302

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082302 has been extended to 31 JAN 2025.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
GLP-1 (active) Calibrator	A0080195	31 Jan 2024	31 JAN 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		06 NOV 2023

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus GLP-1 Active Kit
Kit Lot Number: K0082302
Expiration Date: 31 Jan 2024
Kit Catalog Numbers: K1503O-Series

Precision and Accuracy Specifications:

		Calibrator		GLP-1 (active) Control 1		GLP-1 (active) Control 2		GLP-1 (active) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	GLP-1 (Active)	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 (Active)	0.300	80% - 120%	< 20%	120	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	GLP-1 (Active)	267	pM

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (active) SECTOR Plate	Z0057379	2-8°C	30 Nov 2025
GLP-1 (active) Antibody	D0081849	2-8°C	30 Sep 2025
GLP-1 (active) Calibrator	A0080195	2-8°C	31 Jan 2024
Diluent 12	M0120080	≤ -10°C	31 Aug 2025
Diluent 13	M0130057	≤ -10°C	31 Aug 2025

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		27 JUL 2023