



## 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:** K0082056

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082056 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) Plate	Z0057191	30 NOV 2024	31 JAN 2025
GLP-1 (active) Calibrator	A0080195	31 JAN 2024	31 JAN 2025
Diluent 13	M0130055	30 SEP 2024	31 JAN 2025
Diluent 12	M0120074	31 OCT 2024	31 JAN 2025

Function	Name	Signature	Date
Quality	Xiomara Talavera		05 APR 2024

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:** K0082056  
**Expiration Date:** 31 Jan 2024  
**Kit Catalog Numbers:** K1503O-Series

**Precision and Accuracy Specifications:**

		Calibrator		GLP-1 (active) Control 1		GLP-1 (active) Control 1		GLP-1 (active) Control 1	
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%

		LL			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) Plate	Z0057191	2-8°C	30 Nov 2024
GLP-1 (active) Antibody	D0081833	2-8°C	31 Aug 2025
GLP-1 (active) Calibrator	A0080195	2-8°C	31 Jan 2024
Diluent 13	M0130055	≤ -10°C	30 Sep 2024
Diluent 12	M0120074	≤ -10°C	31 Oct 2024

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		10 AUG 2022