



### CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus GLP-1 Active Kit  
**Lot Number:** K0082561  
**Expiration Date:** 31 JAN 2026  
**Catalog Number:** K1503O-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (active) Antibody	D0082355	2-8°C	30 APR 2028
GLP-1 (active) Calibrator	A0080398	2-8°C	31 MAY 2029
GLP-1 (active) SECTOR Plate	Z0057508	2-8°C	31 MAR 2027
Diluent 12	M0120082	≤ -10°C	31 MAR 2026
Diluent 13	M0130058	≤ -10°C	31 JAN 2026

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 (active) Control 1		GLP-1 (active) Control 2		GLP-1 (active) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	GLP-1 (Active)	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 (Active)	0.3	80% - 120%	< 20%	120	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	GLP-1 (Active)	247	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
<b>Review/Approval</b>	Karen Hamilla	Quality		09 DEC 2024