

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus GLP-1 Active Kit, SECTOR**Lot Number:** K0082923**Expiration Date:** 29 FEB 2028**Catalog Number:** K1503O-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (active) 96-Well Small Spot SECTOR Plate	Z0057676	2-8°C	30 SEP 2028
GLP-1 (active) Antibody	D0082355	2-8°C	30 APR 2028
GLP-1 (active) Calibrator	A0080417	2-8°C	31 JAN 2030
Diluent 12	M0120092	≤ -10°C	29 FEB 2028
Diluent 13	M0130064	≤ -10°C	29 FEB 2028

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)	228	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>	Michelle Likiam	Quality		11 MAY 2026