

### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX PLUS GLP-1 Active Kit  
**Kit Lot Number:** K0082258  
**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	Z0057149	2-8°C	31 Aug 2024
GLP-1 (active) Antibody	D0081833	2-8°C	31 Aug 2025
GLP-1 (active) Calibrator	A0080195	2-8°C	31 Jan 2024
Diluent 12	M0120074	≤ -10°C	31 Oct 2024
Diluent 13	M0130055	≤ -10°C	30 Sep 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		20 APR 2023