



### CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX PLUS Glucagon Kit  
**Lot Number:** K0082466  
**Expiration Date:** 31 AUG 2025  
**Catalog Number:** K150U5-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Glucagon Antibody	D0081964	2-8°C	30 JUN 2026
Glucagon Plate	Z0047978	2-8°C	30 SEP 2026
Metabolic Panel 1 Calibrator Blend	A0080343	2-8°C	30 SEP 2027
Diluent 11	M0110065	≤ -10°C	31 JUL 2026
Diluent 56	M0560005	≤ -10°C	31 AUG 2025

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

**Precision, Accuracy, and Sensitivity Specifications:**

Spot	Name	Calibrator		Metabolic Panel 1 Control 1		Metabolic Panel 1 Control 2		Metabolic Panel 1 Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Glucagon	80% - 120%	< 20%	95% - 135%	< 20%	95% - 135%	< 20%	95% - 135%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Glucagon	0.329	80% - 120%	< 20%	52	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	Glucagon	70	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		26 JUL 2024