



Extension Memo


Page 1 of 1

Product Description: V-PLEX® and V-PLEX PLUS Glucagon Kit
Catalog Number(s): K150U5-Series
Lot Number: K0082466
New Expiration Date: 31 Aug 2026

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 56	M0560005	31 Aug 2025	31 Aug 2026
Diluent 11	M0110065	31 Jul 2026	31 Aug 2026
Glucagon Antibody	D0081964	30 Jun 2026	31 Aug 2026

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 Jun 2025



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:

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


		LLOQ			ULOQ		
Spot	Name	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
Review/Approval	Karen Hamilla	Quality		26 JUL 2024