



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Glucagon Kit, SECTOR
Lot Number: K0082798
Expiration Date: 31 JAN 2028
Catalog Number: K150U5-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Glucagon Antibody	D0082294	2-8°C	31 JAN 2028
Glucagon 96-Well Small Spot SECTOR Plate	Z0048179	2-8°C	31 MAR 2028
Metabolic Panel 1 Calibrator Blend	A0080438	2-8°C	31 JUL 2030
Diluent 11	M0110073	≤ -10°C	29 FEB 2028
Diluent 56	M0560010	≤ -10°C	31 JAN 2028

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%	52.0	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Glucagon	79.4	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		10 NOV 2025