

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Glucagon Kit
Kit Lot Number: K0082147
Expiration Date: 31 Oct 2024
Kit Catalog Numbers: K150U5-Series

Precision and Accuracy 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%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 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	70	pM

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Glucagon Plate	Z0047775	2-8°C	31 Mar 2025
Glucagon Antibody	D0081851	2-8°C	31 Oct 2025
Metabolic Panel 1 Calibrator Blend	A0080343	2-8°C	30 Sep 2027
Diluent 11	M0110056	≤ -10°C	31 Oct 2024
Diluent 56	M0560004	≤ -10°C	31 Jan 2025

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		30 NOV 2022