

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human Insulin Kit  
**Kit Lot Number:** K0082257  
**Expiration Date:** 31 Oct 2024  
**Kit Catalog Numbers:** K151S5-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	Insulin	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	Insulin	0.0690	80% - 120%	< 20%	19.5	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	Insulin	28.9	uIU/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human Insulin Plate	Z0047838	2-8°C	31 Aug 2025
Human Insulin Antibody	D0081847	2-8°C	30 Sep 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		17 MAY 2023