



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human Insulin Kit**Lot Number:** K0082754**Expiration Date:** 31 DEC 2026**Catalog Number:** K151S5-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human Insulin 96-Well Small Spot SECTOR Plate	Z0048149	2-8°C	31 DEC 2027
Human Insulin Antibody	D0082290	2-8°C	31 JAN 2028
Metabolic Panel 1 Calibrator Blend	A0080438	2-8°C	31 JUL 2030
Diluent 11	M0110072	≤ -10°C	30 NOV 2027
Diluent 56	M0560007	≤ -10°C	31 DEC 2026

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	Insulin	80% - 120%	< 20%	95% - 135%	< 20%	95% - 135%	< 20%	95% - 135%	< 20%


Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Insulin	0.069	80% - 120%	< 20%	19.5	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Insulin	29.3	uIU/mL

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		23 SEP 2025