



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX Metabolic Panel 1 kit**Lot Number:** K0082733**Expiration Date:** 31 DEC 2026**Catalog Number:** K151X5-Series, K15325-Series, K150T5-Series, K15333-Series, K151Y5-Series, K15334-Series, K15332-Series, K15335-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
GIP (active) Antibody	D0082293	2-8°C	31 JAN 2028
GLP-1 (active) Antibody	D0082292	2-8°C	31 DEC 2027
Glucagon Antibody	D0082294	2-8°C	31 JAN 2028
Human C-peptide Antibody	D0082291	2-8°C	31 DEC 2027
Human Insulin Antibody	D0082290	2-8°C	31 JAN 2028
Human Leptin Antibody	D0082296	2-8°C	29 FEB 2028
Human PP Antibody	D0082295	2-8°C	31 JAN 2028
Metabolic Panel 1 Calibrator Blend	A0080402	2-8°C	31 JUL 2029
Metabolic Panel 1, 96-Well 10-Spot SECTOR Plate	Z0047999	2-8°C	31 DEC 2026
Diluent 11	M0110071	≤ -10°C	31 JUL 2027
Diluent 56	M0560009	≤ -10°C	30 JUN 2027

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%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	C-Peptide	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	GLP-1 (Active)	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	GIP (active)	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	Glucagon	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	PP	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	Leptin	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Insulin	0.069	80% - 120%	< 20%	19.5	80% - 120%	< 20%
2	C-Peptide	19.0	80% - 120%	< 20%	3150	80% - 120%	< 20%
3	GLP-1 (Active)	0.096	80% - 120%	< 20%	39	80% - 120%	< 20%
7	GIP (active)	3.18	80% - 120%	< 20%	390	80% - 120%	< 20%
8	Glucagon	0.329	80% - 120%	< 20%	52	80% - 120%	< 20%
9	PP	0.41	80% - 120%	< 20%	325	80% - 120%	< 20%
10	Leptin	37.7	80% - 120%	< 20%	7150	80% - 120%	< 20%



CERTIFICATE OF ANALYSIS

Non-Specific Binding Specifications:

Spot	Name	Spec
1	Insulin	< 1.0%
2	C-peptide	< 1.0%
3	GLP-1 (active)	< 1.0%
7	GIP (active)	< 1.0%
8	Glucagon	< 1.0%
9	PP	< 1.0%
10	Leptin	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Insulin	28.9	uIU/mL
2	C-Peptide	10800	pg/mL
3	GLP-1 (Active)	65.4	pM
7	GIP (active)	570	pg/mL
8	Glucagon	81	pM
9	PP	445	pg/mL
10	Leptin	11400	pg/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		25 JUN 2025