



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Metabolic Panel 1 Kit, QuickPlex**Lot Number:** K0082703**Expiration Date:** 31 MAY 2026**Catalog Number:** K15325-Series; K151S5-Series; K151X5-Series; K150T5-Series; K150U5-Series; K151Y5-Series; K151V5-Series; K15332-Series; K15333-Series; K15334-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	D0082355	2-8°C	30 APR 2028
Glucagon Antibody	D0082073	2-8°C	31 OCT 2026
Human C-peptide Antibody	D0082070	2-8°C	30 NOV 2026
Human Insulin Antibody	D0082069	2-8°C	31 OCT 2026
Human Leptin Antibody	D0082075	2-8°C	31 DEC 2026
Human PP Antibody	D0082074	2-8°C	30 NOV 2026
Metabolic Panel 1 Calibrator Blend	A0080402	2-8°C	31 JUL 2029
Metabolic Panel 1, 96-Well 10-Spot QuickPlex Ultra Plate	Z0048117	2-8°C	31 OCT 2027
Diluent 11	M0110069	≤ -10°C	28 FEB 2027
Diluent 56	M0560006	≤ -10°C	31 MAY 2026

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	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%	70% - 130%	< 20%	70% - 130%	< 20%	70% - 130%	< 20%
7	GIP (active)	80% - 120%	< 20%	70% - 130%	< 20%	70% - 130%	< 20%	70% - 130%	< 20%
8	Glucagon	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	PP	80% - 120%	< 20%	70% - 130%*	< 20%	70% - 130%*	< 20%	70% - 130%*	< 20%
10	Leptin	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

*See Additional Comments

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	Insulin	0.069	80% - 120%	< 20%	19.5	80% - 120%	< 20%
2	C-Peptide	19	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	0.405	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	Insulin	C-peptide	GLP-1 (active)	GIP (active)	Glucagon	PP	Leptin
1	Insulin	NA	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%
2	C-peptide	< 1.0%	NA	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%
3	GLP-1 (active)	< 1.0%	< 1.0%	NA	< 1.0%	< 1.0%	< 1.0%	< 1.0%
7	GIP (active)	< 1.0%	< 1.0%	< 1.0%	NA	< 1.0%	< 1.0%	< 1.0%
8	Glucagon	< 1.0%	< 1.0%	< 1.5%	< 1.0%	NA	< 1.0%	< 1.0%
9	PP	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	NA	< 1.0%
10	Leptin	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	NA

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

Additional Comments:

Spot	Name	Specification	Comment
9	PP	Control Conc.	There is a 15% bias for the PP assay relative to SECTOR.

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		02 SEP 2025