

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX and V-PLEX Plus Metabolic Panel 1 (human) Kit, SECTOR
Lot Number: K0082813
Expiration Date: 31 JAN 2028
Catalog Number: K15325-Series, K151X5-Series, K150T5-Series, K151Y5-Series, K15332-Series, K15333-Series, K15334-Series, K15335-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
GIP (active) Antibody	D0082542	2-8°C	30 JUN 2029
GLP-1 (active) Antibody	D0082541	2-8°C	31 MAR 2029
Glucagon Antibody	D0082543	2-8°C	31 MAY 2029
Human C-peptide Antibody	D0082540	2-8°C	31 MAY 2029
Human Insulin Antibody	D0082539	2-8°C	30 JUN 2029
Human Leptin Antibody	D0082545	2-8°C	31 MAY 2029
Human PP Antibody	D0082544	2-8°C	31 MAY 2029
Metabolic Panel 1 Calibrator Blend	A0080438	2-8°C	31 JUL 2030
Metabolic Panel 1, 96-Well 10-Spot SECTOR Plate	Z0048187	2-8°C	31 MAY 2028
Diluent 11	M0110073	≤ -10°C	29 FEB 2028
Diluent 56	M0560010	≤ -10°C	31 JAN 2028

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	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	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	29.3	uIU/mL
2	C-Peptide	11100	pg/mL
3	GLP-1 (Active)	68	pM
7	GIP (active)	563	pg/mL
8	Glucagon	79.4	pM
9	PP	430	pg/mL
10	Leptin	9000	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		12 FEB 2026