

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Metabolic Panel 1 Control Pack
Catalog Number: C4325-1
Lot Number: CP000182
Expiration Date: 31 AUG 2030

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Metabolic Panel 1 Control 1	A00C1026	2-8°C	31 AUG 2030
Metabolic Panel 1 Control 2	A00C1027	2-8°C	31 AUG 2030
Metabolic Panel 1 Control 3	A00C1028	2-8°C	31 AUG 2030

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Insulin	11.0	1.40	0.244	μIU/mL
C-peptide	4130	382	86.8	pg/mL
GLP-1 (active)	21.4	2.66	0.336	pM
GIP (active)	240	23.9	11.1	pg/mL
Glucagon	27.6	4.23	1.07	pM
PP	188	22.3	0.863	pg/mL
Leptin	3480	497	119	pg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		14OCT 2025