



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Metabolic Panel 1 Control Pack

Catalog Number: C4325-1

Lot Number: CP000130

Expiration Date: 31 JUL 2026

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Insulin	13.7	1.86	0.293	μIU/mL
C-Peptide	5560	487	90.3	pg/mL
GLP-1 (active)	23.9	2.98	0.41	pM
GIP (active)	242	27	12.8	pg/mL
Glucagon	31.2	4.1	1.23	pM
PP	170	25.9	1.41	pg/mL
Leptin	3210	448	127	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Metabolic Panel 1 Control 1	A00C0801	2-8°C	31 JUL 2026
Metabolic Panel 1 Control 2	A00C0802	2-8°C	31 JUL 2026
Metabolic Panel 1 Control 3	A00C0803	2-8°C	31 JUL 2026

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		29 APR 2022