



## CERTIFICATE OF ANALYSIS

**Control Pack Description:** Metabolic Panel 1 Control Pack**Catalog Number:** C4325-1**Lot Number:** CP000144**Expiration Date:** 29 FEB 2028**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Insulin	10.7	1.36	0.244	μIU/mL
C-Peptide	4770	450	83.6	pg/mL
GLP-1 (active)	21.4	2.60	0.336	pM
GIP (active)	221	24.4	10.5	pg/mL
Glucagon	26.4	3.40	1.05	pM
PP	154	21.4	1.21	pg/mL
Leptin	3390	438	120	pg/mL

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Metabolic Panel 1 Control 1	A00C0883	2-8°C	29 FEB 2028
Metabolic Panel 1 Control 2	A00C0884	2-8°C	29 FEB 2028
Metabolic Panel 1 Control 3	A00C0885	2-8°C	29 FEB 2028

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		04 MAY 2023