



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Control Pack Description:** Metabolic Panel 1 Control Pack

**Catalog Number:** C4325-1

**Lot Number:** CP000096

**Expiration Date:** 31 OCT 2024

### Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Insulin	12.0	1.65	0.293	uIU/mL
C-Peptide	4990	581	82.3	pg/mL
GLP-1 (active)	22.7	2.78	0.336	pM
GIP (active)	227	22.5	9.61	pg/mL
Glucagon	28.2	3.76	1.11	pM
PP	157	22.2	1.42	pg/mL
Leptin	3290	458	104	pg/mL

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Metabolic Panel 1 Control 1	A00C0673	2-8°C	31 OCT 2024
Metabolic Panel 1 Control 2	A00C0674	2-8°C	31 OCT 2024
Metabolic Panel 1 Control 3	A00C0697	2-8°C	30 NOV 2024

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		06 DEC 2019