



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Metabolic Panel 1 Calibrator Blend

Catalog Number: C0325-2

Lot Number: A0080343

Expiration Date: 30 SEP 2027

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Insulin	28.9	μIU/mL
C-Peptide	10600	pg/mL
GLP-1 (active)	67.2	pM
GIP (active)	578	pg/mL
Glucagon	70	pM
PP	435	pg/mL
Leptin	9720	pg/mL

Reconstitution Volume: 1000 μL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 NOV 2022