



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Metabolic Panel 1 Calibrator Blend

Catalog Number: C0325-2

Lot Number: A0080230

Expiration Date: 30 SEP 2024

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Insulin	30.6	μIU/mL
C-Peptide	11300	pg/mL
GLP-1 (active)	67.7	pM
GIP (active)	551	pg/mL
Glucagon	77.4	pM
PP	443	pg/mL
Leptin	9930	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		22 OCT 2019