



CERTIFICATE OF ANALYSIS

Product Description: Metabolic Panel 1 Calibrator Blend
Catalog Number: C0325-2
Lot Number: A0080438
Expiration Date: 31 JUL 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Insulin	29.3	μIU/mL
C-peptide	11100	pg/mL
GLP-1 (active)	68	pM
GIP (active)	563	pg/mL
Glucagon	79.4	pM
PP	430	pg/mL
Leptin	9000	pg/mL

Reconstitution Volume: 1000 μL

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 SEP 2025