



CERTIFICATE OF ANALYSIS

Product Description: Metabolic Panel 1 Calibrator Blend
Catalog Number: C0325-2
Lot Number: A0080467
Expiration Date: 30 JUN 2031
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Insulin	30	uIU/mL
C-peptide	9470	pg/mL
GLP-1 (active)	67.3	pM
GIP (active)	575	pg/mL
Glucagon	80.5	pM
PP	425	pg/mL
Leptin	8580	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		13 AUG 2026