



CERTIFICATE OF ANALYSIS

Control Pack Description: GLP-1 (total) Control Pack
Catalog Number: C403P-1
Lot Number: CP000169
Expiration Date: 31 OCT 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) Control 1	A00C0969	2-8°C	31 OCT 2029
GLP-1 (total) Control 2	A00C0970	2-8°C	31 OCT 2029
GLP-1 (total) Control 3	A00C0971	2-8°C	31 OCT 2029

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
GLP-1 (total)	83.7	17.5	2.57	pM

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 JAN 2025