



### CERTIFICATE OF ANALYSIS

**Control Pack Description:** GLP-1 (active) Control Pack  
**Catalog Number:** C4266-1  
**Lot Number:** CP000160  
**Expiration Date:** 30 APR 2029

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (active) Control 1	A00C0943	2-8°C	30 APR 2029
GLP-1 (active) Control 2	A00C0944	2-8°C	30 APR 2029
GLP-1 (active) Control 3	A00C0945	2-8°C	30 APR 2029

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
GLP-1 (active)	128	26.2	2.96	pM

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		13 JUN 2024