



## Extension Memo

Page 1 of 1

**Product Description:** GLP-1 (total) Control Pack  
**Catalog Number(s):** C403P-1  
**Lot Number:** CP000085  
**New Expiration Date:** 30 Sep 2025

Based on additional testing, the original date of expiration for this control pack has been extended by 20 months.

The expiration dates for the following controls have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
GLP-1 (total) Control 1	A00C0634	31 Dec 2023	30 Sep 2025
GLP-1 (total) Control 2	A00C0635	31 Dec 2023	30 Sep 2025
GLP-1 (total) Control 3	A00C0636	31 Dec 2023	30 Sep 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		07 Oct 2024



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** GLP-1 (total) Control Pack

**Catalog Number:** C403P-1

**Lot Number:** CP000085

**Expiration Date:** 31 DEC 2023

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

Units = pM	Assay
	<b>GLP (total)</b>
Control 1 (A00C0634)	91.9
Control 2 (A00C0635)	16.7
Control 3 (A00C0636)	3.00

**Reconstitution Volume:** 250 µL

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) Control 1	A00C0634	2-8°C	31 DEC 2023
GLP-1 (total) Control 2	A00C0635	2-8°C	31 DEC 2023
GLP-1 (total) Control 3	A00C0636	2-8°C	31 DEC 2023

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		07 MAR 2019