



EXTENSION MEMO

Page 1 of 1

Product Description(s): GLP-1 (active) Control Pack


Catalog Number(s) / Size(s): C4266-1

Lot Number: CP000063

Based on additional testing, the original date of expiration for this control pack has been extended by 24 months. The expiration date of CP000063 has been extended to 30 Nov 2024.

The expiration dates for the following controls have been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
GLP-1 (active) Control 1	A00C0535	30 NOV 2022	30 NOV 2024
GLP-1 (active) Control 2	A00C0536	30 NOV 2022	30 NOV 2024
GLP-1 (active) Control 3	A00C0545	30 NOV 2022	30 NOV 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		02 NOV 2023

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: GLP-1 (active) Control Pack

Catalog Number: C4266-1

Lot Number: CP000063

Expiration Date: 30 NOV 2022

Storage Temperature: 2-8°C

Concentration of Materials Tested:

Units = pM	Assay
	GLP-1 (active)
Control 1 (A00C0535)	137
Control 2 (A00C0536)	34.2
Control 3 (A00C0545)	3.56

Reconstitution Volume: 250 µL

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (active) Control 1	A00C0535	2-8°C	30 NOV 2022
GLP-1 (active) Control 2	A00C0536	2-8°C	30 NOV 2022
GLP-1 (active) Control 3	A00C0545	2-8°C	30 NOV 2022

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	11 DEC 2017