



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: GLP-1 (total) Control Pack

Catalog Number: C403P-1

Lot Number: CP000067

Expiration Date: 31 JAN 2023

Storage Temperature: 2-8°C

Concentration of Materials Tested:

Units = pM	Assay
	GLP-1 (total)
Control 1 (A00C0562)	96.6
Control 2 (A00C0563)	18.2
Control 3 (A00C0564)	2.91

Reconstitution Volume: 250 µL

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
GLP-1 (total) Control 1	A00C0562	2-8°C	31 JAN 2023
GLP-1 (total) Control 2	A00C0563	2-8°C	31 JAN 2023
GLP-1 (total) Control 3	A00C0564	2-8°C	31 JAN 2023

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>	23 MAR 2018