



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** GLP-1 (total) Calibrator  
**Catalog Number:** C003P-2  
**Lot Number:** A0080316  
**Expiration Date:** 31 DEC 2026  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GLP-1 (total)	195	pM

**Reconstitution Volume:** 1000 µL

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	<i>Karen Hamilla</i>	16 MAR 2022