



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: GLP-1 (total) Calibrator

Catalog Number: C003P-2

Lot Number: A0080330

Expiration Date: 31 MAY 2027


Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GLP-1 (total)	189	pM

Reconstitution Volume: 1000 µL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 JUL 2022