



CERTIFICATE OF ANALYSIS

Product Description: GLP-1 (total) Calibrator
Catalog Number: C003P-2
Lot Number: A0080404
Expiration Date: 31 AUG 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GLP-1 (total)	213	pM

Reconstitution Volume: 1000 µL

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Hamilla, Karen	Quality		30 SEP 2024