



CERTIFICATE OF ANALYSIS

Product Description: GLP-1 (active) Calibrator
Catalog Number: C0266-1
Lot Number: A0080398
Expiration Date: 31 MAY 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GLP-1 (active)	247	pM

Reconstitution Volume: 1000 μ L

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	Karen Hamilla	Quality		22 JUL 2024