



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: GLP-1 (active) Calibrator

Catalog Number: C016L-2

Lot Number: A00U0184

Expiration Date: 30 NOV 2025

Storage Temperature: 2-8°C

Intended Assay Species: Human

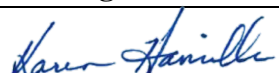
Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GLP-1 (active)	0.165	pmol	658	pM

Reconstitution Volume: 250 µL

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 MAR 2021