



## CERTIFICATE OF ANALYSIS

**Product Description:** GLP-1 (active) Calibrator  
**Catalog Number:** C016L-2  
**Lot Number:** A00U0250  
**Expiration Date:** 31 JAN 2027  
**Storage Temperature:** 2-8°C  
**Intended Assay Species:** Human, Mouse, Rat


**Concentration Assignment:**

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GLP-1 (active)	0.157	pmol	629	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
<b>Review/Approval</b>	Karen Hamilla	Quality		23 APR 2024