



CERTIFICATE OF ANALYSIS

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, Mouse, Rat

Concentration Assignment:

| Analyte | Amount of Analyte in Calibrator | Unit of Measure | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|----------------|---------------------------------|-----------------|---|-----------------|
| GLP-1 (active) | 0.165 | 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 | | 23 APR 2024 |