

**CERTIFICATE OF ANALYSIS**

**Control Pack Description:** GLP-1 (total) Control Pack  
**Catalog Number:** C403P-1  
**Lot Number:** CP000195  
**Expiration Date:** 31 MAY 2031

**Components:**

| Description             | Lot Number | Storage Temperature | Expiration Date |
|-------------------------|------------|---------------------|-----------------|
| GLP-1 (total) Control 1 | A00C1063   | 2-8°C               | 31 MAY 2031     |
| GLP-1 (total) Control 2 | A00C1064   | 2-8°C               | 31 MAY 2031     |
| GLP-1 (total) Control 3 | A00C1065   | 2-8°C               | 31 MAY 2031     |

**Concentration Assignment:**

| Analyte       | Control 1 | Control 2 | Control 3 | Unit of Measure |
|---------------|-----------|-----------|-----------|-----------------|
| GLP-1 (total) | 86.3      | 16.3      | 2.49      | pM              |

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  |  | 01 SEP 2026 |