

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus GLP-1 Active Kit
Lot Number: K0082655
Expiration Date: 31 MAR 2027
Catalog Number: K1503O-Series

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| GLP-1 (active) 96-Well Small Spot SECTOR Plate | Z0048087 | 2-8°C | 31 JUL 2027 |
| GLP-1 (active) Antibody | D0082355 | 2-8°C | 30 APR 2028 |
| GLP-1 (active) Calibrator | A0080417 | 2-8°C | 31 JAN 2030 |
| Diluent 12 | M0120088 | ≤ -10°C | 30 APR 2027 |
| Diluent 13 | M0130061 | ≤ -10°C | 31 MAR 2027 |

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

| Spot | Name | Calibrator | | GLP-1 (active) Control 1 | | GLP-1 (active) Control 2 | | GLP-1 (active) Control 3 | |
|------|----------------|------------|-------|--------------------------|-------|--------------------------|-------|--------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | GLP-1 (Active) | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|----------------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | GLP-1 (Active) | 0.30 | 80% - 120% | < 20% | 120 | 80% - 120% | < 20% |

Stock Concentration of Calibrator:

| Spot | Name | Conc. | Units |
|------|----------------|-------|-------|
| 1 | GLP-1 (Active) | 228 | pM |

Dilution to top of Curve: 1X

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

| | Name | Function | Signature | Date |
|-----------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality | | 10 APR 2025 |