



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human Leptin Kit**Lot Number:** K0082749**Expiration Date:** 30 JUN 2027**Catalog Number:** K151V5-Series**Lot-Specific Components:**

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Human Leptin 96-Well Small Spot SECTOR Plate | Z0048146 | 2-8°C | 31 DEC 2027 |
| Human Leptin Antibody | D0082499 | 2-8°C | 31 DEC 2028 |
| Metabolic Panel 1 Calibrator Blend | A0080402 | 2-8°C | 31 JUL 2029 |
| Diluent 11 | M0110072 | ≤ -10°C | 30 NOV 2027 |
| Diluent 56 | M0560009 | ≤ -10°C | 30 JUN 2027 |

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

Precision, Accuracy, and Sensitivity Specifications:

| Spot | Name | Calibrator | | Metabolic Panel 1 Control 1 | | Metabolic Panel 1 Control 2 | | Metabolic Panel 1 Control 3 | |
|------|--------|------------|-------|-----------------------------|-------|-----------------------------|-------|-----------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | Leptin | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |


| Spot | Name | LLOQ | | | ULOQ | | |
|------|--------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | Leptin | 37.7 | 80% - 120% | < 20% | 7150 | 80% - 120% | < 20% |

Stock Concentration of Calibrator:

| Spot | Name | Conc. | Units |
|------|--------|-------|-------|
| 1 | Leptin | 11400 | pg/mL |

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 |  | 05 AUG 2025 |