



### CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX PLUS Human Leptin Kit  
**Lot Number:** K0082518  
**Expiration Date:** 31 MAY 2026  
**Catalog Number:** K151V5-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human Leptin Antibody	D0082254	2-8°C	31 OCT 2027
Human Leptin Plate	Z0047983	2-8°C	31 OCT 2026
Metabolic Panel 1 Calibrator Blend	A0080343	2-8°C	30 SEP 2027
Diluent 11	M0110067	≤ -10°C	30 NOV 2026
Diluent 56	M0560006	≤ -10°C	31 MAY 2026

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	9720	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
<b>Review/Approval</b>	Hamilla, Karen	Quality		19 SEP 2024