

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human Leptin Kit

**Kit Lot Number:** K0082194

**Expiration Date:** 31 Jan 2025

**Kit Catalog Numbers:** K151V5-Series

**Precision and Accuracy Specifications:**

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

		LLOQ			ULOQ		
Spot	Name	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

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human Leptin Plate	Z0047798	2-8°C	31 May 2025
Human Leptin Antibody	D0081853	2-8°C	30 Nov 2025
Metabolic Panel 1 Calibrator Blend	A0080343	2-8°C	30 Sep 2027
Diluent 11	M0110059	≤ -10°C	30 Apr 2025
Diluent 56	M0560004	≤ -10°C	31 Jan 2025

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		18 JAN 2023