



Meso Scale Diagnostics, LLC.

### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human Leptin Kit  
**Kit Lot Number:** K0082319  
**Expiration Date:** 31 Aug 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 SECTOR Plate	Z0047882	2-8°C	30 Nov 2025
Human Leptin Antibody	D0082093	2-8°C	31 Dec 2026
Metabolic Panel 1 Calibrator Blend	A0080343	2-8°C	30 Sep 2027
Diluent 11	M0110061	≤ -10°C	31 Aug 2025
Diluent 56	M0560005	≤ -10°C	31 Aug 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		02 AUG 2023