

User Guide

# Human Leptin ELISA Kit

96-Well Plate

## EZHL-80SK

<b>Intended Use</b> .....	<b>2</b>	<b>Calculations</b> .....	<b>16</b>
<b>Principles of Assay</b> .....	<b>2</b>	<b>Interpretation</b> .....	<b>16</b>
<b>Reagents Supplied</b> .....	<b>3</b>	<b>Normal Range</b> .....	<b>16</b>
<b>Storage and Stability</b> .....	<b>4</b>	<b>Standard Curve</b> .....	<b>17</b>
<b>Reagent Precautions</b> .....	<b>4</b>	<b>Assay Characteristics</b> .....	<b>18</b>
Sodium Azide.....	4	Sensitivity .....	18
Hydrochloric Acid .....	4	Specificity .....	18
Symbol Definitions.....	4	Precision .....	19
<b>Materials Required</b> .....	<b>5</b>	Spike Recovery of Human Leptin in Assay Samples.....	20
<b>Sample Collection and Storage</b> ...	<b>5</b>	Spike Recovery of Human Leptin in Sensitive Assay Samples.....	21
<b>Human Leptin ELISA Assay Procedure I</b> .....	<b>6</b>	Linearity of Sample Dilution .....	22
<b>Assay Procedure I for Human Leptin ELISA Kit</b> .....	<b>8</b>	Linearity of Sample Dilution (Sensitive Assay) .....	23
<b>Microtiter Plate Arrangement</b> .....	<b>9</b>	<b>Correlation Graph</b> .....	<b>24</b>
<b>Human Leptin ELISA (Sensitive) Assay Procedure II</b> .....	<b>10</b>	Human Leptin Correlation RIA vs. ELISA .....	24
Human Leptin (Sensitive)		<b>Quality Controls</b> .....	<b>25</b>
Standard Preparation .....	10	<b>Troubleshooting</b> .....	<b>25</b>
Human Leptin (Sensitive) Quality Control 1 and 2 Preparation .....	11	<b>Product Ordering</b> .....	<b>26</b>
Human Leptin (Sensitive) Assay .	12	Replacement Reagents.....	26
<b>Assay Procedure II for Human Leptin (Sensitive) ELISA kit</b> ....	<b>14</b>	<b>Notice</b> .....	<b>27</b>
<b>Microtiter Plate Arrangement II</b> .....	<b>15</b>	Technical Assistance .....	27
		Terms and Conditions of Sale.....	27
		Contact Information.....	27

For research use only. Not for use in diagnostic procedures.

---

## Intended Use

This Human Leptin ELISA kit is used for the non-radioactive quantification of human leptin in serum, plasma and other biological media. One kit is sufficient to measure 37 unknown samples in duplicate.

**This kit is for research use only. Not for use in diagnostic procedures.**

## Principles of Assay

This assay is a Sandwich ELISA based, sequentially, on:

- Capture of human leptin by a polyclonal rabbit anti-human leptin antibody immobilized on a 96-well microtiter plate
- Washing of unbound materials from samples
- Binding of a biotinylated monoclonal antibody to the captured human leptin
- Washing of unbound materials from samples
- Binding of streptavidin horseradish peroxidase to the immobilized biotinylated antibodies
- Washing of excess free enzyme conjugates
- Quantification of bound streptavidin-horseradish peroxidase activities in the presence of the substrate 3,3',5,5'-tetramethylbenzidine

The enzyme activity is measured spectrophotometrically by the increased absorbance at 450 nm–590 nm after acidification of formed products. Since the increase in absorbance is directly proportional to the amount of captured human leptin in the unknown sample, the latter can be derived by interpolation from a reference curve generated in the same assay with reference standards of known concentrations of human leptin.

## Reagents Supplied

Each kit is sufficient to run one 96-well plate and contains the following reagents:

**Note:** Store all reagents at 2-8 °C

Reagents Supplied	Volume	Quantity	Cat. No.
Human Leptin ELISA Plate with 2 plate sealers	-	1 plate 2 plates	EP81
10X HRP Wash Buffer Concentrate	50 mL	2 bottles	EWB-HRP
Human Leptin ELISA Standards Human Leptin in Assay Buffer: 0.78, 1.56, 3.125, 6.25, 12.5, 25, 50, and 100 ng/mL <b>Note:</b> The standard(s) in this kit have been calibrated to an International Reference standard, NIBSC code # 97/594.	1 mL	8 vials	E8081-K
Quality Controls 1 and 2	1 mL	1 vial	E6081-K
Assay Buffer	10 mL	1 vial	EABTR
Human Leptin Detection Antibody	11 mL	1 vial	E1081
Enzyme Solution	12 mL	1 vial	EHRP-3
Stop Solution 0.3 M HCl ( <b>Caution:</b> Corrosive material)	12 mL	1 vial	ET-TMB
Substrate Solution 3,3',5,5'-tetramethylbenzidine in buffer (light-sensitive, avoid unnecessary exposure to light)	12 mL	1 vial	ESS-TMB

For research use only. Not for use in diagnostic procedures.

## Storage and Stability

Recommended storage for kit components is 2-8 °C.

All components are shipped and stored at 2-8 °C. Once opened, liquid standards and controls can be stored up to 30 days at 2-8 °C. Refer to expiration dates on all reagents prior to use. Do not mix reagents from different kits unless they have the same lot numbers.

## Reagent Precautions

### Sodium Azide

Sodium azide or ProClin™ has been added to some reagents as a preservative. Although the concentrations are low, Sodium azide and ProClin™ may react with lead and copper plumbing to form highly explosive metal azides. Dispose of unused contents and waste in accordance with international, federal, state, and local regulations.

### Hydrochloric Acid

Hydrochloric acid is corrosive and can cause eye and skin burns. It is harmful if swallowed and can cause respiratory and digestive tract burns. Avoid contact with skin and eyes. Do not swallow or ingest.

## Symbol Definitions

Ingredient	Cat. No.	Full Label
Stop Solution	ET-TMB	 <b>Warning.</b> May be corrosive to metals.
10X HRP Wash Buffer Concentrate	EWB-HRP	 <b>Warning.</b> May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.
Human Leptin Detection Antibody	E1081	 <b>Warning.</b> Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.

## Materials Required (Not Provided)

- Multi-channel Pipettes and pipette tips: 50  $\mu$ L-300  $\mu$ L
- Pipettes with tips: 10  $\mu$ L-200  $\mu$ L
- Buffer and Reagent Reservoirs
- Vortex Mixer
- De-ionized water
- Microtiter Plate Reader capable of reading absorbency at 450 nm
- Orbital Microtiter Plate Shaker
- Absorbent Paper or Cloth

## Sample Collection and Storage

1. To prepare serum, whole blood is directly drawn into a Vacutainer<sup>®</sup> serum tube that contains no anti-coagulant. Let blood clot at room temperature for 30 minutes.
2. Promptly centrifuge the clotted blood at 2000 to 3000 x *g* for 15 minutes at  $4 \pm 2$  °C.
3. Transfer and store serum samples in separate tubes. Date and identify each sample.
4. Use freshly prepared serum or store samples at  $< 0$  or  $-20$  C for later use. Avoid multiple ( $> 5$ ) freeze/thaw cycles.
5. To prepare plasma samples, whole blood should be collected into Vacutainer<sup>®</sup> EDTA plasma tubes and centrifuged immediately after collection. Observe same precautions in the preparation of serum samples.
6. If heparin is to be used as an anti-coagulant, the effect on the assay outcome at the dose of heparin used should be pre-determined.
7. Avoid using samples with gross hemolysis or lipemia.

## Human Leptin ELISA Assay Procedure I

Sensitivity: 0.78 ng/mL–100 ng/mL

Pre-warm all reagents to room temperature immediately before setting up the assay.

1. Dilute the concentrated Wash Buffer 10-fold by adding the entire contents of both bottles of buffer to 900 mL de-ionized or glass distilled water.
2. Remove required number of strips from the Microtiter Assay Plate. Unused strips should be resealed in the foil pouch with the desiccant provided and stored at 2–8 °C. Assemble strips in an empty plate holder and add 300  $\mu$ L of diluted Wash Buffer to each well. Incubate at room temperature for 5 minutes. Decant wash buffer and remove the residual amount from all wells by inverting the plate and tapping it smartly onto absorbent towels several times. Do not let wells dry before proceeding to the next step. If an automated machine is used for the assay, follow the manufacturer's instructions for all washing steps described in this protocol.
3. Add 75  $\mu$ L Assay Buffer into all wells.
4. Add in duplicate 25  $\mu$ L Assay Buffer to blank wells. Refer to [Microtiter Plate Arrangement I](#) for suggested well Orientations.
5. Add in duplicate 25  $\mu$ L Human Leptin Standards in order of ascending concentration to the appropriate wells. Add in duplicate 25  $\mu$ L QC1 and 25  $\mu$ L QC2 to the appropriate wells. Add sequentially 25  $\mu$ L of samples in duplicate to the remaining wells. **For best results all additions should be completed within one hour.**
6. Cover the plate with plate sealer and incubate at room temperature for 2 hours on an orbital microtiter plate shaker set to rotate at moderate speed, about 400 to 500 rpm.
7. Remove plate sealer and decant solutions from the plate. Tap as before to remove residual solutions in the wells.
8. Wash wells 3 times with diluted Wash Buffer, 300  $\mu$ L per well per wash. Decant and tap after each wash to remove residual buffer.
9. Add 100  $\mu$ L Detection Antibody to each well. Cover the plate with sealer and incubate at room temperature for 30 minutes on the microtiter plate shaker.
10. Remove sealer and decant solution from the plate. Tap as before to remove residual solutions in the wells.
11. Add 100  $\mu$ L Enzyme Solution to each well. Cover plate with sealer and incubate with moderate shaking at room temperature for 30 minutes on the microtiter plate shaker.
12. Remove sealer, decant solution from the plate, and tap plate to remove the residual fluid.
13. Wash wells 5 times with diluted Wash Buffer, 300  $\mu$ L per well per wash. Decant and tap firmly after each wash to remove residual buffer.

For research use only. Not for use in diagnostic procedures.

- 
14. Add 100  $\mu$ L of Substrate Solution to each well, cover plate with sealer and shake on the plate shaker for 5-20 minutes. Blue color should be formed in wells of Leptin standards with intensity proportional to increasing concentrations of Leptin.

**Note:** Please be aware that the color may develop more quickly or more slowly than the recommended incubation time depending on the localized room temperature. Please visually monitor the color development to optimize the incubation time. One can monitor color development using 370 nm filter, if available on the spectrophotometer. When the absorbance is between 1.2 and 1.8 at 370 nm, the stop solution can be added to terminate the color development.

15. Remove sealer and add 100  $\mu$ L of Stop Solution (**Caution:** Corrosive solution) and shake plate by hand to ensure complete mixing of solution in all wells. The blue color should turn to yellow after acidification. Wipe the bottom of the microtiter plate to remove any residue prior to reading on plate reader. Read absorbance at 450 nm and 590 nm in a plate reader within 5 minutes and ensure that there are no air bubbles in any well. Record the difference in absorbance units.

# Assay Procedure I for Human Leptin ELISA Kit

	Step 1	Step 2	Step 3-4	Step 5	Step 6-8	Step 9	Step 9-10	Step 11	Step 11-13	Step 14	Step 14	Step 15	Step 15
Well #			Assay Buffer	Standards/ Controls/ Samples		Detection Antibody		Enzyme Solution		Substrate		Stop Solution	
A1, B1	Dilute 2 bottles of 10X Wash Buffer with 900 mL Deionized Water.  Wash plate with 300 µL Wash Buffer and incubate at room temperature for 5 minutes. Remove residual buffer by tapping smartly on absorbent towels.	75 µL	100 µL	--	Seal, Agitate, Incubate 2 hours at Room Temperature. Wash 3X with 300 µL Wash Buffer.	100 µL	Seal, Agitate, Incubate 30 minutes at Room Temperature. Remove residual buffer by tapping smartly on absorbent towels	100 µL	Seal, Agitate, Incubate 30 minutes at Room Temperature on a plate shaker. Wash 5X with 300 µL Wash Buffer.	100 µL	Seal, Agitate, Incubate for 5-20 minutes at Room Temperature.	100 µL	Read Absorbance at 450 nm and 590 nm.
C1, D1			25 µL of 0.78 ng/mL Standard	25 µL		100 µL							
E1, F1			25 µL of 1.56 ng/mL Standard										
G1, H1			25 µL of 3.125 ng/mL Standard										
A2, B2			25 µL of 6.25 ng/mL Standard										
C2, D2			25 µL of 12.5 ng/mL Standard										
E2, F2			25 µL of 25 ng/mL Standard										
G2, H2			25 µL of 50 ng/mL Standard										
A3, B3			25 µL of 100 ng/mL Standard										
C3, D3			25 µL of QC I										
E3, F3			25 µL of QC II										
G3, H3.			25 µL of Sample										
A4, B4	25 µL of Sample												

For research use only. Not for use in diagnostic procedures.

## Microtiter Plate Arrangement

Human Leptin ELISA

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blank	6.25 ng/mL	100 ng/mL	Sample 2								
B	Blank	6.25 ng/mL	100 ng/mL	Sample 2								
C	0.78 ng/mL	12.5 ng/mL	QC 1	Etc.								
D	0.78 ng/mL	12.5 ng/mL	QC 1									
E	1.56 ng/mL	25 ng/mL	QC 2									
F	1.56 ng/mL	25 ng/mL	QC 2									
G	3.125 ng/mL	50 ng/mL	Sample 1									
H	3.125 ng/mL	50 ng/mL	Sample 1									

For research use only. Not for use in diagnostic procedures.

## Human Leptin ELISA (Sensitive) Assay Procedure II

Sensitivity: 0.195 ng/mL–25 ng/mL

### Human Leptin (Sensitive) Standard Preparation

**Note:** The standard curve range recommended for samples with low Leptin levels is 0.195 ng/mL to 25 ng/mL. By performing 1:1 serial dilutions of the 0.78 ng/mL kit standard with Assay Buffer, the 0.39 ng/mL and 0.195 ng/mL standard points can be created for use in the Leptin Sensitive Assay.

1. Label two tubes 0.39 and 0.195 ng/mL. Add 0.2 mL Assay Buffer to each of the two tubes. Prepare serial dilutions by adding 0.2 mL of the 0.78 ng/mL standard to the 0.39 ng/mL tube, mix well and transfer 0.2 mL of the 0.39 ng/mL standard to the 0.195 ng/mL tube and mix well.

**Note:** Do not use a Repeater pipette. Change tip for every dilution. Wet tip with Standard before dispensing. Unused portions of standard should be stored at 2–8 °C. Avoid multiple freeze/thaw cycles.

<b>Standard Concentration ng/mL</b>	<b>Volume of Assay Buffer to Add</b>	<b>Volume of Standard to Add</b>
0.39	0.2 mL	0.2 mL of 0.78 ng/mL
0.195	0.2 mL	0.2 mL of 0.39 ng/mL

## Human Leptin (Sensitive) Quality Control 1 and 2 Preparation

1. Prepare 1:3 dilutions of Quality Control 1 and Quality Control 2 for use in the Human Leptin (Sensitive) assay. Label two tubes QC1 and QC2. Add 0.4 mL Assay Buffer to each of the tubes. Add 0.2 mL Quality Control 1 to the QC1 tube and mix well. Add 0.2 mL Quality Control 2 to the QC2 tube and mix well.

**Note:** Do not use a Repeater pipette. Change tip for every dilution. Wet tip with Quality Control before dispensing. Unused portions of Quality Control should be stored at 2–8 °C. Avoid multiple freeze/thaw cycles.

	<b>Volume of Assay Buffer to Add</b>	<b>Volume of Quality Control to Add</b>
<b>QC1</b>	0.4 mL	0.2 mL of Quality Control 1
<b>QC2</b>	0.4 mL	0.2 mL of Quality Control 2

## Human Leptin (Sensitive) Assay

(Sensitivity: 0.195 ng/mL–25 ng/mL)

Pre-warm all reagents to room temperature immediately before setting up the assay.

1. Dilute the concentrated Wash Buffer 10-fold by adding the entire contents of both bottles of buffer to 900 mL de-ionized or glass distilled water.
2. Remove required number of strips from the Microtiter Assay Plate. Unused strips should be resealed in the foil pouch with the desiccant provided and stored at 2–8 °C. Assemble strips in an empty plate holder and add 300 µL of diluted Wash Buffer to each well. Incubate at room temperature for 5 minutes. Decant wash buffer and remove the residual amount from all wells by inverting the plate and tapping it smartly onto absorbent towels several times. Do not let wells dry before proceeding to the next step. If an automated machine is used for the assay, follow the manufacturer's instructions for all washing steps described in this protocol.
3. Add 50 µL Assay Buffer into all wells.
4. Add in duplicate 50 µL Assay Buffer to blank wells. (Refer to [Microtiter Plate Arrangement II](#) for suggested well orientations.)
5. Add in duplicate 50 µL Human Leptin Standards in order of ascending concentration to the appropriate wells. Add in duplicate 50 µL QC1 and 50 µL QC2 to the appropriate wells. Add sequentially 50 µL of samples in duplicate to the remaining wells. **For best results all additions should be completed within one hour.**
6. Cover the plate with plate sealer and incubate at room temperature for 2 hours on an orbital microtiter plate shaker set to rotate at moderate speed, about 400 to 500 rpm.
7. Remove plate sealer and decant solutions from the plate. Tap as before to remove residual solutions in the wells.
8. Wash wells 3 times with diluted Wash Buffer, 300 µL per well per wash. Decant and tap after each wash to remove residual buffer.
9. Add 100 µL Detection Antibody to each well. Cover the plate with sealer and incubate at room temperature for 30 minutes on the microtiter plate shaker.
10. Remove sealer and decant solution from the plate. Tap as before to remove residual solutions in the wells.
11. Add 100 µL Enzyme Solution to each well. Cover plate with sealer and incubate with moderate shaking at room temperature for 30 minutes on the microtiter plate shaker.
12. Remove sealer, decant solution from the plate, and tap plate to remove the residual fluid.
13. Wash wells 5 times with diluted Wash Buffer, 300 µL per well per wash. Decant and tap firmly after each wash to remove residual buffer.

For research use only. Not for use in diagnostic procedures.

- 
14. Add 100  $\mu\text{L}$  of Substrate Solution to each well, cover plate with sealer and shake on the plate shaker for approximately 5-20 minutes. Blue color should be formed in wells of Leptin standards with intensity proportional to increasing concentrations of Leptin.

**Note:** Please be aware that the color may develop more quickly or more slowly than the recommended incubation time depending on the localized room temperature. Please visually monitor the color development to optimize the incubation time. One can monitor color development using 370 nm filter, if available on the spectrophotometer. When the absorbance is between 1.2 and 1.8 at 370 nm, the stop solution can be added to terminate the color development.

15. Remove sealer and add 100  $\mu\text{L}$  of Stop Solution (**Caution: Corrosive solution**) and shake plate by hand to ensure complete mixing of solution in all wells. The blue color should turn to yellow after acidification. Wipe the bottom of the microtiter plate to remove any residue prior to reading on plate reader. Read absorbance at 450nm and 590nm in a plate reader within 5 minutes and ensure that there are no air bubbles in any well. Record the difference in absorbance units.

# Assay Procedure II for Human Leptin (Sensitive) ELISA kit

	Step 1	Step 2	Step 3-4	Step 5	Step 6-8	Step 9	Step 9-10	Step 11	Step 11-13	Step 14	Step 14	Step 15	Step 15
Well #			Assay Buffer	Standards/ Controls/ Samples		Detection Antibody		Enzyme Solution		Substrate		Stop Solution	
A1, B1			100 µL	--		100 µL		100 µL		100 µL		100 µL	
C1, D1			50 µL	50 µL of 0.195 ng/mL Standard		↓		↓		↓		↓	
E1, F1				50 µL of 0.39 ng/mL Standard									
G1, H1				50 µL of 0.78 ng/mL Standard									
A2, B2				50 µL of 1.56 ng/mL Standard									
C2, D2				50 µL of 3.12 ng/mL Standard									
E2, F2				50 µL of 6.25 ng/mL Standard									
G2, H2				50 µL of 12.5 ng/mL Standard									
A3, B3				50 µL of 25 ng/mL Standard									
C3, D3				50 µL of diluted QC I									
E3, F3				50 µL of diluted QC II									
G3, H3				50 µL of Sample									
A4, B4				50 µL of Sample									
	Dilute 2 bottle of 10X Wash Buffer with 900 mL Deionized Water.		Wash plate with 300 µL Wash Buffer and incubate at room temperature for 5 minutes. Remove residual buffer by tapping smartly on absorbent towels.		Seal, Agitate, Incubate 2 hours at Room Temperature. Wash 3X with 300 µL Wash Buffer.		Seal, Agitate, Incubate 30 minutes at Room Temperature. Remove residual buffer by tapping smartly on absorbent towels		Seal, Agitate, Incubate 30 minutes at Room Temperature. Wash 5X with 300 µL Wash Buffer.		Seal, Agitate, Incubate 5–20 minutes at Room Temperature.		Read Absorbance at 450 nm and 590 nm

For research use only. Not for use in diagnostic procedures.

## Microtiter Plate Arrangement II

### Human Leptin (Sensitive) ELISA

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blank	1.56 ng/mL	25.0 ng/mL	Sample 2								
B	Blank	1.56 ng/mL	25.0 ng/mL	Sample 2								
C	0.195 ng/mL	3.12 ng/mL	QC 1	Etc.								
D	0.195 ng/mL	3.12 ng/mL	QC 1									
E	0.39 ng/mL	6.25 ng/mL	QC 2									
F	0.39 ng/mL	6.25 ng/mL	QC 2									
G	0.78 ng/mL	12.5 ng/mL	Sample 1									
H	0.78 ng/mL	12.5 ng/mL	Sample 1									

For research use only. Not for use in diagnostic procedures.

---

## Calculations

The dose-response curve of this assay fits best to a sigmoidal 5-parameter logistic equation. The results of unknown samples can be calculated with any computer program having a 5-parameter logistic function.

**Note:** When sample volumes assayed differ from 25  $\mu\text{L}$  (in normal assay), an appropriate mathematical adjustment must be made to accommodate for the dilution factor (for example, if 12.5  $\mu\text{L}$  of sample is used, then calculated data must be multiplied by 2).

## Interpretation

- The assay should be rejected if one of the two QCs falls outside of 2 standard deviations of the applicable mean.
- The limit of sensitivity of this assay is 0.2 ng/mL + 2 SD human leptin (25  $\mu\text{L}$  sample size).
- The appropriate range of this assay is 0.78 to 100 ng/mL human leptin (25  $\mu\text{L}$  sample size). Any result greater than 100 ng/mL in a 25  $\mu\text{L}$  sample assayed should be diluted and repeated using assay buffer as diluent until it falls within range.

## Normal Range

Normal range: Leptin levels are directly correlated with degree of adiposity.

Mean Serum Leptin levels from normal lean individuals (BMI ranges 18-25):

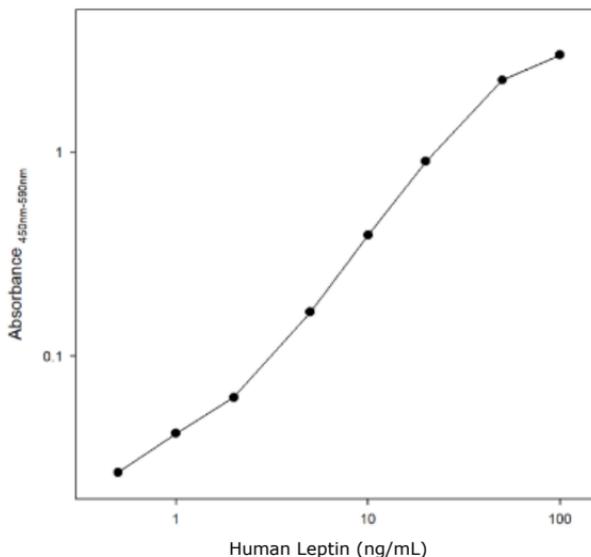
Lean Men:  $3.8 \pm 1.8$  ng/mL

Lean Women:  $7.4 \pm 3.7$  ng/mL

Serum levels are approximately 2.5 times higher in women per unit BMI as compared to men.

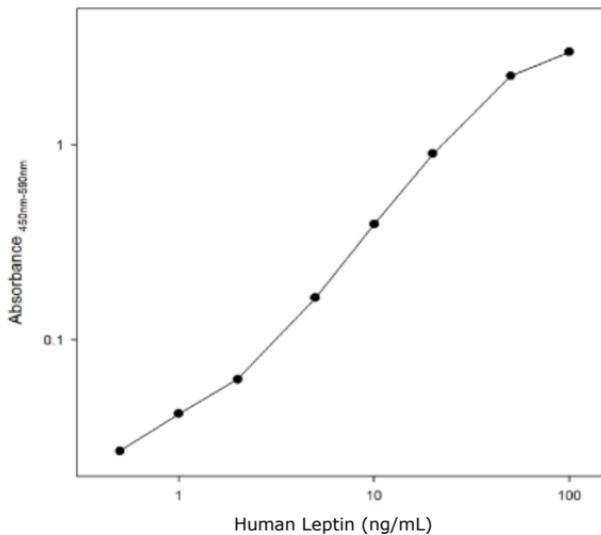
## Standard Curve

### Human Leptin ELISA



Typical Standard Curve – Not to be used to calculate data

### Sensitive Human Leptin



Typical Standard Curve – Not to be used to calculate data

For research use only. Not for use in diagnostic procedures.

## Assay Characteristics

### Sensitivity

The lowest level of human leptin that can be detected by this assay is 0.2 ng/mL + 2 SD (25 µL sample size). For the sensitive assay, the lowest level of human leptin that can be detected is 0.2 ng/mL + 2 SD (50 µL sample size).

### Specificity

The antibody pair used in this assay is specific human leptin and does not significantly cross-react to the following molecules/hormones tested:

Human Leptin	100%
Chicken Leptin	n.d.*
Mouse Leptin	n.d.*
Ovine Leptin	n.d.*
Porcine Leptin	n.d.*
Rat Leptin	n.d.*
Bovine Proinsulin	n.d.*
Human Proinsulin	n.d.*
Porcine Proinsulin	n.d.*
Human Insulin	n.d.*
Human IGF-I	n.d.***
Human IGF-II	n.d.***
Glucagon	n.d.*

n.d.: Not detectable at concentrations up to \* 625 nM; \*\* 400 nM and \*\*\* 200 nM.

## Precision

	<b>Mean Leptin Levels (ng/mL)</b>	<b>Intra-Assay % CV</b>	<b>Inter-Assay % CV</b>
1	2.34	4.6	6.2
2	9.53	3.4	4.2
3	22.08	4.1	3.0
4	28.9	2.6	2.6

## Sensitive Assay

	<b>Mean Leptin Levels (ng/mL)</b>	<b>Intra-Assay % CV</b>	<b>Inter-Assay % CV</b>
1	0.86	4.9	8.6
2	4.59	2.2	3.4
3	11.26	1.4	1.7
4	14.08	1.9	1.3

The assay variations of Human Leptin ELISA kits were studied on four human serum samples with varying concentrations of endogenous leptin. The mean within variations of four duplicate determinations in each assay of the indicated samples were calculated from results of three separate assays. The mean between variation of each sample was calculated from results of three separate assays with four duplicate samples in each assay.

## Spike Recovery of Human Leptin in Assay Samples

Varying amounts of human leptin were added to three human serum samples and the leptin content was determined in three separate assays.

The % of recovery = observed leptin concentrations/expected leptin concentrations x 100%.

<b>Sample</b>	<b>Leptin Added (ng/mL)</b>	<b>Expected (ng/mL)</b>	<b>Observed (ng/mL)</b>	<b>Recovery</b>
1	0	4.4	4.4	100
	5	9.4	9.84	104.7
	20	24.4	27.2	111.5
	50	54.4	51.89	95.4
2	0	12.02	12.02	100
	5	17.02	18.91	111.1
	20	32.02	36.54	114.1
	50	62.02	61.61	99.3
3	0	32.79	32.79	100
	5	37.79	39.95	105.7
	20	52.79	52.79	100
	50	82.79	94.33	113.9

## Spike Recovery of Human Leptin in Sensitive Assay Samples

<b>Sample</b>	<b>Leptin Added (ng/mL)</b>	<b>Expected (ng/mL)</b>	<b>Observed (ng/mL)</b>	<b>Recovery</b>
1	0	0.87	0.87	100.0
	1	1.87	1.8	96.3
	5	5.87	5.7	97.1
	10	10.87	10.4	95.7
2	0	3.59	3.59	100.0
	1	4.59	4.58	99.8
	5	8.59	8.78	102.2
	10	13.59	12.98	95.5
3	0	7.21	7.21	100.0
	1	8.21	8.15	99.3
	5	12.21	11.53	94.4
	10	17.21	14.83	86.2

## Linearity of Sample Dilution

### Effect of Serum Dilution

Three human serum samples with the indicated sample volumes were assayed in three separate experiments. Required amounts of assay buffer were added to compensate for lost volumes below 25  $\mu\text{L}$ .

$\% \text{ expected} = (\text{observed}/\text{expected}) \times 100\%$

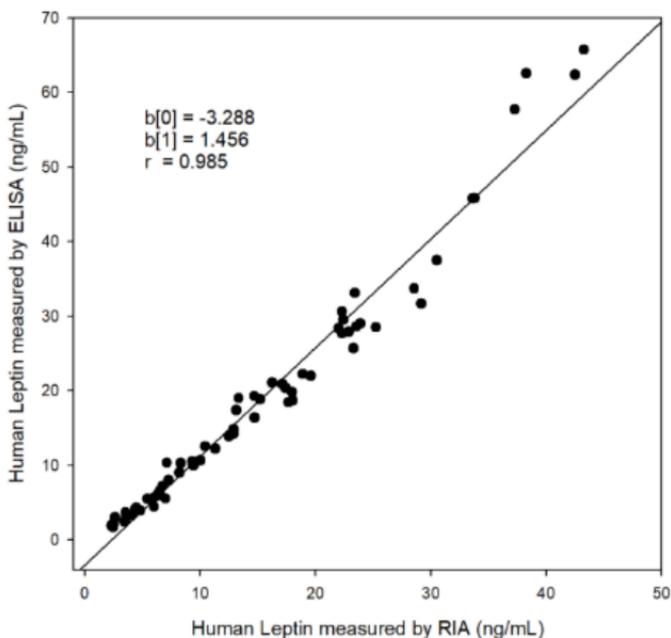
<b>Sample</b>	<b>Volume (<math>\mu\text{L}</math>)</b>	<b>Expected (ng/mL)</b>	<b>Observed (ng/mL)</b>	<b>Expected</b>
1	25	61.31	61.31	100
	12.5	-	32.22	105.1
	5	-	11.73	95.7
	2.5	-	5.62	91.7
2	25	33.05	33.05	100
	12.5	-	15	90.8
	5	-	5.55	80.4
	2.5	-	2.65	80.2
3	25	15.21	15.21	100
	12.5	-	7.35	96.6
	5	-	2.6	85.5
	2.5	-	1.23	80.9

## Linearity of Sample Dilution (Sensitive Assay)

Sample	Volume (µL)	Expected (ng/mL)	Observed (ng/mL)	Expected
1	50	3.50	3.50	100
	25	-	1.88	107.3
	12.5	-	1.00	114.4
	6.25	-	0.51	117.2
	3.125	-	0.24	108.3
	1.56	-	0.12	106.9
2	50	5.08	5.08	100
	25	-	2.70	106.1
	12.5	-	1.40	109.9
	6.25	-	0.78	122.1
	3.125	-	0.39	122.9
	1.56	-	0.16	100.8
3	50	6.54	6.54	100
	25	-	3.15	96.3
	12.5	-	1.73	105.6
	6.25	-	0.97	118.0
	3.125	-	0.47	114.5
	1.56	-	0.26	125.3

## Correlation Graph

### Human Leptin Correlation RIA vs. ELISA



66 serum samples collected from 66 human subjects were assayed for leptin content using both the Human Leptin RIA Kit (Cat. No. EZHL-81K) and the Human Leptin ELISA Kit (Cat. No. EZHL-80SK). Correlation of the two assay kits derived by linear regression analysis of paired results from each sample.

---

## Quality Controls

The ranges for each analyte in Quality Control 1 and 2 are provided on the card insert, or available at our website [SigmaAldrich.com](http://SigmaAldrich.com).

## Troubleshooting

- To obtain reliable and reproducible results the operator should carefully read this manual and fully understand all aspects of each assay step before attempting to run the assay.
- Throughout the assay the operator should adhere strictly to the procedures with good laboratory practice.
- Have all necessary reagents and equipment ready on hand before starting. Once the assay has been started all steps should be completed with precise timing and without interruption.
- Avoid cross contamination of any reagents or samples to be used in the assay.
- Make sure all reagents and samples are added to the bottom of each well.
- Careful and complete mixing of solutions in the well is critical. Poor assay precision will result from incomplete mixing or cross well contamination due to inappropriate mixing.
- Remove any air bubbles formed in the well after acidification of substrate solution because bubbles interfere with spectrophotometric readings.
- Do not let the absorbency reading of the highest standard reach 2.0 units or higher before adding the stop solution.
- High absorbance in the background or blank wells could be due to:
  - Well cross contamination by standard solution or sample.
  - Inadequate washing of the well with HRP.

## Product Ordering

Products are available for online ordering at [SigmaAldrich.com](https://www.sigmaaldrich.com).

### Replacement Reagents

<b>Reagents</b>	<b>Cat. No.</b>
Human Leptin ELISA Plate	EP81
10X HRP Wash Buffer Concentrate	EWB-HRP
Human Leptin ELISA Standards	E8081-K
Quality Controls 1 and 2	E6081-K
Assay Buffer for assay	EABTR
Human Leptin Detection Antibody	E1081
Enzyme Solution	EHRP-3
Substrate	ESS-TMB
Stop Solution	ET-TMB
10 pack of Human Leptin ELISA Kits	EZHL-80BK

---

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

## Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://SigmaAldrich.com/techservice).

## Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://SigmaAldrich.com/terms).

## Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://SigmaAldrich.com/offices).

---

The life science business of Merck operates  
as MilliporeSigma in the U.S. and Canada.

Merck, Millipore and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2007-2024 Merck KGaA, Darmstadt, Germany and/or its affiliates.  
All Rights Reserved.

For research use only. Not for use in diagnostic procedures.

The Merck logo is displayed in a bold, blue, sans-serif font. The letters are closely spaced and have a slightly irregular, hand-drawn appearance.