

User Guide

# Rat/Mouse Ghrelin (Active) ELISA Kit

## 96-Well Plate

### EZRGRA-90K

Intended Use .....	2	Graph of Typical Reference Curve ..	14
Principles of Assay .....	2	Assay Characteristics .....	15
Reagents Supplied .....	3	Sensitivity .....	15
Storage and Stability .....	4	Specificity .....	15
Reagent Precautions .....	4	Precision .....	16
Sodium Azide .....	4	Spike Recovery of Rat/Mouse Ghrelin (Active) in Assay Samples .....	17
Hydrochloric Acid .....	4	Linearity of Sample Dilution .....	18
Materials Required .....	6	Normal Range of Active Ghrelin Levels in Rat/Mouse Blood .....	19
Sample Collection and Storage .....	7	Post Prandial Attenuation of Active Ghrelin in Blood .....	20
Reagent Preparation .....	8	Effects of Fasting on Serum/Plasma Ghrelin (Active) Levels .....	20
Rat/Mouse Ghrelin (Active) Standard Preparation .....	8	RIA Cat. No. GHRA-88HK vs. ELISA Cat. No. EZRGRA-90K .....	21
Quality Control 1 and 2 Preparation .....	9	Quality Controls .....	22
Preparation of Capture and Detection Antibody Mixture .....	9	Troubleshooting .....	22
Rat/Mouse Ghrelin (Active) ELISA Assay Procedure .....	9	Product Ordering .....	23
Assay Procedure for Rat/Mouse Ghrelin (Active) ELISA Kit .....	11	Replacement Reagents .....	23
Microtiter Plate Arrangement .....	12	Notice .....	24
Calculations .....	13	Technical Assistance .....	24
Interpretation .....	13	Terms and Conditions of Sale .....	24
		Contact Information .....	24

---

## Intended Use

This kit is used for the non-radioactive quantification of rat/mouse ghrelin (active) in serum and plasma. There is no cross reactivity to des octanoyl-ghrelin. Circulating ghrelin is a multifunctional hormone produced primarily by the stomach. It consists of 28 amino acids and the n-octanoylation of serine3 position in the molecule is necessary for its bioactivity. Originally found as an endogenous ligand for the growth hormone secretagogue receptor in the pituitary gland, it distinguishes itself from the hypothalamic growth hormone-releasing hormone as another potent stimulator for growth hormone secretion. It is also an important orexigenic hormone in the regulation of energy homeostasis. One kit is sufficient to measure 38 unknown samples in duplicate.

**For Research Use Only. Not for Use in Diagnostic Procedures.**

## Principles of Assay

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

- Capture of ghrelin molecules (active form) in the sample by anti-ghrelin IgG and immobilization of the resulting complex to the wells of a microtiter plated coated by a pre-titered amount of anchor antibodies
- Binding of a second biotinylated antibody to the captured molecules
- Washing of unbound materials from samples
- Conjugation of horseradish peroxidase to the immobilized biotinylated antibodies
- Washing of excess free enzyme conjugates
- Quantification of immobilized antibody-enzyme conjugates by monitoring 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, corrected from the absorbency at 590 nm, after acidification of formed products. Since the increase in absorbance is directly proportional to the amount rat/mouse ghrelin (active form) in the unknown sample, the concentration of active ghrelin can be derived by interpolation from a reference curve generated in the same assay with reference standards of known concentrations of rat/mouse ghrelin.

# 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.
Microtiter Plate with 2 plate sealers coated with pre-titered anchor antibodies. <b>Note:</b> Unused strips should be resealed in the foil pouch with the desiccant provided and stored at 2-8 °C	-	1 plate 2 sealers	EPDAR
Rat/Mouse Ghrelin (Active) Standard	2 mL/vial Lyophilized	1 vial	E8090-K
Quality Controls 1 and 2	0.5 mL/vial upon hydration Lyophilized	1 vial each	E6090-K
Matrix Solution	1 mL	1 vial	EMTX-GA
Assay Buffer	15 mL	1 bottle	EABGR
10X HRP Wash Buffer Concentrate.	50 mL	2 bottles	EWB-HRP
Rat/Mouse Ghrelin (Active) Detection Antibody	3 mL	1 bottle	E1090-D
Rat/Mouse Ghrelin (Active) Capture Antibody	3 mL	1 bottle	E1090-C
Enzyme Solution	12 mL	1 bottle	EHRP-88
Substrate Solution <b>Note:</b> Minimize light exposure.	12 mL	1 bottle	ESS-TMB2
Stop Solution <b>Caution:</b> Corrosive Solution	12 mL	1 bottle	ET-TMB

---

## Storage and Stability

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

All components are shipped and stored at 2-8 °C. Reconstituted standards and controls can be frozen for future use but repeated freeze/thaw cycles should be avoided. 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. Harmful if swallowed. Causes respiratory and digestive tract burns. Avoid contact with skin and eye. Do not swallow or ingest.

**Note:** See full labels of Hazardous Components on the next page.

## Symbol Definitions

Ingredient	Cat. No.	Full Label
Rat/Mouse Ghrelin (Active) Capture Antibody	E1090-C	 <p><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.</p>
Rat/Mouse Ghrelin (Active) Detection Antibody	E1090-D	 <p><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.</p>
Rat/Mouse Ghrelin (Active) Quality Controls 1 & 2	E6090-K	  <p><b>Danger:</b> Harmful if swallowed. Causes serious eye damage. Harmful to aquatic life with long lasting effects. Avoid release to the environment. Wear eye protection. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/attention.</p>
Rat/Mouse Ghrelin (Active) Standard	E8090-K	  <p><b>Danger:</b> Harmful if swallowed. Causes serious eye damage. Harmful to aquatic life with long lasting effects. Avoid release to the environment. Wear eye protection. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/attention.</p>
Assay Buffer	EABGR	 <p><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.</p>
Stop Solution	ET-TMB	 <p><b>Warning:</b> May be corrosive to metals.</p>
10X HRP Wash Buffer Concentrate	EWB-HRP	 <p><b>Warning.</b> May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.</p>

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

---

## Materials Required (Not Provided)

- Multi-channel Pipettes and pipette tips: 5  $\mu$ L-50  $\mu$ L and 50  $\mu$ L-300  $\mu$ L
- Pipettes and pipette tips: 10  $\mu$ L-20  $\mu$ L or 20  $\mu$ L-100  $\mu$ L
- Buffer and Reagent Reservoirs
- Vortex Mixer
- De-ionized water
- Microtiter Plate Reader capable of reading absorbency at 450 nm and 590 nm
- Orbital Microtiter Plate Shaker
- Absorbent Paper or Cloth
- Pefabloc® or AEBSF [4-(2-Aminoethyl)-benzenesulfonyl fluoride], 100 mg/mL aqueous stock solution (store at -20 °C, minimize multiple freeze/thaw cycles) is recommended for Sample Collection and Storage.
- 5 N HCl, recommended for Sample Collection and Storage.

---

## Sample Collection and Storage

The active ghrelin molecule is extremely unstable in serum/plasma and should be rigorously protected during blood sample collection. Ideally all samples should be processed as quickly as possible and kept on ice to retard the breakdown of active ghrelin. For maximum protection, we recommend addition of Pefabloc® or AEBSF and acidification of all samples. Acidification will result in noticeable protein precipitation but does not affect the assay. However, if the presence of precipitates interferes with the sample pipetting accuracy, the sample should be centrifuged and the supernatant used for assay.

1. To prepare serum, whole blood is directly drawn into a centrifuge tube that contains no anti-coagulant. Immediately add enough Pefabloc® or AEBSF to a final concentration of 1 mg/mL. Let blood clot at room temperature for 30 min.
2. Promptly centrifuge the clotted blood at 2,000 to 3,000 x g for 15 minutes at 4 ± 2 °C.
3. Transfer serum samples in separate tubes and acidify with HCl to a final concentration of 0.05 N. Aliquot acidified serum in small quantities. Date and identify each sample.
4. Use freshly prepared serum or store samples at -20 ± 5 °C for later use. Avoid multiple (> 5) freeze/thaw cycles.
5. To prepare plasma sample, whole blood should be collected into a centrifuge tube containing enough K<sub>3</sub>EDTA to achieve a final concentration of 1.735 mg/mL and treated with Pefabloc® or AEBSF as described for serum, followed by immediate centrifugation. Acidify plasma samples with HCl to a final concentration of 0.05 N. Observe same precautions in the preparation of serum samples.
6. If heparin is to be used as 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.

## Reagent Preparation

### Rat/Mouse Ghrelin (Active) Standard Preparation

1. Use care in opening the lyophilized Standard vial. Using a pipette, reconstitute the Rat/Mouse Ghrelin (Active) Standard with 2 mL of deionized water. Please refer to the analysis sheet for exact concentration. Invert and mix gently until completely in solution.
2. Label six tubes 1, 2, 3, 4, 5, and 6. Add Assay Buffer to each of the six tubes according to the volumes outlined in the chart below. Dilute the reconstituted standard stock according to the chart below. Vortex each tube briefly to ensure complete mixing.

**Note:** Change tip for every dilution. Wet tip with standard before dispensing. Unused portions of reconstituted standard should be stored in small aliquots at  $\leq -20^{\circ}\text{C}$ . Avoid multiple freeze/thaw cycles.

Volume of Deionized Water to Add	Volume of Standard to Add	Standard Stock Concentration
2 mL	0	X (refer to analysis sheet for exact concentration)

Tube #	Volume of Assay Buffer to Add	Volume of Standard to Add	Standard Concentration (pg/mL)
1	500 $\mu\text{L}$	500 $\mu\text{L}$ of reconstituted standard	X/2
2	500 $\mu\text{L}$	500 $\mu\text{L}$ of Tube 1	X/4
3	500 $\mu\text{L}$	500 $\mu\text{L}$ of Tube 2	X/8
4	500 $\mu\text{L}$	500 $\mu\text{L}$ of Tube 3	X/16
5	500 $\mu\text{L}$	500 $\mu\text{L}$ of Tube 4	X/32
6	500 $\mu\text{L}$	500 $\mu\text{L}$ of Tube 5	X/64



## Quality Control 1 and 2 Preparation

Use care in opening the lyophilized Quality Control vials. Reconstitute each Quality Control 1 and Quality Control 2 with 0.5 mL distilled or de-ionized water and gently invert to ensure complete hydration. Unused portions of the reconstituted Quality Controls should be stored in small aliquots at  $\leq -20^{\circ}\text{C}$ . Avoid further freeze/thaw cycles.

## Preparation of Capture and Detection Antibody Mixture

Prior to use, combine the entire contents of Rat/Mouse Ghrelin (Active) Capture Antibody (3 mL) and Rat/Mouse Ghrelin (Active) Detection Antibody (3 mL), or at a 1:1 ratio, and invert to mix thoroughly.

## Rat/Mouse Ghrelin (Active) ELISA Assay Procedure

Warm all reagents to room temperature before setting up the assay.

1. Dilute 10X concentrated HRP wash buffer 10-fold by mixing the entire contents of both buffer bottles with 900 mL de-ionized or glass distilled water.
2. Remove the required number of strips from the Microtiter Assay Plate. Unused strips should be resealed in the foil pouch and stored at 2-8 °C. Assemble the strips in an empty plate holder. Add 300  $\mu\text{L}$  diluted Wash Buffer to each well of the plate. Decant Wash Buffer and remove the residual volume by inverting the plate and tapping it smartly onto absorbent towels several times. Repeat wash procedure 2 additional 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 20  $\mu\text{L}$  of appropriate Matrix Solution to Blank, Standards and Quality Control wells (refer to [Microtiter Plate Arrangement](#)).
4. Add 30  $\mu\text{L}$  Assay Buffer to each of the Blank and Sample wells.
5. Add 10  $\mu\text{L}$  Assay Buffer to each of the Standard and Quality Control wells.
6. Add in duplicate 20  $\mu\text{L}$  Ghrelin Standards in order of ascending concentrations to the appropriate wells.
7. Add in duplicate 20  $\mu\text{L}$  QC1 and 20  $\mu\text{L}$  QC2 to the appropriate wells.
8. Add sequentially 20  $\mu\text{L}$  of the unknown samples in duplicate to the remaining wells.
9. Transfer the Antibody Solution Mixture (1:1 mixture of capture and detection antibody) to a buffer or reagent reservoir and add 50  $\mu\text{L}$  to each well with a multi-channel pipette.
10. 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.

- 
11. Remove plate sealer and decant solutions from the plate. Tap as before to remove residual solutions in well.
  12. 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.
  13. Add 100  $\mu$ L Enzyme Solution to each well. Cover plate with sealer and incubate with moderate shaking at room temperature for 30 min on the micro-titer plate shaker.
  - 14.** Remove sealer, decant solutions from the plate and tap plate to remove the residual fluid.
  15. Wash wells 6 times with diluted Wash Buffer, 300  $\mu$ L per well per wash. Decant and tap after each wash to remove residual buffer.
  16. Add 100  $\mu$ L of Substrate Solution to each well, cover plate with sealer and shake in the plate shaker for approximately 5-20 minutes. Blue color should be formed in wells of Ghrelin standards with intensity proportional to increasing concentrations of Ghrelin.  
  
**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.
  17. Remove sealer and add 100  $\mu$ L Stop Solution (**Caution:** Corrosive Solution) and shake plate by hand to ensure complete mixing of solution in all wells. The blue color should turn into 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.

# Assay Procedure for Rat/Mouse Ghrelin (Active) ELISA Kit

	Step 1	Step 2	Step 3	Step 4-5	Step 6-8	Step 9	Step 10-12	Step 13	Step 14-15	Step 16	Step 17
Well #	Dilute both bottles of 10X HRP Wash Buffer with 900 mL de-ionized water.	Wash plate 3X with 300 µL Wash Buffer. Remove residual buffer by tapping smartly on absorbent towels.	Matrix Solution	Assay Buffer	Standards/ QCs/Samples	Capture/Detecti on Ab Mixture	Seal, Agitate, Incubate 2 hours at Room Temperature. Wash 3X with 300 µL Wash Buffer.	Enzyme Solution	Seal, Agitate, Incubate 30 minutes at Room Temperature. Wash 6X with 300 µL Wash Buffer.	Substrate	Seal, Agitate, Incubate 15 minutes at Room Temperature.  Read absorbance at 450 nm and 590 nm.
A1, B1			20 µL	30 µL	-	50 µL		100 µL		100 µL	
C1, D1			20 µL	10 µL	20 µL of Tube 6	↓		↓		↓	
E1, F1			20 µL	10 µL	20 µL of Tube 5						
G1, H1			20 µL	10 µL	20 µL of Tube 4						
A2, B2			20 µL	10 µL	20 µL of Tube 3						
C2, D2			20 µL	10 µL	20 µL of Tube 2						
E2, F2			20 µL	10 µL	20 µL of Tube 1						
G2, H2			20 µL	10 µL	20 µL Reconstituted standard						
A3, B3			20 µL	10 µL	20 µL of QC 1						
C3, D3			20 µL	10 µL	20 µL of QC 2						
E3, F3			-	30 µL	20 µL of Sample 1						
G3, H3, etc.	-	30 µL	20 µL of Sample 2								

## Microtiter Plate Arrangement

### Rat/Mouse Ghrelin (Active) ELISA

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blank	Tube 3 Standard	QC 1	Etc.								
B	Blank	Tube 3 Standard	QC 1	Etc.								
C	Tube 6 Standard	Tube 2 Standard	QC 2									
D	Tube 6 Standard	Tube 2 Standard	QC 2									
E	Tube 5 Standard	Tube 1 Standard	Sample 1									
F	Tube 5 Standard	Tube 1 Standard	Sample 1									
G	Tube 4 Standard	Reconstituted Standard	Sample 2									
H	Tube 4 Standard	Reconstituted Standard	Sample 2									

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

---

## Calculations

Graph a reference curve by plotting the absorbance unit of 450 nm, less unit at 590 nm, on the Y-axis against the concentrations of Ghrelin standard on the X-axis. The dose-response curve of this assay fits best to a sigmoidal 4- or 5-parameter logistic equation. The results of unknown samples can be calculated with any computer program having a 4- or 5-parameter logistic function.

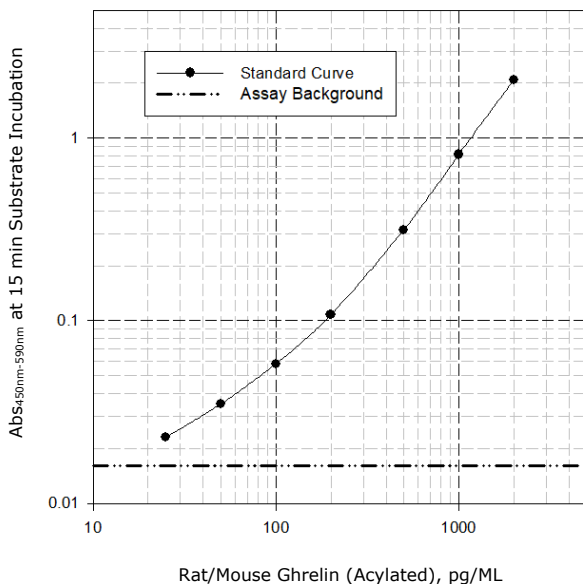
**Note:** When sample volumes assayed differ from 20  $\mu$ L, an appropriate mathematical adjustment must be made to accommodate for the dilution factor (for example, if 10  $\mu$ L of sample is used, then calculated data must be multiplied by 2). When sample volume assayed is less than 20  $\mu$ L, compensate the volume deficit with matrix solution.

## Interpretation

1. The assay will be considered accepted when all Quality Control values fall within the calculated QC range. If any QCs fall outside of the control range, review results with a supervisor.
2. If the difference between duplicate results of a sample is > 15% CV, repeat the sample.
3. The theoretical minimal detecting concentration of this assay is 8 pg/mL Active Ghrelin (20  $\mu$ L sample size).
4. The appropriate range of this assay is 25 pg/mL to 2,000 pg/mL Active Ghrelin (20  $\mu$ L sample size). Any result greater than 2,000 pg/mL in a 20  $\mu$ L sample should be diluted using matrix solution and the assay repeated until the results fall within range.

## Graph of Typical Reference Curve

Typical Standard Curve, not to be used to calculate data.



For demonstration only. Do not use for calculations

# Assay Characteristics

## Sensitivity

The Minimum Detectable Concentration (MinDC) of Active Ghrelin is 8 pg/mL when using a 20 µL sample size.

## Specificity

Rat/Mouse Ghrelin (Active)	100%
Des-Octanoyl Rat/Mouse Ghrelin	0%
Canine Ghrelin (Active)	111%
Porcine Ghrelin (Active)	98%
Human Ghrelin (Active)	53%
Des-Octanoyl Human Ghreline	0%
Motilin Related Peptide (Human, Rat/Mouse)	0%
PYY 3-36 (Human, Mouse, Rat)	0%
NPY (Human/Rat)	0%
Pancreatic Polypeptide (Human, Rat)	0%
Human GIP (1-42)	0%

## Precision

### Intra- and Inter-Assay Variation

<b>Samples</b>	<b>Mean Active Ghrelin (pg/mL)</b>	<b>Intra-Assay %CV</b>	<b>Inter-Assay %CV</b>
Rat Serum 1	61	4.90	9.81
Rat Serum 2	206	1.08	2.10
Rat Serum 3	1,046	1.00	1.40
Mouse Serum 1	54	7.09	9.26
Mouse Serum 2	195	1.60	3.41
Mouse Serum 3	1,036	0.27	1.56
Rat Plasma 1	90	0.57	4.90
Rat Plasma 2	231	1.27	3.27
Rat Plasma 3	982	0.65	1.15
Mouse Plasma 1	58	5.56	4.70
Mouse Plasma 2	210	2.70	4.08
Mouse Plasma 3	1,035	0.86	1.23

Serum or plasma samples from rats and mice are pooled and treated with AEBSF and HCl, then divided into 3 aliquots each. Various amounts of rat/mouse ghrelin are added to the aliquots to create low, intermediate and high levels of ghrelin samples for precision tests. Intra-assay variations were calculated from results of six duplicate determinations in one assay. Inter-assay variations were calculated from results of six separate assays with duplicate samples in each assay.



## Spike Recovery of Rat/Mouse Ghrelin (Active) in Assay Samples

Sample I.D.	Active Ghrelin Spiked (pg/mL)	Serum Ghrelin		Plasma Ghrelin	
		pg/mL	Recovery Rate	pg/mL	Recovery Rate
Rat	0 (Basal)	0	-	53	-
	50	60	96%	106	106%
	200	200	94%	252	100%
	1,000	885	102%	1,019	97%
Rat	0 (Basal)	0	-	103	-
	50	36	72%	154	102%
	200	172	86%	285	91%
	1,000	1,070	90%	973	87%
Rat	0 (Basal)	168	--	181	-
	50	218	100%	233	104%
	200	356	94%	374	97%
	1,000	1,070	90%	1,100	92%
Mouse	0 (Basal)	92	-	53	-
	50	129	74%	109	112%
	200	269	89%	263	105%
	1,000	965	87%	1,073	102%
Mouse	0 (Basal)	67	-	116	-
	50	125	116%	174	116%
	200	296	115%	317	101%
	1,000	1,156	109%	1,079	96%
Mouse	0 (Basal)	136	-	213	-
	50	193	114%	259	92%
	200	359	112%	396	92%
	1,000	1,166	103%	1,112	90%
Mean S.D. (n = 6)	50	-	95.3 ±19%	-	105.3 ±8.4%
	200	-	98.3 ±12%	-	97.7 ±5.4%
	1,000	-	95.3 ±11%	-	94.0 ±5.4%

Varying amounts of rat/mouse ghrelin were added to individual serum and plasma samples from rat and mouse and the resulting ghrelin content of each sample was assayed by Rat/Mouse Ghrelin (active) ELISA. The recovery = [(observed ghrelin concentration after spike – Basal Ghrelin level) / spiked ghrelin concentration] X 100%.

## Linearity of Sample Dilution

Sample	Volume Assayed	Serum Ghrelin		Plasma Ghrelin	
		pg/mL	% of expected	pg/mL	% of Expected
Rat	20 µL	212	100%	221	100%
	15 µL	153	96%	159	96%
	10 µL	100	94%	110	100%
	5 µL	54	102%	57	103%
Rat	20 µL	640	100%	607	100%
	15 µL	460	96%	448	98%
	10 µL	304	95%	300	99%
	5 µL	159	99%	153	101%
Rat	20 µL	1288	100%	1220	100%
	15 µL	920	95%	912	100%
	10 µL	606	94%	611	100%
	5 µL	302	94%	309	101%
Mouse	20 µL	192	100%	219	100%
	15 µL	140	97%	159	97%
	10 µL	98	102%	104	95%
	5 µL	50	104%	45	82%
Mouse	20 µL	567	100%	621	100%
	15 µL	140	97%	159	97%
	10 µL	98	102%	104	95%
	5 µL	50	104%	45	83%
Mouse	20 µL	1,067	100%	1,297	100%
	15 µL	808	101%	961	99%
	10 µL	547	103%	635	98%
	5 µL	286	108%	323	100%
Mean S.D. (n = 6)	20 µL	-	100%	-	100%
	15 µL	-	97.2 ±2.1%	-	98.0 ±1.4%
	10 µL	-	98.2 ±4.3%	-	98.3 ±1.9%
	5 µL	-	102.3 ±5.2%	-	96.5 ±8.1%

Spiked and pooled samples were assayed at 20, 15, 10 and 5 µL each for active ghrelin by ELISA. Measured ghrelin levels are corrected for various dilution factors and then divided by levels found at 20 µL sample size to obtain the % of expected values.

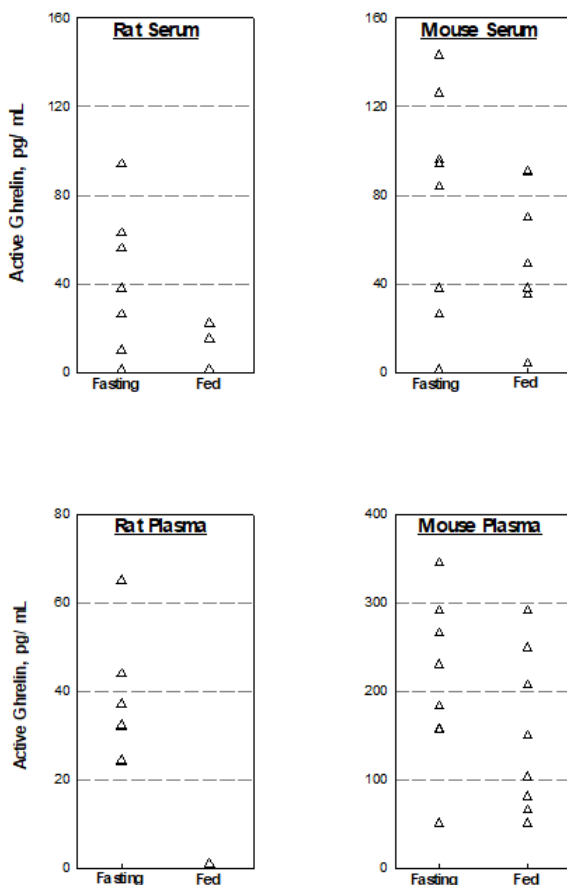
---

## Normal Range of Active Ghrelin Levels in Rat/Mouse Blood

The normal range of active ghrelin in 24-hour fasted rat (Sprague Dawley) and mouse (CD-1) blood is ~ 120 pg/mL and 600 pg/mL, respectively. The levels are lower in non-fasted animals.

# Post Prandial Attenuation of Active Ghrelin in Blood

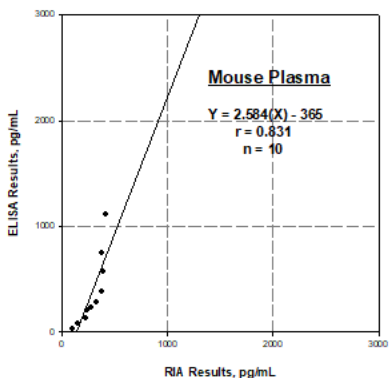
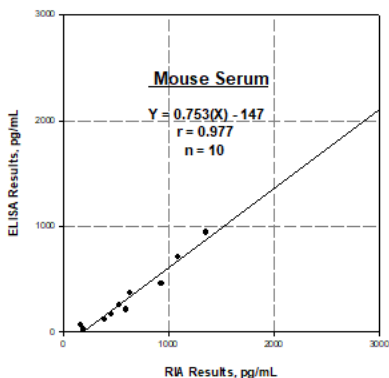
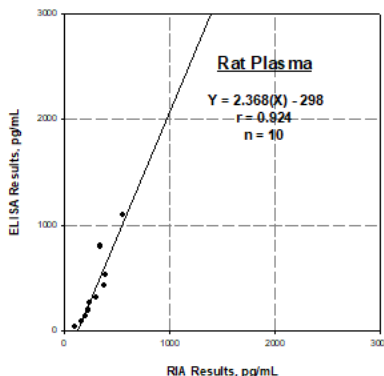
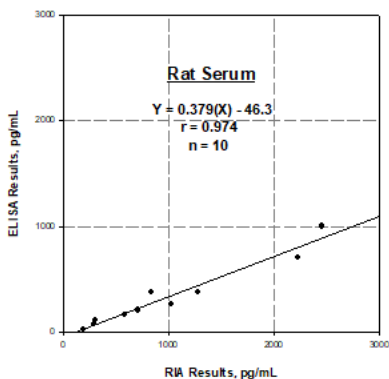
## Effects of Fasting on Serum/Plasma Ghrelin (Active) Levels



Each group contains 8 animals, either fed *ad lib* or 24-hour fasted before blood collection. All blood samples were treated immediately with 1 mg/mL AEBSF and processed for serum/plasma isolation. Resulting serum/plasma samples are acidified to 0.05 N HCl and stirred at -20 °C before ELISA assay.

## RIA Cat. No. GHRA-88HK vs. ELISA Cat. No. EZRGRA-90K

Pooled neat serum or plasma sample from rats/mice are thawed, treated with 1 mg/mL AESBF followed by acidification with 0.05 N HCl, spiked with rat/mouse ghrelin at levels from 50–1,000 pg/mL, and then assayed for active ghrelin by ELISA and RIA. Paired results are analyzed using linear regression.



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

---

## 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](https://www.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.
- High absorbance in background or blank wells could be due to
  - cross well contamination by standard solution or sample, or
  - inadequate washing of wells with Wash Buffer, or
  - overexposure to light after substrate has been added.

## Product Ordering

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

### Replacement Reagents

Reagents	Cat. No.
Microtiter Plate	EPDAR
10X HRP Wash Buffer Concentrate	EWB-HPR
Rat/Mouse Ghrelin (Active) Standard	E8090-K
Rat/Mouse Ghrelin (Active) Quality Controls 1 & 2	E6090-K
Matrix Solution	EMTX-GA
Assay Buffer	EABGR
Rat/Mouse Ghrelin (Active) Capture Antibody	E1090-C
Rat/Mouse Ghrelin (Active) Detection Antibody	E1090-D
Enzyme Solution	EHRP-88
Substrate	ESS-TMB2
Stop Solution	ET-TMB

---

## 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.

© 2021-2024 Merck KGaA, Darmstadt, Germany and/or its affiliates.  
All Rights Reserved.

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

IFU-EZRGRA-90K Rev 03/24

The Merck logo, consisting of the word "MERCK" in a bold, blue, sans-serif font.