

Rat Leptin ELISA Kit ab100773

★★★★★ 1 Abreviews 22 References 4 Images

Overview

Product name	Rat Leptin ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 30 pg/ml
Range	10.97 pg/ml - 8000 pg/ml
Recovery	84 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	75.43	68% - 89%
Serum	92.32	83% - 102%
Plasma	84.25	75% - 94%

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Rat

Product overview Abcam’s Leptin Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of rat Leptin in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for rat Leptin coated on a 96-well plate. Standards and samples are pipetted into the wells and Leptin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-rat Leptin antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Leptin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Microplate

Properties

**Storage instructions**

Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Biotinylated anti-Rat Leptin	2 vials
Leptin Microplate (12 x 8 wells)	1 unit
Recombinant rat Leptin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function**

May function as part of a signaling pathway that acts to regulate the size of the body fat depot. An increase in the level of LEP may act directly or indirectly on the CNS to inhibit food intake and/or regulate energy expenditure as part of a homeostatic mechanism to maintain constancy of the adipose mass.

**Involvement in disease**

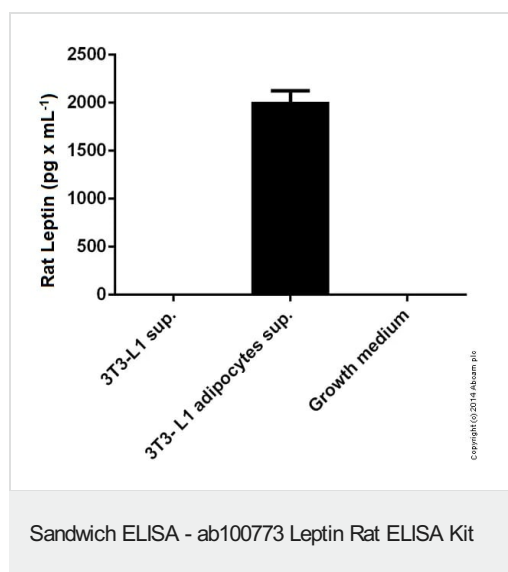
Defects in LEP may be a cause of obesity (OBESITY) [MIM:601665]. It is a condition characterized by an increase of body weight beyond the limitation of skeletal and physical requirements, as the result of excessive accumulation of body fat.

**Sequence similarities**

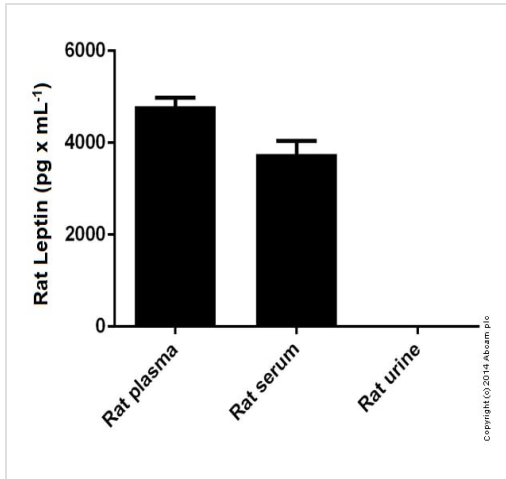
Belongs to the leptin family.

**Cellular localization**

Secreted.

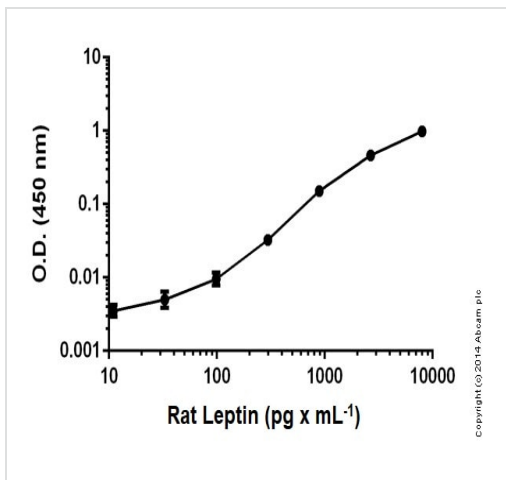
**Images**

Rat Leptin measured in supernatanses showing quantity (ng) per mL of tested sample



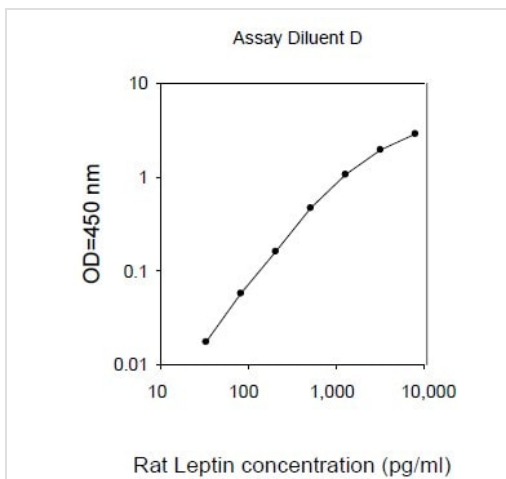
Rat Leptin measured in biofluids showing quantity (ng) per mL of tested sample

Sandwich ELISA - ab100773 Leptin Rat ELISA Kit



Standard curve: mean of duplicates (+/-SD) with background readings subtracted

Sandwich ELISA - ab100773 Leptin Rat ELISA Kit



ELISA - Leptin Rat ELISA Kit (ab100773)

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### **Our Abpromise to you: Quality guaranteed and expert technical support**

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
  
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors