

Mouse Leptin ELISA Kit ab100718

★★★★☆ 2 Abreviews 29 References 4 Images

Overview

Product name	Mouse Leptin ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 4 pg/ml
Range	4.1 pg/ml - 1000 pg/ml
Recovery	94 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	94.47	83% - 103%
Serum	93.29	82% - 102%
Plasma	95.38	84% - 104%

Assay duration	Multiple steps standard assay
Species reactivity	<b>Reacts with:</b> Mouse
Product overview	Mouse Leptin ELISA kit is designed for the quantitative measurement of Mouse Leptin in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Mouse 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-Mouse 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
120X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse Leptin	2 vials
Leptin Microplate (12 strips x 8 wells)	1 unit
Recombinant Mouse 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.

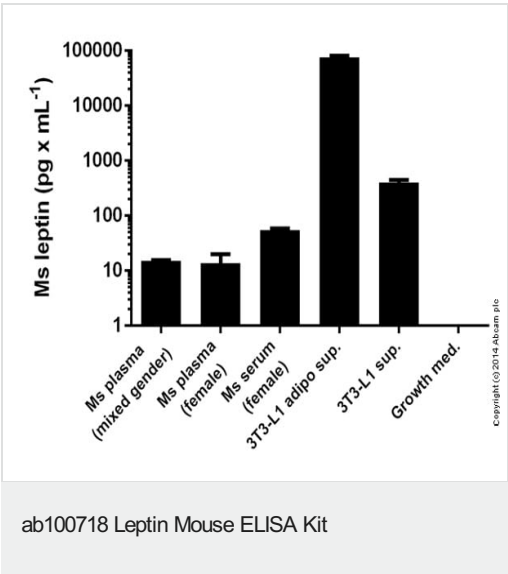
Sequence similarities

Belongs to the leptin family.

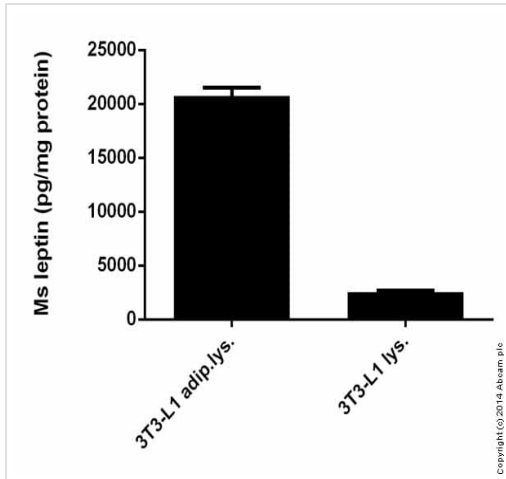
Cellular localization

Secreted.

Images

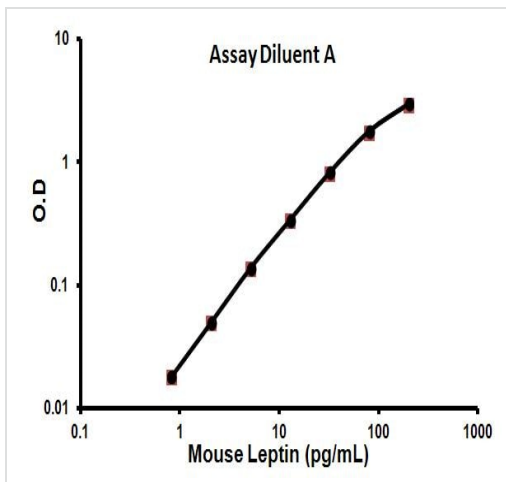


Ms Leptin measured in biological fluids showing quantity (pg) per mL of tested sample



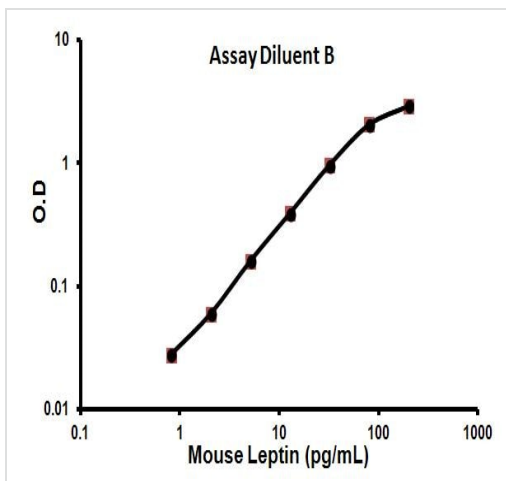
Ms Leptin measured in cell lysates showing quantity (pg) per mg protein

ab100718 Leptin Mouse ELISA Kit



Representative Standard Curve using ab100718

Typical Standard Curve



Representative Standard Curve using ab100718

Typical Standard Curve

**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