

Mouse Leptin ELISA Kit ab100718

★★★★☆ [2 Abreviews](#) [29 References](#) [4 图像](#)

概述

产品名称	小鼠Leptin ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 4 pg/ml
范围	4.1 pg/ml - 1000 pg/ml
回收率	94 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	94.47	83% - 103%
Serum	93.29	82% - 102%
Plasma	95.38	84% - 104%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Mouse
产品概述	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.

平台	Microplate
----	------------

性能

存放说明

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

组件	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

功能

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.

疾病相关

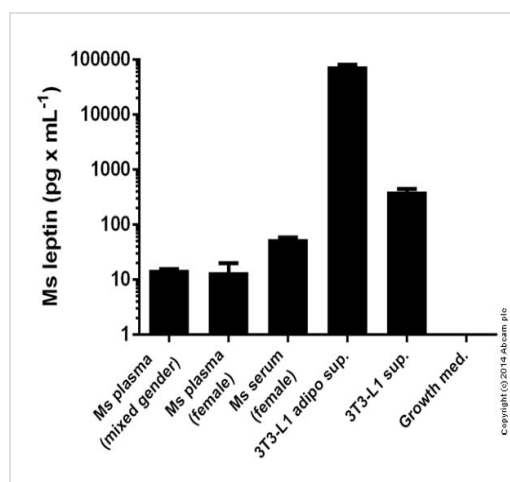
Defects in LEP may be a cause of obesity (OBESITY) [MIM:601665]. It is a condition characterized by an increase of body weight beyond the limitation of skeletal and physical requirements, as the result of excessive accumulation of body fat.

序列相似性

Belongs to the leptin family.

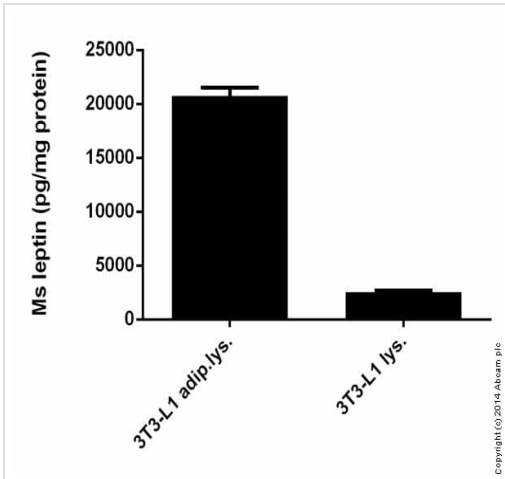
细胞定位

Secreted.

图片

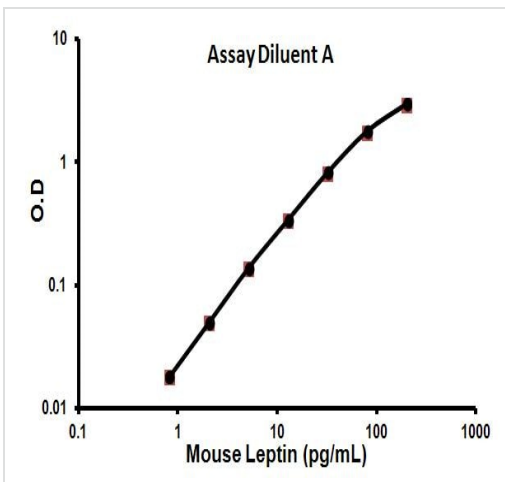
Ms Leptin measured in biological fluids showing quantity (pg) per mL of tested sample

ab100718 Leptin Mouse ELISA Kit



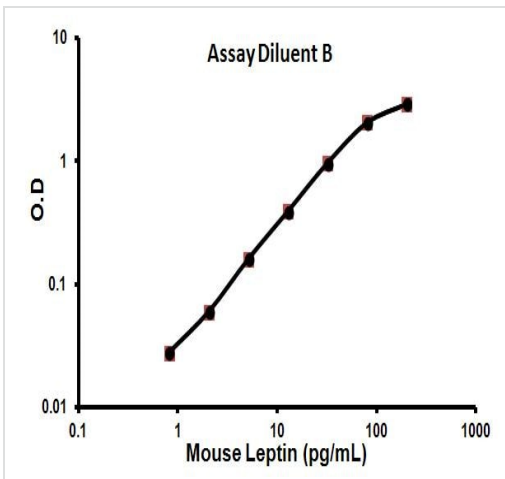
Ms Leptin measured in cell lysates showing quantity (pg) per mg protein

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.cn/abpromise> or contact our technical team.

Terms and conditions

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