

Mouse Leptin ELISA Kit ab108881

1 References 1 图像

概述

产品名称

小鼠Leptin ELISA试剂盒

检测方法

Colorimetric

精确度

批次内

| 样品      | n | Mean | SD | CV%  |
|---------|---|------|----|------|
| Overall |   |      |    | 4.5% |

批次间

| 样品      | n | Mean | SD | CV%  |
|---------|---|------|----|------|
| Overall |   |      |    | 9.1% |

样品类型

Cell culture supernatant, Serum, Plasma

检测类型

Sandwich (quantitative)

灵敏度

= 0.3 ng/ml

范围

0.375 ng/ml - 24 ng/ml

检测时间

5h 00m

实验步骤

Multiple steps standard assay

种属反应性

与反应: Mouse

产品概述

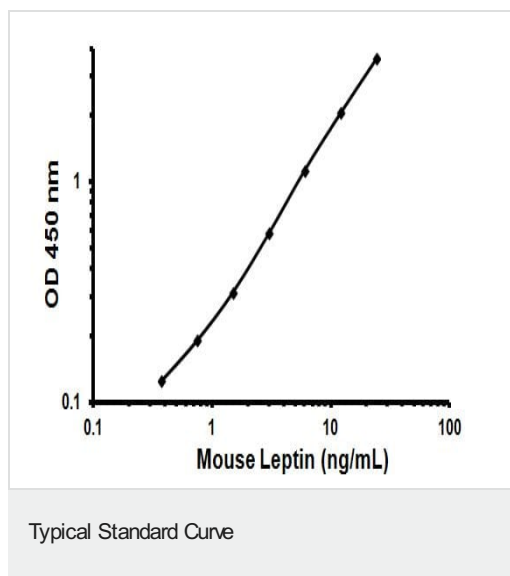
Abcam's Leptin mouse *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of leptin levels in plasma, serum and cell culture supernatants.

A Leptin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Leptin specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Leptin captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

|  |   |
|--|---|
| 平台                                     | Microplate  |
| 性能                                     |   |
| 存放说明                                   | Store at -20°C. Please refer to protocols.  |
| 组件                                     | 1 x 96 tests  |
| 100X Streptavidin-Peroxidase Conjugate | 1 x 80µl  |
| 10X Diluent N Concentrate              | 1 x 30ml  |
| 1X Standard Diluent                    | 1 x 2ml   |
| 20X Wash Buffer Concentrate            | 2 x 30ml  |
| 70X Biotinylated Mouse Leptin Antibody | 1 x 90µl  |
| Chromogen Substrate                    | 1 x 7ml   |
| Leptin Microplate (12 x 8 well strips) | 1 unit  |
| Leptin Standard                        | 1 x 2 vials   |
| Sealing Tapes                          | 3 units   |
| Stop Solution                          | 1 x 11ml  |
| 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



Representative Standard Curve using ab108881

**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