

# Leptin (rat), ELISA kit

The Leptin (rat), ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The Leptin (rat), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 3 hours.

The adipokine leptin is a hormone secreted predominantly by adipose tissue that signals through leptin receptors in the hypothalamus to decrease appetite and increase energy expenditure. Binding of leptin to the long-form of the leptin receptor in the hypothalamus reduces neuropeptide Y (NPY) and agouti-related protein (AgRP) activity, while stimulating anorexigenic pro-opiomelanocortin (POMC) neuron activity to reduce appetite. In peripheral tissues, leptin antagonizes insulin signaling, increases fatty acid oxidation, decreases insulin production in pancreatic beta cells, and promotes fertility. Its expression in adipocytes can be regulated in a paracrine fashion by other adipokines such as IL-6 (stimulation) and TNF- $\alpha$  (inhibition). Leptin deficient mice display severe insulin resistance, obesity, and decreased fertility, all of which are reversible by administration of exogenous leptin. In contrast, elevated circulating levels of leptin are associated with increased obesity, indicative of an acclimated state of leptin resistance which is not well understood.

Citations: 21

[View Online »](#)

## Ordering Information

[Order Online »](#)

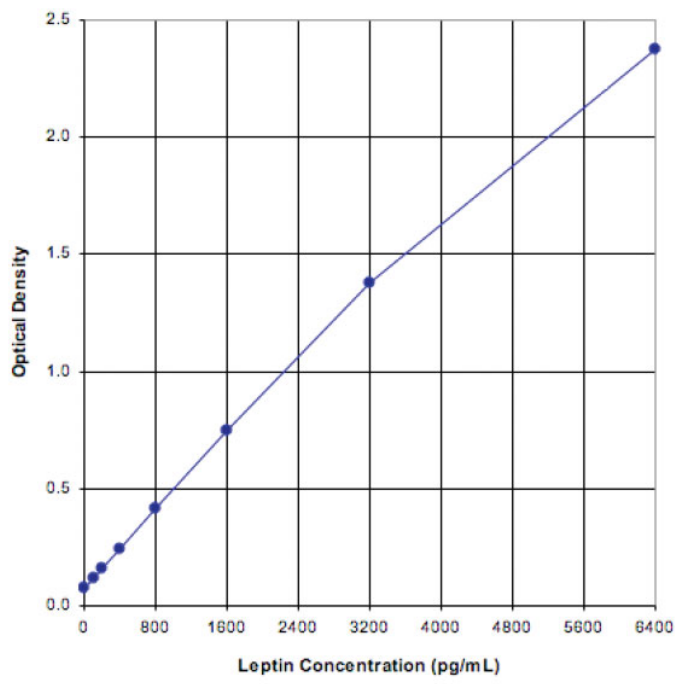
ADI-900-015A

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- No edge effect - all incubations at room temperature eliminate risk of evaporation at 37°C
- Broad dynamic range - easily quantify high and low levels of leptin
- Convenient- store all reagents at 4°C
- Fast - fully quantitative results in just 3 hours from up to 39 samples in duplicate
- Pre-coated plates and ready-to-use reagents save you time



## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name OB gene product

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of rat leptin in culture supernatants, plasma, and serum.

Assay Time 3 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter plate, Standard, Antibody, Conjugate, Assay buffer 28, TMB Substrate, Stop solution 2, Wash buffer concentrate

Sensitivity 67.2 pg/ml (range 100 – 6,400 pg/ml)

Species Reactivity Rat

Technical Info / Product Notes **Cited samples:**  
[Cytokine ELISAs Cited Samples](#)

UniProt ID P50596

Wavelength 450 nm



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)