

Leptin (human), ELISA kit

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

The Leptin ELISA 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- α (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: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

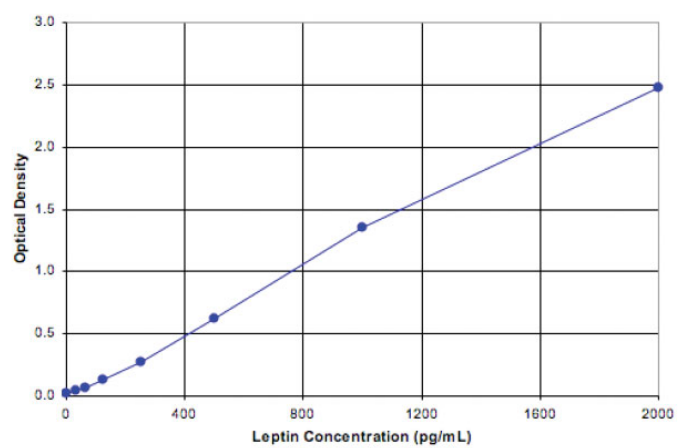
ADI-900-028A

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Sensitive - accurately quantify as little as 23.4 pg/mL
- Save time - obtain results from 39 samples in duplicate in just 3 hours!
- Convenient- store all reagents at 4°C
- 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 human 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 17, TMB Substrate, Stop solution 2, Wash buffer concentrate

Sensitivity 23.4 pg/ml (range 31.3 – 2,000 pg/ml)

Species Reactivity Human

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

UniProt ID P41159

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)