

Leptin (mouse), (recombinant)

Leptin, the product of the *ob* (obese) gene, is a 16kDa protein consisting of 146 amino acid residues. Leptin is produced in the adipose tissue, and is considered to play an important role in appetite control, fat metabolism and regulation of body weight. Studies have shown that it may also influence reproductive function. Mice with *ob/ob* genotype develop a form of diabetes similar to type II in humans.

Ordering Information

[Order Online »](#)

ALX-201-035-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Reconstituted protein is stable for 4 weeks when stored at +4°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	OB gene product
Biological Activity	Induces proliferation of BAF/3 cells stably transfected with the long form of human leptin receptor.
CAS	181030-10-4
Endotoxin Content	<0.1EU/μg purified protein.
Formulation	Lyophilized from 0.0045mM sodium bicarbonate.
MI	14: 5443
MW	~16kDa.
Purity	≥98% (SDS-PAGE)
Reconstitution	Reconstitute with sterile water or 0.4% sodium bicarbonate, pH 8-9. Do not reconstitute to less than 0.1mg/ml. Further dilutions should be made with medium containing a carrier protein.
Source	Produced in <i>E. coli</i> . Homologous to mouse serum leptin.
UniProt ID	P41160



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-
fr@enzolifesciences.com](mailto:info-fr@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)