

# Leptin (human), (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. It targets the central nervous system, particularly hypothalamus, affecting food intake. Leptin levels are high in most obese individuals. Studies have shown that it may also influence reproductive function.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-201-034-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 freeze in liquid nitrogen.
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	177404-21-6
Endotoxin Content	<0.1EU/μg purified protein.
Formulation	Lyophilized from 0.0045mM sodium bicarbonate.
MI	14: 5443
MW	~16kDa.
Purity	≥95% (SDS-PAGE; reducing and non-reducing)
Purity Detail	The purified recombinant human leptin consists of >93% monomers as determined by gel-filtration chromatography.
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 0.1% HSA or BSA.

**Source** Produced in *E. coli*. Homologous to human serum leptin.



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

P41159

European Sales Office  
ENZO LIFE SCIENCES  
(UK) 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)