

# Obestatin (mouse, rat)

Food intake modulator

Obestatin derives like ghrelin from proghrelin, but exhibits opposite physiological effects. Obestatin decreases food intake and body weight in rodents. Moreover, obestatin also counteracts ghrelin's effect by decelerating gastric emptying and decreasing intestinal contractility. Obestatin binds to orphan G protein-coupled receptor GPR39. Amidation of the C-terminus seems to be important for obestatin's activity.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-164-005-M001	1mg
------------------	-----

**Manuals, SDS & CofA**

**[View Online »](#)**

# Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

# Product Details

Formula	C <sub>114</sub> H <sub>174</sub> N <sub>34</sub> O <sub>31</sub>
Formulation	Lyophilized.
MW	2516.8
Peptide Content	≥80%
Purity	≥97% (HPLC)
Sequence	H-Phe-Asn-Ala-Pro-Phe-Asp-Val-Gly-Ile-Lys-Leu-Ser-Gly-Ala-Gln-Tyr-Gln-Gln-His-Gly-Arg-Ala-Leu-NH <sub>2</sub>
Solubility	Soluble in DMSO.

Last modified: May 29, 2024

Image not found or type unknown  
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)