

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 $C_{114}H_{174}N_{34}O_{31}$

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-NH₂

Solubility Soluble in DMSO.

Last modified: May 29, 2024



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)