

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 »](#)

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-NH ₂
Solubility	Soluble in DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com