

Ghrelin (human)

Endogenous GHS-R ligand

Ghrelin, a 28 aa peptide predominantly produced by the stomach, is an endogenous ligand for the growth hormone secretagogue receptor (GHS-R) that potently stimulates growth hormone release. Studies suggest that peripherally or centrally administered ghrelin, independent of growth hormone, decreases fat oxidation and increases food intake and adiposity. In addition, plasma ghrelin levels are lower in obese human subjects. Ghrelin promotes the production of orexigenic neuropeptides (neuropeptide Y and agouti-related protein) in the hypothalamic arcuate nuclei and activates the neurons which produce these orexigenic neuropeptides, resulting in an increase in feeding and body weight and thereby opposing the effects of leptin. Moreover, ghrelin shows regulatory effects on the cardiovascular and immune system.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-157-021-MC05	0.5mg
ALX-157-021-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

CAS 258279-04-8

Formula $C_{149}H_{249}N_{47}O_{42}$

Formulation Lyophilized.

MW 3370.9

Purity ≥98%

Sequence H-Gly-Ser-Ser(octanoyl)-Phe-Leu-Ser-Pro-Glu-His-Gln-Arg-Val-Gln-Gln-Arg-Lys-Glu-Ser-Lys-Lys-Pro-Pro-Ala-Lys-Leu-Gln-Pro-Arg-OH

Solubility Soluble in distilled water or 5% acetic acid.



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)