

# 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

## 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.

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-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)