

PRODUCT INFORMATION



Ghrelin (human) (trifluoroacetate salt)

Item No. 15072

Formal Name: glycy-L-seryl-O-(1-oxooctyl)-L-seryl-L-phenylalanyl-L-leucyl-L-seryl-L-prolyl-L- α -glutamyl-L-histidyl-L-glutamyl-L-arginyl-L-valyl-L-glutamyl-L-glutamyl-L-arginyl-L-lysyl-L- α -glutamyl-L-seryl-L-lysyl-L-lysyl-L-prolyl-L-prolyl-L-alanyl-L-lysyl-L-leucyl-L-glutamyl-L-prolyl-L-arginine, trifluoroacetate salt



Synonym: Lenomorelin

MF: $\text{C}_{149}\text{H}_{249}\text{N}_{47}\text{O}_{42} \cdot \text{XCF}_3\text{COOH}$

FW: 3,370.9

$\cdot \text{XCF}_3\text{COOH}$

Purity: $\geq 98\%$

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥ 4 years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ghrelin (human) (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the ghrelin (human) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Ghrelin, an endogenous ligand for the growth hormone secretagogue receptor, is synthesized principally in the stomach.¹ It stimulates food intake and transduces signals to hypothalamic regulatory nuclei that control energy homeostasis.² It is a peptide of 28 amino acids, in which the serine 3 residue is n-octanoylated, which is necessary for biological activity.^{1,3}

References

1. Kojima, M., Hosoda, H., Date, Y., *et al.* Ghrelin is a growth-hormone-releasing acylated peptide from stomach. *Nature* **402(6762)**, 656-660 (1999).
2. Murphy, K.G. and Bloom, S.R. Gut hormones in the control of appetite. *Exp. Physiol.* **89(5)**, 507-516 (2004).
3. Gutierrez, J.A., Solenberg, P.J., Perkins, D.R., *et al.* Ghrelin octanoylation mediated by an orphan lipid transferase. *Proc. Nat. Acad. Sci. USA* **105(17)**, 6320-6325 (2008).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/31/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM