

PRODUCT INFORMATION



Somatorelin (1-44) amide (human) (trifluoroacetate salt)

Item No. 33942

Formal Name: somatoliberin (human pancreatic islet),
trifluoroacetate salt

Synonyms: Growth Hormone-releasing Factor (1-44) amide,
hpGRF-44, Human Pancreas GRF-44,
Somatoliberin

MF: $C_{215}H_{358}N_{72}O_{66}S \cdot XCF_3COOH$

FW: 5,039.7

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Somatorelin (1-44) amide (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the somatorelin (1-44) amide (human) (trifluoroacetate salt) in water. The solubility of somatorelin (1-44) amide (human) (trifluoroacetate salt) in water is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Somatorelin (1-44) amide is a synthetic peptide that corresponds to the full 44-amino acid sequence of growth hormone-releasing hormone (GHRH), a hormone that stimulates the production of growth hormone from the pituitary gland and is involved in the regulation of cell proliferation and differentiation.^{1,2} Somatorelin induces growth hormone release similarly to GHRH *in vitro* and *in vivo*.¹

References

1. Guillemin, R., Brazeau, P., Böhlen, P., *et al.* Growth hormone-releasing factor from a human pancreatic tumor that caused acromegaly. *Science* **218(4572)**, 585-587 (1982).
2. Kiaris, H., Chatzistamou, I., Papavassiliou, A.G., *et al.* Growth hormone-releasing hormone: Not only a neurohormone. *Trends Endocrinol. Metab.* **22(8)**, 311-317 (2011).

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, 11/21/2022

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