

PRODUCT INFORMATION



Alexamorelin (trifluoroacetate salt)

Item No. 36586

Formal Name: L-alanyl-L-histidyl-2-methyl-D-tryptophyl-L-alanyl-L-tryptophyl-D-phenylalanyl-L-lysineamide, trifluoroacetate salt

Synonym: Ala-His-D-2-Methyl-Trp-Ala-Trp-D-Phe-Lys-NH₂

MF: C₅₀H₆₃N₁₃O₇ • XCF₃COOH

FW: 958.1

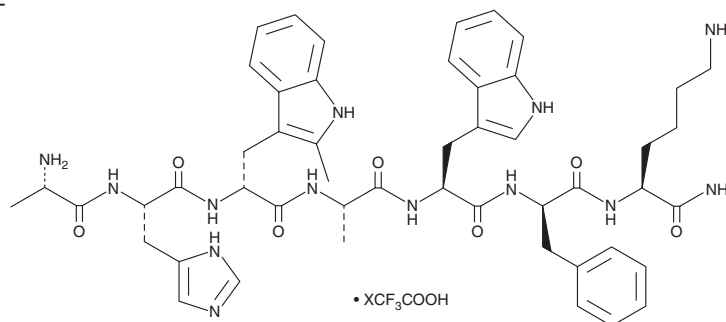
Purity: ≥98%

UV/Vis.: λ_{max}: 222 nm

Supplied as: A 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

Alexamorelin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the alexamorelin (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Alexamorelin (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

Alexamorelin is a synthetic heptapeptide agonist of the growth hormone secretagogue receptor (GHS-R).¹

Reference

1. Broglio, F., Benso, A., Gottero, C., *et al.* Endocrine activities of alexamorelin (Ala-His-D-2-methyl-Trp-Ala-Trp-D-Phe-Lys-NH₂), a synthetic GH secretagogue, in humans. *Eur. J. Endocrinol.* **143(3)**, 419-425 (2000).

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, 05/08/2023

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