

PRODUCT INFORMATION



Alytesin (trifluoroacetate salt)

Item No. 34428

Formal Name: 5-oxo-L-prolylglycyl-L-arginyl-L-leucylglycyl-L-threonyl-L-glutamyl-L-tryptophyl-L-alanyl-L-valylglycyl-L-histidyl-L-leucyl-L-methioninamide, trifluoroacetate salt

MF: C₆₈H₁₀₆N₂₂O₁₇S • XCF₃COOH

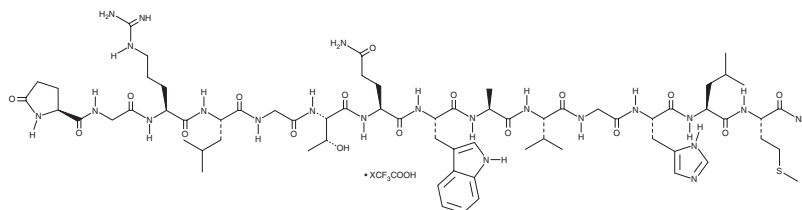
FW: 1,535.8

Purity: ≥98%

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

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

Description

Alytesin is a neuropeptide originally isolated from frog (*A. obstetricans*) skin that has diverse biological activities.¹⁻³ It induces hypertension with tachyphylaxis and increases blood glucose levels in dogs.¹ Alytesin reduces gastric acid secretion in rats.² It also induces anorexigenic effects in neonatal chicks.³

References

1. Erspamer, V., Erpamer, G.F., and Inselvini, M. Some pharmacological actions of alytesin and bombesin. *J. Pharm. Pharmacol.* **22(11)**, 875-876 (1970).
2. Guglietta, A., Strunk, C.L., Irons, B.J., et al. Central neuromodulation of gastric acid secretion by bombesin-like peptides. *Peptides* **6 Suppl 3**, 75-81 (1985).
3. Cline, M.A., Fouse, D.N., and Prall, B.C. Central and peripheral alytesin cause short-term anorexigenic effects in neonatal chicks. *Neuropeptides* **42(3)**, 283-291 (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, 12/06/2022

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