

PRODUCT INFORMATION



Allatostatin I (trifluoroacetate salt)

Item No. 43824

Formal Name: L-alanyl-L-prolyl-L-serylglycyl-L-alanyl-L-glutaminy-L-arginyl-L-leucyl-L-tyrosylglycyl-L-phenylalanylglycyl-L-leucinamide, trifluoroacetate salt

Peptide Sequence: APSGAQRLYGFGGL-NH₂

MF: C₆₁H₉₄N₁₈O₁₆ • XCF₃COOH

FW: 1,335.5

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

H-Ala-Pro-Ser-Gly-Ala-Gln-Arg-Leu-Tyr-Gly-Phe-Gly-Leu-NH₂ • XCF₃COOH

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

Laboratory Procedures

Allatostatin I (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the allatostatin I (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Allatostatin I (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO and a 1:1 solution of acetonitrile:water.

Description

Allatostatin I is an insect neuropeptide.¹ It is found primarily in the insect corpora allata, which are endocrine glands involved in juvenile hormone biosynthesis. Allatostatin I inhibits juvenile hormone biosynthesis in female cockroaches (*D. punctata*) at a variety of life stages.² It reduces hindgut motility, but not juvenile hormone production, in 2-day-old ring-legged earwigs (*E. annulipes*).¹

References

1. Rankin, S.M., Stay, B., Chan, K., *et al.* Cockroach allatostatin-immunoreactive neurons and effects of cockroach allatostatin in earwigs. *Arch. Insect Biochem. Physiol.* **38(4)**, 155-165 (1998).
2. Pratt, G.E., Farnsworth, D.E., and Feyerisen, R. Changes in the sensitivity of adult cockroach corpora allata to a brain allatostatin. *Mol. Cell. Endocrinol.* **70(2)**, 185-195 (1990).

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, 09/15/2025

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