

PRODUCT INFORMATION



Myomodulin A (*Aplysia*, *Lymnaea*) (trifluoroacetate salt)

Item No. 40151

Formal Name: L-prolyl-L-methionyl-L-seryl-L-methionyl-L-leucyl-L-arginyl-L-leucinamide, trifluoroacetate salt

Synonyms: MMA, Pro-Met-Ser-Met-Leu-Arg-Leu-NH₂

Peptide Sequence: PMSMLRL-NH₂

MF: C₃₆H₆₇N₁₁O₈S₂ • XCF₃COOH

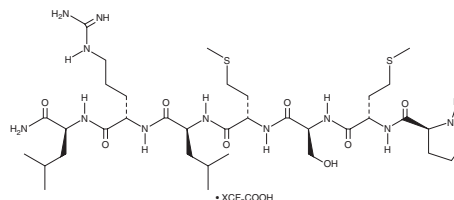
FW: 846.1

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

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

Description

Myomodulin A is a gastropod neuropeptide.¹ It is found in gastropod ventral lobe neurons, neural elements of the accessory radula closer (ARC) muscle, and penis nerve.¹⁻³ Myomodulin A is biosynthesized by ARC motor neurons in *A. californica* and potentiates motor neuron-elicited contractions of the *A. californica* buccal mass ARC.² Myomodulin A also increases the relaxation rate of electrically evoked muscle contractions in *L. stagnalis* penis retractor muscle.³

References

1. Miller, M.W., Beushausen, S., Vitek, A., *et al.* The myomodulin-related neuropeptides: Characterization of a gene encoding a family of peptide cotransmitters in *Aplysia*. *J. Neurosci.* **13**(8), 3358-3367 (1993).
2. Cropper, E.C., Tenenbaum, R., Kolks, M.A., *et al.* Myomodulin: a bioactive neuropeptide present in an identified cholinergic buccal motor neuron of *Aplysia*. *Proc. Natl. Acad. Sci. USA* **84**(15), 5483-5486 (1987).
3. Li, K.W., van Golen, F.A., van Minnen, J., *et al.* Structural identification, neuronal synthesis, and role in male copulation of myomodulin-A of *Lymnaea*: A study involving direct peptide profiling of nervous tissue by mass spectrometry. *Brain Res. Mol. Brain Res.* **25**(3-4), 355-358 (1994).

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, 02/29/2024

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