

PRODUCT INFORMATION



Buccalin A (trifluoroacetate salt)

Item No. 39728

Formal Name: (3S,6S,9S,12S,15S,18S,27S)-3-((S)-2-(2-aminoacetamido)-4-(methylthio)butanamido)-15-benzyl-27-carbamoyl-6,18-bis(hydroxymethyl)-9-isobutyl-12,29-dimethyl-4,7,10,13,16,19,22,25-octaoxo-5,8,11,14,17,20,23,26-octazatriacontanoic acid, trifluoroacetate salt

Peptide Sequence: H-Gly-Met-Asp-Ser-Leu-Ala-Phe-Ser-Gly-Gly-Leu-NH₂

MF: C₄₅H₇₂N₁₂O₁₅S • XCF₃COOH

FW: 1,053.2

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

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

Description

Buccalin A is a neuropeptide that is found in gastropod motor neurons and is involved in neuromuscular signaling.¹ It reduces current-induced muscle contraction amplitude but not muscle relaxation rate in isolated and washed *A. californica* accessory radula closer (ARC) muscle fibers when used at a concentration of 10 nM. Buccalin A (10 μM) decreases the amplitude of inhibitory post-synaptic currents (IPSCs) in isolated and perfused voltage-clamped *A. californica* buccal ganglia.²

References

1. Cropper, E.C., Miller, M.W., Tenenbaum, R., *et al.* Structure and action of buccalin: A modulatory neuropeptide localized to an identified small cardioactive peptide-containing cholinergic motor neuron of *Aplysia californica*. *Proc. Natl. Acad. Sci. USA* **85(16)**, 6177-6181 (1988).
2. Baux, G., Fossier, P., Trudeau, L.E., *et al.* Transmitter release and calcium currents at an *Aplysia* buccal ganglion synapse--II. Modulation by presynaptic receptors. *Neuroscience* **53(2)**, 581-593 (1993).

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, 02/13/2024

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