

PRODUCT INFORMATION

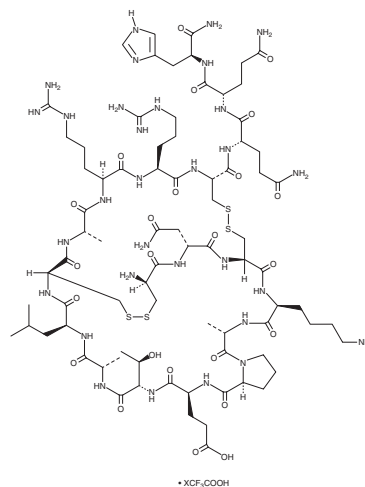


Apamin (trifluoroacetate salt)

Item No. 17082

CAS Registry No.: 24345-16-2
Formal Name: L-cysteinyl-L-asparaginyl-L-cysteinyl-L-lysyl-L-alanyl-L-prolyl-L- α -glutamyl-L-threonyl-L-alanyl-L-leucyl-L-cysteinyl-L-alanyl-L-arginyl-L-arginyl-L-cysteinyl-L-glutamyl-L-glutamyl-L-histidinamide, cyclic (1 \rightarrow 11),(3 \rightarrow 15)-bis(disulfide), trifluoroacetate salt

Synonym: Ro 23-6721
MF: C₇₉H₁₃₁N₃₁O₂₄S₄ • XCF₃COOH
FW: 2,027.3
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Apamin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the apamin (trifluoroacetate salt) in water. The solubility of apamin (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Apamin is an 18-amino acid peptide toxin that has been found in honey bee (*A. mellifera*) venom.^{1,5} It is an inhibitor of small conductance calcium-activated potassium (K_{Ca}2.2/SK2) channels (IC₅₀s = 0.06-0.4 nM).¹ It is selective for K_{Ca}2.2/SK2 channels over K_{Ca}2.1/SK1 and K_{Ca}2.3/SK3 channels (IC₅₀s = 1-12 and 1-13 nM, respectively), as well as the intermediate-conductance calcium-activated potassium channel K_{Ca}3.1/SK4 at 1 μ M but does not inhibit voltage-gated potassium channel 1.3 (K_v1.3; IC₅₀ = 13 nM).^{1,2,6} Apamin (50 μ g/kg) reduces CD8⁺ cell infiltration into the CNS and the severity of experimental autoimmune encephalomyelitis (EAE), as well as increased peripheral blood levels of CD62L⁻, CD44L⁻, and CD44⁺ memory T cells in mice.⁷ It has commonly been used to determine the role of K_{Ca}/SK channels in the brain, vascular endothelium, and bladder.^{3-5,8}

References

1. Wulff, H. and Köhler, R. *J. Cardiovasc. Pharmacol.* **61(2)**, 102-112 (2013).
2. Maylie, J., Bond, C.T., Herson, P.S., et al. *J. Physiol.* **554(Pt 2)**, 255-261 (2004).
3. Triggle, C.R., Hollenberg, M., Anderson, T.J., et al. *J. Smooth Muscle Res.* **39(6)**, 249-267 (2003).
4. Pedarzani, P. and Stocker, M. *Cell. Mol. Life Sci.* **65(20)**, 3196-3217 (2008).
5. Gu, H., Han, S.M., and Park, K.-K. *Toxins (Basel)* **12(3)**, 195 (2020).
6. Voos, P., Yazar, M., Lautenschläger, R., et al. *Eur. Biophys. J.* **46(6)**, 517-523 (2017).
7. Fattahi, H., Esmaeil, N., and Aliomrani, M. *Neurotox. Res.* **39(6)**, 1881-1891 (2021).
8. Thorneloe, K.S., Knorn, A.M., Doetsch, P.E., et al. *Am. J. Physiol. Regul. Integr. Comp. Physiol.* **294(5)**, R1737-R1743 (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, 11/17/2022

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