

PRODUCT INFORMATION



Kassinin (frog) (trifluoroacetate salt)

Item No. 39727

Peptide Sequence: DVPKSDQFVGLM-NH₂ H-Asp-Val-Pro-Lys-Ser-Asp-Gln-Phe-Val-Gly-
MF: C₅₉H₉₅N₁₅O₁₈S • XCF₃COOH Leu-Met-NH₂
FW: 1,334.6 • XCF₃COOH
Purity: ≥95%
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

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

Description

Kassinin is a tachykinin neuropeptide and agonist of the neurokinin-2 (NK₂) and NK₃ receptors.¹⁻⁴ It is selective for the NK₂ and NK₃ receptors over the NK₁ receptor.⁴ Kassinin (1 μM) increases the short-circuit current in isolated frog skin.⁵ It induces contractions in isolated guinea pig ileum and isolated rabbit jejunum in a concentration-dependent manner.³ Kassinin decreases blood pressure in anesthetized guinea pigs (ED₅₀ = 1,832 pmol/kg).²

References

1. Anastasi, A., Montecucchi, P., Erspamer, V., *et al.* Amino acid composition and sequence of kassinin, a tachykinin dodecapeptide from the skin of the African frog *Kassina senegalensis*. *Experientia* **33(7)**, 857-858 (1977).
2. Hancock, J.C. and Hoover, D.B. Effect of substance P and other tachykinins on arterial pressure in guinea-pigs. *J. Auton. Pharmacol.* **5(1)**, 25-30 (1985).
3. Holzer-Petsche, U., Schimek, E., Amann, R., *et al.* *In vivo* and *in vitro* actions of mammalian tachykinins. *Naunyn Schmiedebergs Arch. Pharmacol.* **330(2)**, 130-135 (1985).
4. Regoli, D., Drapeau, G., Dion, S., *et al.* New selective agonists for neurokinin receptors: pharmacological tools for receptor characterization. *Trends Pharmacol. Sci.* **9(8)**, 290-295 (1988).
5. Lippe, C., Bellantuono, V., Ardizzone, C., *et al.* Eledoisin and Kassinin, but not Enterokassinin, stimulate ion transport in frog skin. *Peptides* **25(11)**, 1971-1975 (2004).

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, 02/20/2024

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