

PRODUCT INFORMATION



Neurokinin A (trifluoroacetate salt)

Item No. 24221

Formal Name: L-histidyl-L-lysyl-L-threonyl-L- α -aspartyl-L-seryl-L-phenylalanyl-L-valylglycyl-L-leucyl-L-methioninamide, trifluoroacetate salt

Synonyms: Neurokinin α , Neuromedin L, Substance K

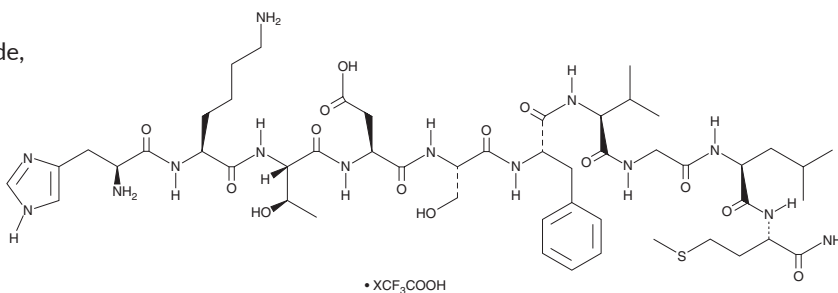
MF: C₅₀H₈₀N₁₄O₁₄S • XCF₃COOH
FW: 1,133.3

Purity: \geq 95%

Supplied as: A lyophilized powder

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

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

Description

Neurokinin A is a peptide agonist of the neurokinin-2 (NK₂) receptor.¹⁻³ It potentiates electrically-induced contractions of isolated rat vas deferens and increases vasodilation in the rat hind paw.² Neurokinin A (0.3-30 nmol/kg, i.v.) induces bronchoconstriction in guinea pigs in a dose-dependent manner (EC₅₀ = 29 nM), an effect which can be reversed by the NK₂ receptor antagonist MEN 10,207.^{1,3}

References

1. Maggi, C.A., Giuliani, S., Ballati, L., *et al.* *In vivo* evidence for tachykininergic transmission using a new NK-2 receptor-selective antagonist, MEN 10,376. *J. Pharmacol. Exp. Ther.* **257**(3), 1172-1178 (1991).
2. 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).
3. Laneuville, O., Couture, R., and Pace-Asciak, C.R. Neurokinin A-induced contraction of guinea-pig isolated trachea: Potentiation by hepoxilins. *Br. J. Pharmacol.* **107**(3), 808-812 (1992).

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, 11/16/2022

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