

PRODUCT INFORMATION



Spantide II (trifluoroacetate salt)

Item No. 43054

Formal Name: N⁶-(3-pyridinylcarbonyl)-D-lysyl-L-prolyl-3-(3-pyridinyl)-L-alanyl-L-prolyl-3,4-dichloro-D-phenylalanyl-L-asparaginyl-D-tryptophyl-L-phenylalanyl-D-tryptophyl-L-leucyl-L-norleucinamide, trifluoroacetate salt
Synonym: [D-NicLys¹,3-Pal³,D-Cl₂Phe⁵,Asn⁶,D-Trp^{7,9},Nle¹¹]-Substance P

Peptide Sequence: x₁PX₂PX₃NwFwLX₄-NH₂ (X₁ = N⁶-(nicotinylyl)-D-lysine; X₂ = 3-pyridinyl-L-alanine; X₃ = 3,4-dichloro-D-phenylalanine; X₄ = L-norleucine)

MF: C₈₆H₁₀₄Cl₂N₁₈O₁₃ • XCF₃COOH

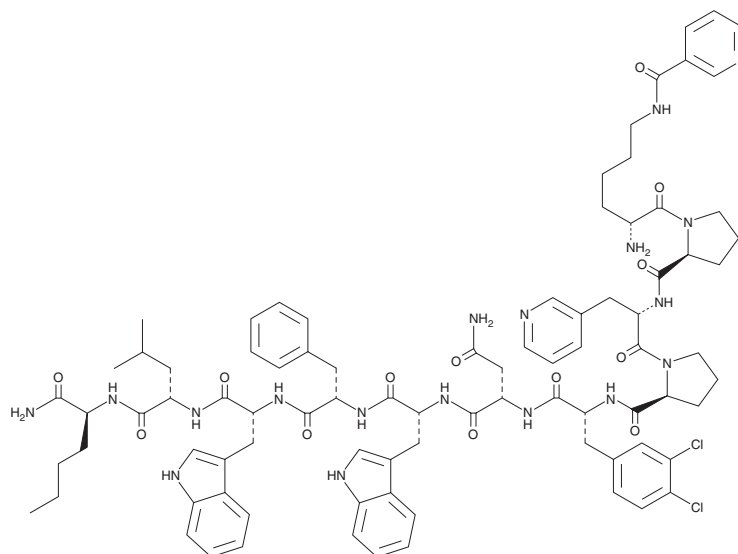
FW: 1,668.8

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

Spantide II (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the spantide II (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Spantide II (trifluoroacetate salt) is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Spantide II is a peptide neurokinin-1 (NK₁) and NK₂ receptor antagonist (K_s = 2.4 and 19.05 nM, respectively).¹ It induces histamine release from isolated rat mast cells (EC₅₀ = 3.2 μM).² Spantide II (10 μM) inhibits bombesin-induced contractions in isolated and perfused guinea pig taenia coli when used at a concentration of 10 μM. Topical application of spantide II (0.05-1% (w/v) three times per day) decreases ear swelling in a mouse model of allergic contact dermatitis induced by 2,4-dinitro-1-fluorobenzene.³

References

1. Janecka, A., Poels, J., Fichna, J., et al. *Regul. Pept.* **131(1-3)**, 23-28 (2005).
2. McLean, S., Ganong, A., Seymour, P.A., et al. *J. Pharmacol. Exp. Ther.* **267(1)**, 472-479 (1993).
3. Kikwai, L., Babu, R.J., Prado, R., et al. *AAPS PharmSciTech* **6(4)**, E565-E572 (2005).

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, 05/06/2025

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