

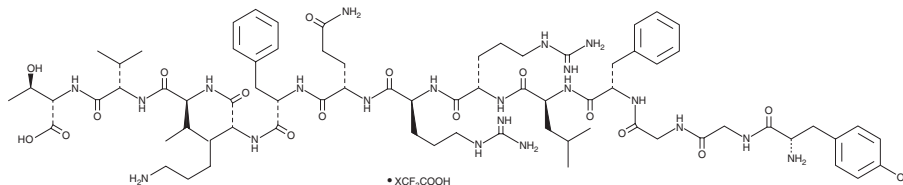
PRODUCT INFORMATION



Dynorphin B (trifluoroacetate salt)

Item No. 18178

Formal Name: dynorphin B (swine), trifluoroacetate salt
Synonyms: Dynorphin B-13, Rimorphin
MF: C₇₄H₁₁₅N₂₁O₁₇ • XCF₃COOH
FW: 1,570.8
Purity: ≥95%
UV/Vis.: λ_{max}: 278 nm
Supplied as: A crystalline 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

Dynorphin B (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the dynorphin B (trifluoroacetate salt) in the solvent of choice. Dynorphin B (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of dynorphin B (trifluoroacetate salt) in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of dynorphin B (trifluoroacetate salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of dynorphin B (trifluoroacetate salt) in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Dynorphin B is a 13-residue opioid peptide released from the cleavage of prodynorphin and found widely distributed in the central nervous system.^{1,2} It acts as a κ_{1b}-opioid receptor agonist (K_i = 1.1 nM) and has been implicated in antinociceptive functions.³ Dynorphin B has also been used to prime cardiogenesis in pluripotent embryonic stem cells.⁴

References

1. Berman, Y., Juliano, L., and Devi, L.A. Specificity of the dynorphin-processing endoprotease: Comparison with prohormone convertases. *J. Neurochem.* **72**, 2120-2196 (1999).
2. Janecka, A., Fichna, J., and Janecki, T. Opioid receptors and their ligands. *Curr. Top. Med. Chem.* **4**, 1-17 (2004).
3. Dhawan, B.N., Cesselin, F., Raghbir, R., *et al.* International Union of Pharmacology. XII. Classification of opioid receptors. *Pharm. Res.* **48(4)**, 567-592 (1996).
4. Ventura, C., Zinellu, E., Maninchedda, E., *et al.* Dynorphin B is an agonist of nuclear opioid receptors coupling nuclear protein kinase C activation to the transcription of cardiogenic genes in GTR1 embryonic stem cells. *Circ. Res.* **92**, 623-629 (2003).

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, 08/03/2023

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