

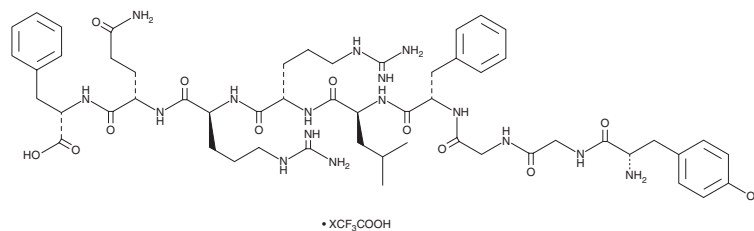
# PRODUCT INFORMATION



## Dynorphin B (1-9) (trifluoroacetate salt)

Item No. 24743

**MF:**  $C_{54}H_{78}N_{16}O_{12} \cdot XCF_3COOH$   
**FW:** 1,143.3  
**Purity:**  $\geq 95\%$   
**Supplied as:** A lyophilized powder  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Dynorphin B (1-9) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the dynorphin B (1-9) (trifluoroacetate salt) in water. The solubility of dynorphin B (1-9) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Dynorphin B (1-9) is a neuropeptide and N-terminal cleavage product of dynorphin B, an endogenous opioid.<sup>1</sup> Formation of dynorphin B (1-9) is blocked by N-ethylmaleimide (NEM), a non-selective inhibitor of cysteine peptidases. Dynorphin B (1-9) is a competitive inhibitor of dynorphin A (1-9), binding to the 60 kDa N-terminal fragment of puromycin-sensitive aminopeptidase (ApPS<sup>60</sup>).<sup>2</sup>

### References

1. Bivehed, E., Strömvall, R., Bergguist, J., *et al.* Region-specific bioconversion of dynorphin neuropeptide detected by *in situ* histochemistry and MALDI imaging mass spectrometry. *Peptides* **87**, 20-27 (2017).
2. Ma, Z., Daquin, A., Yao, J., *et al.* Proteolytic cleavage of the puromycin-sensitive aminopeptidase generates a substrate binding domain. *Arch. Biochem. Biophys.* **415**(1), 80-86 (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, 11/11/2022

#### 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](http://WWW.CAYMANCHEM.COM)