

PRODUCT INFORMATION



Pentapeptide-18 (acetate)

Item No. 35198

Formal Name: L-tyrosyl-D-alanylglycyl-L-phenylalanyl-L-leucine, acetate

Synonyms: [Ala²,Leu⁵]-Enkephalin,
[D-Ala²]Leucine-Enkephalin,
Tyr-D-Ala-Gly-Phe-Leu-OH

MF: C₂₉H₃₉N₅O₇ • XC₂H₄O₂

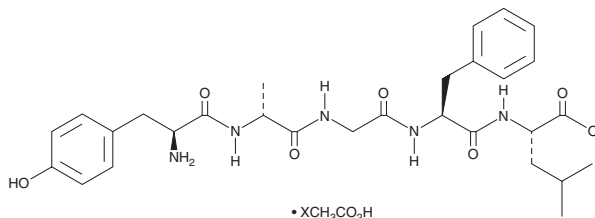
FW: 569.7

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

Pentapeptide-18 (acetate) is supplied as a solid. A stock solution may be made by dissolving the pentapeptide-18 (acetate) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Pentapeptide-18 is a peptide derivative of enkephalin.^{1,2} It reduces tail pinch-induced hyperalgesia in rats.² Formulations containing pentapeptide-18 have been used in cosmeceuticals.

References

1. Beddell, C.R., Clark, R.B., Hardy, G.W., *et al.* Structural requirements for opioid activity of analogues of the enkephalins. *Proc. R. Soc. Long. B. Biol. Sci.* **198(1132)**, 249-265 (1977).
2. Maeda, M., Kawasaki, K., Mayumi, T., *et al.* Amino acids and peptides. XXII. Preparation and antinociceptive effect of [D-Ala²]Leu-enkephalin-poly(ethylene glycol) hybrid. *Biol. Pharm. Bull.* **17(6)**, 823-825 (1994).

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, 12/12/2022

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