

PRODUCT INFORMATION



Acetyl Tetrapeptide-15 (acetate)

Item No. 35180

Formal Name: N-acetyl-L-tyrosyl-L-prolyl-L-phenylalanyl-L-phenylalaninamide, acetate

Synonyms: N-acetyl-L-Tyr-L-Pro-L-Phe-L-Phe-NH₂,
N-acetyl-YPFF-NH₂

MF: C₃₄H₃₉N₅O₆ • XC₂H₄O₂

FW: 613.7

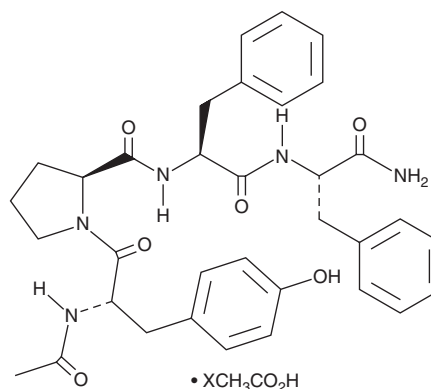
Purity: ≥95%

UV/Vis.: λ_{max}: 210 nm

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

Acetyl tetrapeptide-15 (acetate) is supplied as a solid. A stock solution may be made by dissolving the acetyl tetrapeptide-15 (acetate) in the solvent of choice, which should be purged with an inert gas. Acetyl tetrapeptide-15 (acetate) is soluble in DMSO.

Description

Acetyl tetrapeptide-15 is a peptide derivative of endomorphin 2 (Item No. 23281).¹ Formulations containing acetyl tetrapeptide-15 have been used in cosmeceuticals.

Reference

1. Resende, D.I.S.P., Ferreira, M.S., Sousa-Lobo, J.M., *et al.* Usage of synthetic peptides in cosmetics for sensitive skin. *Pharmaceuticals (Basel)* **14(8)**, 702 (2021).

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