

PRODUCT INFORMATION



Pentapeptide-3 (acetate)

Item No. 35186

Formal Name: glycyl-L-prolyl-L-arginyl-L-prolyl-L-alaninamide, acetate

Synonyms: H-GPRPA-NH₂,
H-Gly-Pro-Arg-Pro-Ala-NH₂

MF: C₂₁H₃₇N₉O₅ • XC₂H₄O₂

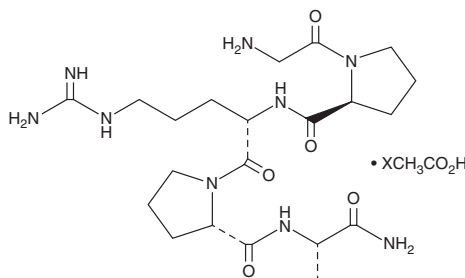
FW: 495.6

Purity: ≥95%

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-3 (acetate) is supplied as a solid. A stock solution may be made by dissolving the pentapeptide-3 (acetate) in the solvent of choice, which should be purged with an inert gas. Pentapeptide-3 (acetate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of pentapeptide-3 (acetate) in these solvents is approximately 10 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 pentapeptide-3 (acetate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of pentapeptide-3 (acetate) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Pentapeptide-3 is a cosmeceutical peptide and pentapeptide fragment of the temple viper (*T. wagleri*) venom peptide Waglerin-1.¹ It is an antagonist of nicotinic acetylcholine receptors (nAChRs).²

References

1. Schagen, S.K. Topical peptide treatments with effective anti-aging results. *Cosmetics* **4(2):16**, (2017).
2. Zhmak, M.N., Utkin, Y.N., Andreeva, T.V., et al. Peptide inhibitors of nicotinic acetylcholine receptor. *Syneuro LLC US 9,550,808 B2*, (2017).

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, 11/02/2023

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