

PRODUCT INFORMATION



Acetyl Hexapeptide-3 (acetate)

Item No. 27756

Formal Name: N-acetyl-L- α -glutamyl-L- α -glutamyl-L-methionyl-L-glutaminyl-L-arginyl-L-argininamide, acetate

Synonyms: Ac-EEMQRR-NH₂, Acetyl-Glu-Glu-Met-Gln-Arg-Arg-NH₂, Acetyl Hexapeptide-8

MF: C₃₄H₆₀N₁₄O₁₂S • XC₂H₄O₂

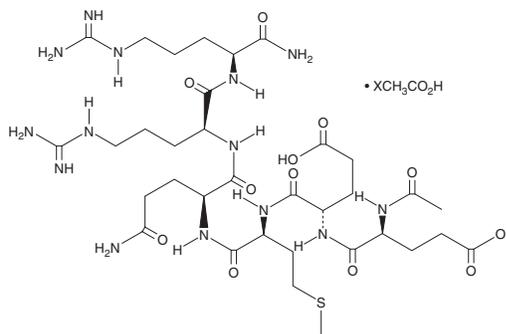
FW: 889.0

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

Acetyl hexapeptide-3 (acetate) is supplied as a solid. Aqueous solutions of acetyl hexapeptide-3 (acetate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of acetyl hexapeptide-3 (acetate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Acetyl hexapeptide-3 is a synthetic acetylated and amidated hexapeptide consisting of amino acid residues 12-17 of synaptosomal-associated protein 25 (SNAP-25).¹ It inhibits formation of the SNARE complex *in vitro* in a concentration-dependent manner. It also inhibits calcium-induced norepinephrine release in permeabilized chromaffin cells (IC₅₀ = 110 μ M). Acetyl hexapeptide-3 (10% in an oil-and-water emulsion) inhibits decreases in type I collagen levels and skin thickness, as well as increases in type III collagen levels, in skin in a mouse model of subacute aging induced by D-galactose (Item No. 20890) when administered topically.²

References

1. Blanes-Mira, C., Clemente, J., Jodas, G., *et al.* A synthetic hexapeptide (argireline) with antiwrinkle activity. *Int. J. Cosmet. Sci.* **24**(5), 303-310 (2002).
2. Wang, Y., Wang, M., Xiao, X.S., *et al.* The anti-wrinkle efficacy of argireline. *J. Cosmet Laser Ther.* **15**(4), 237-241 (2013).

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, 12/02/2022

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