

PRODUCT INFORMATION

Amyloid-β Precursor Protein (96-110) Peptide (cyclized) (human) (trifluoroacetate salt)

Item No. 27419

Formal Name: N²-acetyl-L-asparaginyl-L-tryptophyl-L-cysteinyl-L-lysyl-L-arginylglycyl-L-arginyl-L-lysyl-L-glutaminyll-L-cysteinyl-L-lysyl-L-threonyl-L-histidyl-L-prolyl-L-histidinamide cyclic (3→10)-disulfide, trifluoroacetate salt
Synonyms: Cyclized APP (96-110), Cyclized APP₉₆₋₁₁₀

MF: C₈₁H₁₂₈N₃₂O₁₉S₂ • XCF₃COOH

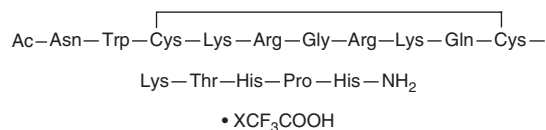
FW: 1,918.2

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

Amyloid-β precursor protein (96-110) peptide (cyclized) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the amyloid-β precursor protein (96-110) peptide (cyclized) (human) (trifluoroacetate salt) in water. The solubility of amyloid-β precursor protein (96-110) peptide (cyclized) (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Amyloid-β precursor protein (96-110) peptide (cyclized) is a synthetic peptide consisting of amino acids 96-110 of amyloid precursor peptide (APP) that is cyclized *via* a bridge between the cysteine residues at positions 3 and 10.¹ It is homologous to the heparin-binding domain of human APP and inhibits binding of [¹²⁵I]APP to heparin *in vitro* with an IC₅₀ value of approximately 100 nM. Amyloid-β precursor protein (96-110) peptide (cyclized) inhibits neurite outgrowth induced by APP and mouse postnatal heparan sulfate proteoglycans (HSPG) in primary mouse hippocampal neurons when used at a concentration of 10 μg/ml.

Reference

- Small, D.H., Nurcombe, V., Reed, G., *et al.* A heparin-binding domain in the amyloid protein precursor of Alzheimer's disease is involved in the regulation of neurite outgrowth. *J. Neurosci.* **14(4)**, 2117-2127 (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, 11/30/2022

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