

PRODUCT INFORMATION



Amyloid- β (22-35) Peptide (trifluoroacetate salt)

Item No. 27417

Formal Name:	L- α -glutamyl-L- α -aspartyl-L-valylglycyl-L-seryl-L-asparaginyl-L-lysylglycyl-L-alanyl-L-isoleucyl-L-isoleucylglycyl-L-leucyl-L-methionine, trifluoroacetate salt	H—Glu—Asp—Val—Gly—Ser—Asn—Lys—Gly—Ala—Ile—
Synonym:	A β (22-35)	Ile—Gly—Leu—Met—OH
MF:	C ₅₉ H ₁₀₂ N ₁₆ O ₂₁ S • XCF ₃ COOH	• XCF ₃ COOH
FW:	1,403.6	
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

Amyloid- β (22-35) peptide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the amyloid- β (22-35) peptide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Amyloid- β (22-35) peptide (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

Amyloid- β (A β) (22-35) is a 13-residue fragment of A β that corresponds to residues 693-705 of the human amyloid precursor protein (APP) full-length sequence.¹ It assembles straight fibrils and bundles that bind thioflavin T *in vitro*. A β (22-35) (40 μ g/ml) induces cell death in rat hippocampal neurons.

Reference

1. Takadera, T., Sakura, N., Mohri, T., *et al.* Toxic effect of a β -amyloid peptide (β 22-35) on the hippocampal neuron and its prevention. *Neurosci. Lett.* **161(1)**, 41-44 (1993).

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, 12/22/2022

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