

PRODUCT INFORMATION



Amyloid- β (40-1) Peptide (human) (trifluoroacetate salt)

Item No. 27112

Formal Name: L-valyl-L-valylglycylglycyl-L-valyl-L-methionyl-L-leucylglycyl-L-isoleucyl-L-isoleucyl-L-alanylglycyl-L-lysyl-L-asparaginyll-L-serylglycyl-L-valyl-L- α -aspartyl-L- α -glutamyl-L-alanyl-L-phenylalanyl-L-phenylalanyl-L-valyl-L-leucyl-L-lysyl-L-glutaminyll-L-histidyl-L-histidyl-L-valyl-L- α -glutamyl-L-tyrosylglycyl-L-seryl-L- α -aspartyl-L-histidyl-L-arginyl-L-phenylalanyl-L- α -glutamyl-L-alanyl-L-aspartic acid, trifluoroacetate salt



Synonym: A β (40-1)

MF: C₁₉₄H₂₉₅N₅₃O₅₈S • XCF₃COOH

FW: 4,329.9

Purity: \geq 95%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

• XCF₃COOH

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Amyloid- β (40-1) (A β (40-1)) peptide (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the A β (40-1) peptide (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. A β (40-1) peptide (human) (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

A β (40-1) peptide is a peptide that contains the reverse sequence of A β (1-40) peptide (Item No. 21617).¹ A β (40-1) does not induce neurotoxicity of rat cortical neurons when administered *via* microinjection into the cerebral cortex at a dose of 3 nmol. It has been used as an inactive control for A β (1-40) peptide.²

References

1. Kowall, N.W., Beal, M.F., Busciglio, J., *et al.* An *in vivo* model for the neurodegenerative effects of β amyloid and protection by substance P. *Proc. Natl. Acad. Sci. U.S.A.* **88**(16), 7247-7251 (1991).
2. Terranova, J.-P., Kan, J.-P., Storme, J.-J., *et al.* Administration of amyloid β -peptides in the rat medial septum causes memory deficits: Reversal by SR 57746A, a non-peptide neurotrophic compound. *Neurosci. Lett.* **213**(2), 79-82 (1996).

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