

PRODUCT INFORMATION



Amyloid-β (1-28) Peptide (human) (trifluoroacetate salt)

Item No. 27107

Formal Name:	L-α-aspartyl-L-alanyl-L-α-glutamyl-L-phenylalanyl-L-arginyl-L-histidyl-L-α-aspartyl-L-serylglycyl-L-tyrosyl-L-α-glutamyl-L-valyl-L-histidyl-L-histidyl-L-glutamyl-L-lysyl-L-leucyl-L-valyl-L-phenylalanyl-L-phenylalanyl-L-alanyl-L-α-glutamyl-L-α-aspartyl-L-valylglycyl-L-seryl-L-asparaginyl-L-lysine, trifluoroacetate salt	H—Asp—Ala—Glu—Phe—Arg—His—Asp—Ser—Gly—Tyr— Glu—Val—His—His—Gln—Lys—Leu—Val—Phe—Phe— Ala—Glu—Asp—Val—Gly—Ser—Asn—Lys—OH
Synonyms:	Aβ (1-28), Aβ28	
MF:	C ₁₄₅ H ₂₀₉ N ₄₁ O ₄₆ • XCF ₃ COOH	• XCF ₃ COOH
FW:	3,262.5	
Purity:	≥95%	
Supplied as:	A crystalline 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 (1-28) peptide (human) (trifluoroacetate salt) is supplied as a crystalline solid. Aqueous solutions of amyloid-β (1-28) peptide (human) (trifluoroacetate salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of amyloid-β (1-28) peptide (human) (trifluoroacetate salt) in PBS, pH 7.2, is approximately 0.15 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Amyloid-β (1-28) (Aβ28) is a synthetic peptide that lacks 14 C-terminal amino acids and is more soluble than Aβ42 (Item No. 20574).¹ Aβ28 forms fibrils and plaques *in vitro* that are similar to those formed by Aβ42 and induces learning deficits in a passive avoidance task in rats.²⁻⁴

References

1. Syme, C.D., Nadal, R.C., Rigby, S.E., *et al.* Copper binding to the amyloid-beta (Aβ) peptide associated with Alzheimer's disease. Folding, coordination geometry, pH dependence, stoichiometry, and affinity of Aβ(1-28): Insights from a range of complementary spectroscopic techniques. *J. Biol. Chem.* **279**(18), 18169-18177 (2004).
2. Kirschner, D.A., Inouye, H., Duffy, L.K., *et al.* Synthetic peptide homologous to β protein from Alzheimer disease forms amyloid-like fibrils *in vitro*. *Proc. Natl. Acad. Sci. USA* **84**(19), 6953-6957 (1987).
3. Burdick, D., Soreghan, B., Kwon, M., *et al.* Assembly and aggregation properties of synthetic Alzheimer's A4/β amyloid peptide analogs. *J. Biol. Chem.* **267**(1), 546-554 (1992).
4. Alvarez, X.A., Miguel-Hidalgo, J.J., Fernández-Novoa, L., *et al.* Intrahippocampal injections of the beta-amyloid 1-28 fragment induces behavioral deficits in rats. *Methods Find. Exp. Clin. Pharmacol.* **19**(7), 471-479 (1997).

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, 12/21/2022

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