

PRODUCT INFORMATION



TAMRA-Amyloid- β (1-42) Peptide (trifluoroacetate salt)

Item No. 27409

Synonyms:	TAMRA-Amyloid β (1-42), TAMRA-A β 42	TAMRA—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—Gly—Ala— Ile—Ile—Gly—Leu—Met—Val—Gly—Gly—Val—Val— Ile—Ala—OH • XCF ₃ COOH
MF:	C ₂₂₈ H ₃₃₁ N ₅₇ O ₆₄ S • XCF ₃ COOH	
FW:	4,926.5	
Purity:	≥95%	
Ex./Em. Max:	543/572 nm	
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

TAMRA-Amyloid- β (1-42) peptide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the TAMRA-amyloid- β (1-42) peptide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. TAMRA-Amyloid- β (1-42) peptide (trifluoroacetate salt) is soluble in the organic solvent formic acid, at a concentration of approximately 1 mg/ml.

Description

TAMRA-Amyloid- β (1-42) peptide is a fluorescently labeled peptide. Amyloid- β (1-42) (A β 42) is a neurotoxic 42-residue protein fragment found in amyloid plaques in postmortem cerebral cortex from patients with Alzheimer's disease.^{1,2} Aggregation of A β 42 results in the formation of neurotoxic fibrils or globular oligomers.¹ TAMRA-Amyloid- β (1-42) peptide is a labeled form of A β 42 containing carboxytetramethyl rhodamine (TAMRA), which displays excitation/emission maxima of 543/572 nm, respectively.

References

1. Wolfe, M.S. Therapeutic strategies for Alzheimer's disease. *Nat. Rev. Drug Discov.* **1**(11), 859-866 (2002).
2. Iwatsubo, T., Odaka, A., Suzuki, N., *et al.* Visualization of A β 42(43) and A β 40 in senile plaques with end-specific A β monoclonals: Evidence that an initially deposited species is A β 42(43). *Neuron* **13**(1), 45-53 (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