

PRODUCT INFORMATION



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

Item No. 27411

Formal Name: N-[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]-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-asparaginy-L-lysylglycyl-L-alanyl-L-isoleucyl-L-isoleucylglycyl-L-leucyl-L-methionyl-L-valylglycylglycyl-L-valyl-L-valine, trifluoroacetate salt

Biotin—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—OH
• XCF₃COOH

Synonyms: Biotin-A β (1-40), Biotin-A β 40
MF: C₂₀₄H₃₀₉N₅₅O₆₀S₂ • XCF₃COOH
FW: 4,556.1
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years

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

Laboratory Procedures

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

Description

Biotin-A β 40 is an affinity probe that allows amyloid- β (1-40) (A β 40) to be detected or immobilized through interaction with the biotin ligand. It has been used to characterize A β 40 distribution amongst the lipoprotein fractions and identify A β 40 interaction partners in human plasma.¹

Reference

1. Koudinov, A., Matsubara, E., Frangione, B., *et al.* The soluble form of Alzheimer's amyloid beta protein is complexed to high density lipoprotein 3 and very high density lipoprotein in normal human plasma. *Biochem. Biophys. Res. Commun.* **205**(2), 1164-1171 (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