

# PRODUCT INFORMATION



## Biotin-Amyloid- $\beta$ (1-42) Peptide (trifluoroacetate salt)

Item No. 27410

**Formal Name:** N-[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]-L- $\alpha$ -aspartyl-L-alanyl-L- $\alpha$ -glutamyl-L-phenylalanyl-L-arginyl-L-histidyl-L- $\alpha$ -aspartyl-L-serylglycyl-L-tyrosyl-L- $\alpha$ -glutamyl-L-valyl-L-histidyl-L-histidyl-L-glutamyl-L-lysyl-L-leucyl-L-valyl-L-phenylalanyl-L-phenylalanyl-L-alanyl-L- $\alpha$ -glutamyl-L- $\alpha$ -aspartyl-L-valylglycyl-L-seryl-L-asparaginyl-L-lysylglycyl-L-alanyl-L-isoleucyl-L-isoleucylglycyl-L-leucyl-L-methionyl-L-valylglycylglycyl-L-valyl-L-valyl-L-isoleucyl-L-alanine, 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  
• XCF<sub>3</sub>COOH

**Synonyms:** Biotin-A $\beta$ (1-42), Biotin-A $\beta$ 42

**MF:** C<sub>213</sub>H<sub>325</sub>N<sub>57</sub>O<sub>62</sub>S<sub>2</sub> • XCF<sub>3</sub>COOH

**FW:** 4,740.4

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

### Description

Biotin-amyloid- $\beta$  (1-42) peptide is an affinity probe that allows amyloid- $\beta$  (1-42) (A $\beta$ 42) to be detected or immobilized through interaction with the biotin ligand. It has been used to identify A $\beta$ 42 interaction partners in rat hippocampal synaptosomal membranes.<sup>1</sup>

### References

1. Jerecic, J. and Krafft, G.A. ADDL receptor polypeptides, polynucleotides and host cells for recombinant production. *Acumen Pharmaceuticals, Inc.* US 9,217,024 B2 (2015).

#### 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, 10/25/2022

#### 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