

Amyloid beta 42, N-Alpha 5-TAMRA labeled

Synthetic peptide derived from human amyloid (1-42)

Catalog # A06-58G

Lot # T2065-7

Product Description

The synthetic Amyloid β 42 Peptide (Modified DAEFRHDSGY EVHHQKLVFF AEDVGSNKGAIIGLMVGGVV IA) is derived from human amyloid β (amino acid 1-42).

Molecular Weight

The theoretical molecular weight is 4926.50

Purity

The purity was determined to be $\geq 95\%$ by HPLC analysis.

Storage and Stability

Store product at -20°C . For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

Amyloid peptides, derived from amyloid precursor protein (APP), are thought to play a role in the development of the senile plaques associated with Alzheimer's Disease. The amyloid hypothesis presupposes that flaws in the processing of APP result in abnormally high levels of the longer, "stickier" forms of beta amyloid, known as A β 42 and A β 43, leading to aggregation of amyloid in the neuronal cell death and ultimately neuronal death. Mutations in the structure of A β 40 and related peptides as well as in some of the enzymes involved in the processing of APP have been shown to alter the processing of APP. The sporadic (i.e., non-genetic) form of the disease, however, is far more common, caused by aging in concert with a number of both genetic and environmental risk factors.

References

1. Esler WP, et al: Stereochemical specificity of Alzheimer's disease beta-peptide assembly. *Biopolymers*. 1999 May;49(6):505-14.
2. Roher AE, et al: beta-Amyloid-(1-42) is a major component of cerebrovascular amyloid deposits: implications for the pathology of Alzheimer disease. *Proc Natl Acad Sci U S A*. 1993 Nov 15;90(22):10836-40.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Formulation

The peptide supplied in trifluoroacetate salt as a lyophilized powder.

Reconstitution Protocol

Dilute peptide in distilled water to make the working solution.

Amyloid beta 42, N-Alpha 5-TAMRA labeled

Synthetic peptide derived from human amyloid (1-42)

Catalog Number	A06-58G
Specific Lot Number	T2065-7
Purity	> 95%
Format	Lyophilized powder
Stability	1 yr At -20°C from date of shipment
Storage & Shipping	Store product at -20°C . For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped at ambient temperature.