

Product Data Sheet

Product Name: β-Amyloid (12-28)

Catalog Number: AS-24229 (0.5 mg) Lot Number: See label on vial

AS-24230 (1 mg)

Sequence: H-Val-His-His-Gln-Lys-Leu-Val-Phe-Phe-Ala-Glu-Asp-Val-Gly-Ser-Asn-

Lys-OH (3-letter code)

VHHQKLVFFAEDVGSNK (1-letter code)

Molecular Weight: 1955.2

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: β-Amyloid (12-28) peptide is freely soluble in water.

Storage: β -Amyloid (12-28) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20° C or lower. Reconstituted peptide can be aliquoted and stored at -20° C or lower.

Description: Aβ (12–28) residues are the binding site for apolipoprotein E (apoE) on Aβ. This sequence encompasses a hydrophobic domain (residues 14–21) and a β-turn (residues 22–28) which place two hydrophobic domains of Aβ 14 to 21 and 29 to 40/42 opposite each other, allowing for the assembly of Aβ peptides into fibrils. The secondary structure of Aβ (12–28), a neutral peptide, is dominated by a-helix and random coil. The interaction of apoE with residues 12 to 28 of Aβ is not just a non-specific hydrophobic interaction but plays a pivotal role in the mechanism of Aβ pathology in Alzheimer's disease (AD). Aβ (11-28) and five other fragments enhanced aggregation of full length Aβ (1-40). All of the peptides that enhance aggregation contained either residues 17 to 20 or 30 to 35, indicating the importance of these regions for promoting aggregation of full-length Aβ. Ref: Sadowski, M. et al. *Am. J. Pathol.* 165, 937 (2004); Liu, R. et al. J. *Neurosci. Res.* 75, 162 (2004).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Aβ₁₋₁₁, Aβ₁₀₋₂₀, Aβ₁₅₋₂₀, Aβ₁₂₋₂₈, Aβ₂₅₋₃₅, Aβ₃₇₋₄₃, Aβ₂₉₋₄₀, biotinated Aβ₁₋₄₂, and FITC-conjugated Aβ₁₋₄₂ were obtained from AnaSpec. o determine whether exposure to exogenous Aβ₄₂ increases Aβ₄₂—α7nAChR association and causes Aβ₄₂-induced α7nAChR and NMDAR dysfunction, ~20 mg of FCX slices from either control subjects or AD individuals were incubated with 0.1 μM Aβ₄₂ at 37°C for 1 h. To test their effects, the following drugs were added immediately after Aβ₄₂: S 24795 (1–100 μM), Aβ₁₂₋₂₈ (10 μM), memantine (30 μM), galantamine (30 μM), PNU 282987 (30 μM), MLA (10 μM), or MLA (10 μM) plus S 24795 (10 μM). Incubation continued for 1 h in the dark to minimize light destruction of the test agents such as S 24795. The incubation mixture in a total incubation volume of 0.5 ml was aerated with 95% $O_2/5\%$ CO_2 every 15 min for

1 min during the incubation. Reaction was terminated by the addition of 1.5 ml of ice-cold Ca²⁺-free K-R. Tissue slices were harvested by brief centrifugation and used as the tissue sources for various assays-Wang, H.Y. et al. *J Neuro* **10**, 10961 (2009).

Published Citations:

Mouedden, M. et al. *J. Neuro.* **145**, 97 (2005). Osada, Y. et al. *JBC* **280**, 8596 (2005). Solorzano-Vargas, RS. et al. *Mole. Immunol.* **45**, 881 (2008). Wang, H.Y. et al. *J Neuro* **10**, 10961 (2009).

For Research Use Only