

## **Product Data Sheet**

Product Name: β-Amyloid (1-42), 5-FAM-labeled

Catalog Number: AS-23526-01 (0.1 mg) Lot Number: See label on vial

AS-23525-05 (0.5 mg)

Sequence: 5-FAM-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 (3-letter code)

5-FAM-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA

(1-letter code)

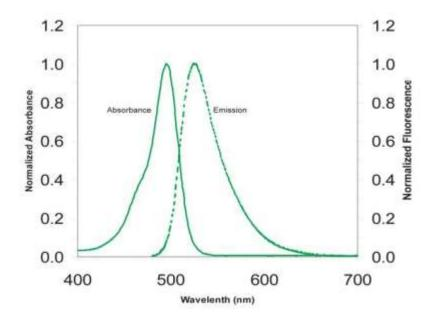
Molecular Weight: 4872.4 % Peak Area by HPLC:≥95

Appearance: Lyophilized yellow powder

Peptide Reconstitution: Use 1.0% NH4OH as the solvent, followed by buffer (i.e.1XPBS). Add 1.0% NH4OH directly to the lyophilized peptide powder (add 35-40  $\mu$ l to 0.5 mg peptide or 70-80  $\mu$ l to 1 mg peptide). The peptide cannot be stored long term in 1.0% NH4OH, and it is therefore important to immediately dilute this solution with 1X PBS or other buffer to a concentration of approximately 1mg/mL or less. Gently vortex to mix.

Storage: Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at – 20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at –20°C or lower. Do not freeze thaw.

Description: This is a fluorescent (FAM)-labeled β-Amyloid peptide, Abs/Em=494/521 nm. Ref: Wei, H. et al. *J. Pharmacol.* **392**, 117 (2000).



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.

Fluorescein-labeled  $\beta$ -Amyloid peptide A $\beta$  1-42 (fluo A $\beta$  1-42) was purchased from AnaSpec(San Jose, CA, USA). Fluo-A $\beta$  (1-42) was provided to sympathetic neurons cultured in 3-compartment cultures or 8-well slides. The internalization of fluo-A $\beta$ (1-42) in living neurons was examined using a Nikon TE300 inverted fluorescence microscope - Saavedra, L. et al. J Biol Chem 10.1074/jbc.M701823200 (2007).

## **Published Citations:**

Choucair, A. et al. *Biochim Biophys Acta* **1768**, 146 (2007). Saavedra, L. et al. *J Biol Chem* 10.1074/jbc.M701823200 (2007). LeVine III, H. *Anal. Biochem* **335**, 81 (2004).

For Research Use Only