



## Product Data Sheet

**Product Name:**  $\beta$ -Amyloid (1-42), 5-FAM-labeled

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

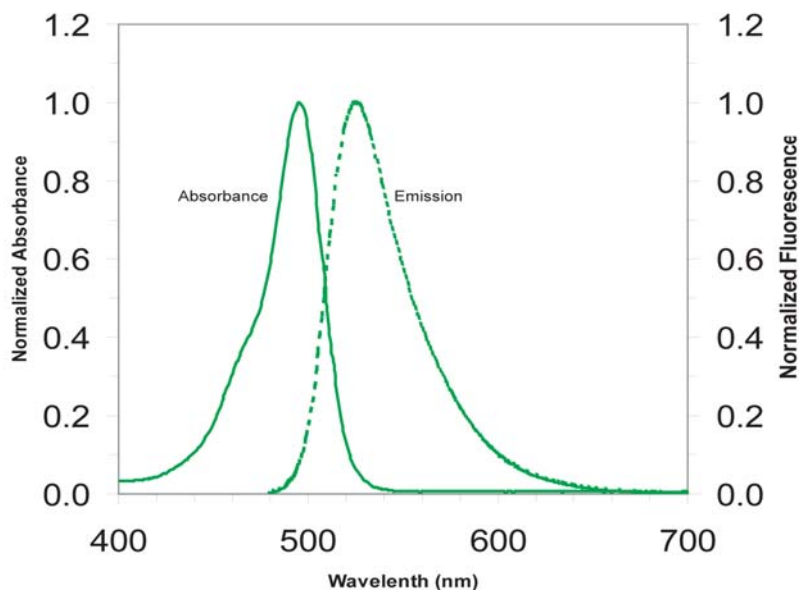
**% Peak Area by HPLC:**  $\geq 95$

**Appearance:** Lyophilized yellow powder

**Peptide Reconstitution:** Use 1.0% NH<sub>4</sub>OH as the solvent, followed by buffer (i.e. 1X PBS). Add 1.0% NH<sub>4</sub>OH 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% NH<sub>4</sub>OH, 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  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at  $-20^{\circ}\text{C}$  or lower. Do not freeze thaw.

**Description:** This is a fluorescent (FAM)-labeled  $\beta$ -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 \(2007\).](#)

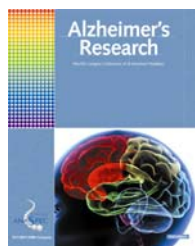
**Published Citations:**

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).

**Related Products:**



Click Alzheimer's Research Brochure to view the full listing.

***For Research Use Only***