

Product Name:	β-Amyloid (1-40), HiLyte Fluor™488-labeled	d
Catalog Number:	AS-60491-01 (0.1 mg)	Lot Number: See label on vial
Sequence:	HiLyte Fluor <sup>™</sup> 488-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 (3-letter code) HiLyte Fluor <sup>™</sup> 488-DAEFRHDSGYEVHHQKLVFFAEDVGSNKG AIIGLMVGGVV (1-letter code)	
Molecular Weight:	4687.3	
% Peak Area by HPLC:	≥ 95	
Appearance:	Lyophilized orange powder	

Peptide Reconstitution: Reconstitute by adding 35-40  $\mu$ I 1%NH<sub>4</sub>OH to 0.1 mg  $\beta$ -Amyloid (1-40). Dilute this peptide solution to approximately 1 mg/ml (or more dilute) with a buffer such as PBS or another buffer; aliquot and store at -20C.

Storage:  $\beta$ -Amyloid (1-40) 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: This is a fluorescent (HiLyte Fluor™ 488)-labeled β-Amyloid peptide, Abs/Em=503/528 nm.

## 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.

 $A\beta_{1-40}$ -HiLyte Fluor (Alexa Fluor 488) peptide was freshly dissolved in PBS and added to complete culture medium (0.5 ml) at a final concentration of 1 µM for 18 h at 37°C. Subsequently, medium was changed and 585-nM MitoTracker Orange (Molecular Probes Inc.) was added for 30 min at 37°C. Cells were incubated for another 15 min in new medium before washing with PBS and fixation in 2% paraformaldehyde for 5 min-<u>Petersen, CAH. et al. *PNAS* **105**, 13145 (2008).</u>

To test whether adsorption of A $\beta_{40}$  at the aqueous/fluorous interface can be minimized, we compared the behavior of A $\beta_{40}$  labeled with HiLyte-488 at the N terminus at two liquid/liquid interfaces: 1) aqueous peptide/fluorocarbon interface and 2) aqueous peptide/R<sub>f</sub>-OEG<sub>3</sub>-protected fluorocarbon interface, where R<sub>f</sub>-OEG<sub>3</sub> is an amphiphilic fluorinated surfactant (see Supporting Information) that is added to the carrier fluid, assembles spontaneously at the aqueous/fluorous interface, presents triethylene glycol groups to the aqueous phase, and thereby prevents protein adsorption-Meier, M. et al. Angew Chem Int Ed Engl. 48, 1487 (2009).

## **Published Citations:**

Petersen, CAH. et al. *PNAS* **105**, 13145 (2008). Meier, M. et al. *Angew Chem Int Ed Engl.* **48**, 1487 (2009).

## **Related Products:**

<b>Name</b> β-Amyloid (1-40), HiLyte Fluor™ 555-labeled	Cat # AS-60492-01	<b>Size</b> 0.1 mg
β-Amyloid (1-40), HiLyte Fluor™ 647-labeled	AS-60493	0.1 mg
β-Amyloid (1-40), FAM-labeled	AS-23514-01 AS-23513-05	0.1 mg 0.5 mg
β-Amyloid (1-40), TAMRA-labeled	AS-60488-01	0.1 mg
SensoLyte® 520 β-Secretase Assay Kit	AS-71144	1 kit

For Research Use Only