



Product Data Sheet

Product Name: β -Amyloid (1-42), HiLyte Fluor™ 488-labeled

Catalog Number: AS-60479-01 (0.1 mg) Lot Number: See label on vial

Sequence: HiLyte Fluor™ 488-labeled-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)
HiLyte Fluor™ 488-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA (1-letter code)

Molecular Weight: 4870.5

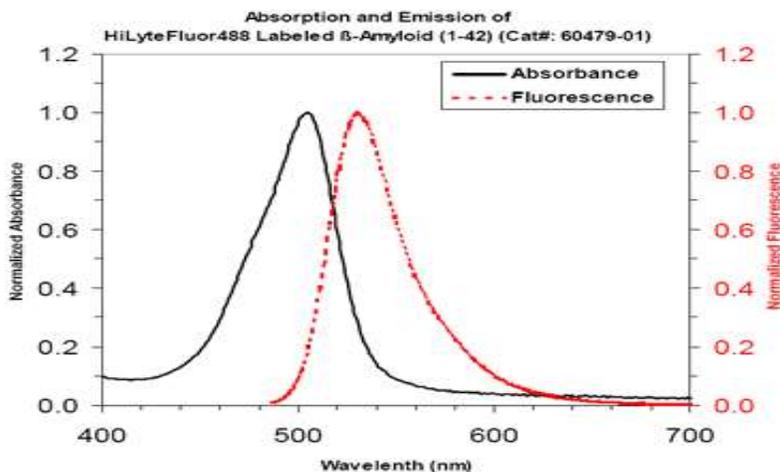
% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized orange-pink color powder

Peptide Reconstitution: Reconstitute by adding 50 μ l 1%NH₄OH to 0.1 mg β -Amyloid (1-42), HiLyte Fluor™ 488-labeled peptide. 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: β -Amyloid (1-42), HiLyte Fluor™ 488-labeled 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. HiLyte 488™ Fluor labeled β -Amyloid (1-42) has a brighter intensity than β -Amyloid (1-42) 5-FAM-labeled.



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.

β -Amyloid (1-42), HiLyte Fluor™ 488-labeled (AnaSpec, San Jose, CA) were added at various timepoints, and cells were washed twice with PBS and then removed from the plate using 0.25% trypsin/EDTA solution- [Nazer, B. et al. *Neurobio Dis.* **30**, 94 \(2008\).](#)

Fluorescence-labeled A β -42 (HiLyte Fluor™488- β -Amyloid(1–42); Anaspec Inc., CA, USA) were prepared 5:1 (w/w) in DMSO at 200 μ M concentration- [Vestergaard, M. et al. *Biochem & Biophys Res Com.* **377**, 725 \(2008\).](#)

Fluorescence-labeled A β -42 (HiLyte Fluor™488- β -Amyloid(1–42); Anaspec Inc., CA, USA) were prepared 5:1 (w/w) in DMSO at 200 μ M concentration. Final working concentration of the A β -42 and probe were 80 and 8 μ M, respectively in 20 mM Tris/HCl buffer, pH 7.4 (TBS) for fluorescence imaging studies. This solution was allowed to spontaneously aggregate in TBS at 37 ± 1 °C for a defined period of time and analysed using various techniques. Unless otherwise stated, all analyses were carried out at RT.-[Vestergaard, M. et al. *Biochem & Biophys Res Com.* **377**, 725 \(2008\).](#)

Published Citations:

[Hickman, SE. et al. *Neurobio. Dis.* **28**, 8354 \(2008\).](#)

[Nazer, B. et al. *Neurobio Dis.* **30**, 94 \(2008\).](#)

[Vestergaard, M. et al. *Biochem & Biophys Res Com.* **377**, 725 \(2008\).](#)

For Research Use Only