



## Product Data Sheet

Product Name:	$\beta$ -Amyloid (1-42)	
Catalog Number:	AS-24224 (0.5 mg)	Lot Number: See label on vial
	AS-20276 (1 mg)	
	AS-20276-5 (5 mg)	
	AS-20276-25 (25 mg)	

Sequence: H-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)  
DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVIA (1-letter code)

Molecular Weight: 4514.1

% Peak Area by HPLC:  $\geq$ 95

Peptide Content:  $\geq$  60%

Appearance: Lyophilized white 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:** A $\beta$  (1-42), a major component of amyloid plaques, accumulates in neurons of Alzheimer's disease brains. Biochemical analysis of the amyloid peptides isolated from Alzheimer's disease brain indicates that A $\beta$  (1-42) is the principal species associated with senile plaque amyloids, while A $\beta$  (1-40) is more abundant in cerebro-vascular amyloid deposits. Ref: Nagele, R. et al. *Nseurosci* **110**, 199 (2002); Garzon-Rodriguez, W. et al. *J Biol Chem* **272**, 21037 (1997).

**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.

Amyloid  $\beta$ (1–42) peptide was purchased from AnaSpec (San Jose, CA, USA). A 1 mg sample of peptide was dissolved in 200  $\mu$ L hexafluoroisopropanol (HFIP) and aliquoted to obtain 0.1 mg stocks (handle HFIP in a chemical fume hood taking the necessary precautions) - [Reinke, AA. et al. Chem Biol Drug Design](#) **70**, 206 (2007).

Synthetic A $\beta$  (1–42) (AnaSpec, San Jose, CA) was prepared for aggregation by resuspending lyophilized A $\beta$  (1–42) in hexafluoroisopropanol, dried under a nitrogen stream, and stored as a film at –20 °C. Immediately prior to use, A $\beta$  (1–42) was resuspended in Me SO to 10<sub>2</sub> mM and sonicated for 10 min. For experiments in which early stages of aggregation were studied, these aliquots were rapidly brought to 25 uM in phosphate-buffered saline (PBS), pH 7.2, and used immediately. Oligomers were prepared by diluting the A $\beta$  (1–42) to 25 uM with phenol red-free DMEM-F12 and incubating for 24 h at 4 °C without shaking. Fibrils were similarly prepared

by incubating 25  $\mu$ M A $\beta$  (1–42) in PBS at 37 °C for 24 h with vigorous shaking - [Evans, CG. et al.](#)  
[J Biol Chem 281, 33182 \(2006\).](#)

The ability to produce the presumably neurotoxic, aggregated  $\beta$ -sheet structure is dependent on many factors, particularly the peptide concentration ionic strength, solvent polarity. For example, the aggregation rate is extremely rapid in aqueous acetonitrile solutions, such as those used for HPLC purification of the peptide. The longer that the A $\beta$  (1-42) peptide remains in aqueous acetonitrile solution, the more likely it will become an aggregated  $\beta$ -sheet structure. Additionally, different commercially prepared batches of HPLC-purified Ab (1-42) peptides can have different starting aggregations states and structures, which will then in turn affect their solubility, aggregation rates, biological activities in solution and the ability to reproduce biophysical measurements. To partly overcome the above complications, we develop a pretreatment method that involves sonicating the dry peptide in conc. TFA before biophysical measurements. TFA breaks up the pre-aggregated peptides and affords monomeric random coil structures. This method ensures that different batches of purified Ab (1-42) peptide will provide reproducible starting points for biophysical and neurotoxicity studies. - [Salomon, AR. et al.](#)  
[Biochem 35, 13568 \(1996\).](#)

Solubility of A $\beta$  (1-42) is pH and concentration dependent. It is significantly insoluble at pH 7.4; it is highly insoluble in aqueous media but are soluble at 40 mg/ml in the  $\alpha$ -helix-promoting solvent, 1,1,1,3,3,3-hexafluoro-2-propanol (HFIP, handle HFIP in a fumehood and take the necessary precautions) - [Burdick, D. et al. J Biol Chem 267, 546 \(1992\).](#)

#### Published Citations:

- [Davis, RC. et al. Mol Neurodegener 6, 10 \(2011\).](#)  
[Dillen, L. et al. Bioanalysis 3, 45 \(2011\).](#)  
[Li, W. et al. Brain Res 10.1016/j.brainres.2011.02.078 \(2011\).](#)  
[Pandya, A. et al. J Mol Neurosci 10.1007/s12031-011-9509-3 \(2011\).](#)  
[Yin, F. et al. Neurochem Int 58, 399 \(2011\).](#)  
[Aksenov, M. et al. Neurosci Lett 475, 174 \(2010\).](#)  
[Helmond, Z. et al. Brain Patho 20, 787 \(2010\).](#)  
[Nagano, T. et al. Brain Res 1323, 11 \(2010\).](#)  
[Ravindran, C. et al. Cell Immunol 260, 113 \(2010\).](#)  
[Ryan, D. et al. J Neurosci Meth 190, 171 \(2010\).](#)  
[Spuch, C. et al. Neurotox Res 17, 421 \(2010\).](#)  
[Storage, D. et al. Arch Neurol 67, 867 \(2010\).](#)  
[Tian, G. et al. J Neurochem 113, 978 \(2010\).](#)  
[Tran, T. et al. Neurobiol Disease 37, 756 \(2010\).](#)  
[Tran, T. et al. Hippocampus 10.1002/hipo.20790 \(2010\).](#)  
[Vargas, T. et al. J Biomed Biotech 10.1155/2010/805405 \(2010\).](#)  
[Wang, Y. et al. J Phys Chem B 114, 7112 \(2010\).](#)  
[Wirths, O. et al. J Neural Transmission 117, 85 \(2010\).](#)  
[Arriagada, C. et al. Neuro Lett 470, 81 \(2009\).](#)  
[Giliberto, L. et al. J Biol Chem 284, 9027 \(2009\).](#)  
[Gu, Z. et al. J Biol Chem 284, 10639 \(2009\).](#)  
[Guix, F.X. et al. Brain 132, 1335 \(2009\).](#)  
[Li, M. et al. J Immunol 182, 5778 \(2009\). Zhao, L. et al. J Neurosci 29, 3603 \(2009\). Gevorkian, G. et al. Neurochem Int 52, 1030 \(2008\). Watanabe, T. et al. J Pharmacol Sci 106, 84 \(2008\).](#)  
[Chafekar, SM. et al. Antioxidants & Redox Signaling 9, 2245 \(2007\).](#)  
[Nelson, T.J. & DL Aklon, J Biol Chem 10.1074/jbc.M705558200 \(2007\).](#)  
[Bales, KR. et al. J Clin Invest 116, 825 \(2006\).](#)  
[Guo, J-P. et al. Proc Natl Acad Sci 103, 1953 \(2006\).](#)  
[Shaked, G. M. et al. FASEB J 20, 1254 \(2006\).](#)  
[Osada, Y. et al. J Biol Chem 280, 8596 \(2005\).](#)  
[Piccini, A. et al. J Biol Chem 280, 34186 \(2005\).](#)  
[Boyd-Kimball, D. et al. Chem Res Toxicol 17, 1743 \(2004\).](#)

- Legleiter, J. et al. *J Mol Bio* **335**, 997 (2004).  
Bard, F. et al. *PNAS* **100**, 2023 (2003).  
Takata, K. et al. *BBRC* **301**, 699 (2003).  
Uryu, S. et al. *BBRC* **303**, 302 (2003).  
Arispe, N. et al. *FASEB J* **16**, 1526 (2002).  
Kansi, J. et al. *Neurotox Res* **4**, 219 (2002).  
Barger, SW. et al. *J Neurochem* **76**, 846 (2001).  
Yao, Z. et al. *Brain Res* **889**, 181 (2001).  
Liang, JJN. *FEBS Lett* **484**, 98 (2000).  
Yatin, SM. et al. *J Alzheimer's Dis* **2**, 123 (2000).  
Yao, ZX. Et al. *Brain Res* **847**, 203 (1999).  
Yatin, SM. et al. *Neurosci Lett* **263**, 17 (1999).  
Bradt, BM. et al. *J Exp Med* **188**, 431 (1998).

*For Research Use Only*