

PRODUCT ANALYSIS SHEET

Catalog Number:	03-136
Lot Number:	????
Quantity:	1 mg
Description:	Major constituent of plaques and tangles that occur in Alzheimer's disease (AD) patients.
Sequence:	H ₂ N-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
Molecular Weight:	4329.9
Purity:	≥95% by HPLC analysis (purity based on peak area)
Amino Acid Analysis and Identity:	Confirms expected sequence
Peptide Content:	????? %
Physical Appearance:	Lyophilized powder
Solubility:	H ₂ O, TFA (trifluoroacetic acid), DMSO, HFIP (1,1,1,3,3,3-hexafluoro-2-propanol). Solubility in H ₂ O: 1 mg/mL.
Storage:	-20°C
Expiration Date:	See product label
References:	Borchelt, D.R., et al. (1997) <i>Neuron</i> 19:939-945. Arendash, G.W., et al. (1999) <i>Neuroscience Letters</i> 268:17-20. Yan, S.D., et al. (1999) <i>J. Biol. Chem.</i> 274:2145-2156. Bradt, B.M., et al. (1999) <i>J. Exp. Med.</i> 188:431-438. Ulery, P.G., et al. (2000) <i>J. Biol. Chem.</i> 275:7410-7415. Eisenhauer, P.B., et al. (2000) <i>J. Neurosci. Res.</i> 60(6):804-810. Lefterov, I.M., et al. (2000) <i>FASEB J.</i> 14:1837-1847. Town, T., et al. (2001) <i>Neuroscience Letters</i> 307:101-104. Viel, J.J., et al. (2001) <i>J. Neurosci. Res.</i> 64(5):454-465. Wang, S., et al. (2001) <i>J. Biol. Chem.</i> 276(45):42027-42034. Paris, D., et al. (1999) <i>Exp. Neurol.</i> 157:211-221. Paris, D., et al. (2002) <i>Atherosclerosis</i> 161(2):293-299. Paris, D., et al. (2002) <i>Prostaglandins and Other Lipid Mediators</i> 70(1-2):1-12. Wei, W.L., et al. (2002) <i>J. Biol. Chem.</i> 277(20):17649-17656. Tian, G.C., et al. (2002) <i>J. Biol. Chem.</i> 277(35):31499-31505. Kung, M.-P., et al. (2002) <i>Brain Research</i> 956:202-210. Agdeppa, E.D., et al. (2003) <i>Neuroscience</i> 117(3):723-730. Luo, X.G., et al. (2003) <i>J. Immunol.</i> 135(1-2):62-71.

For Research Use Only. CAUTION: Not for human or animal therapeutic or diagnostic use.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

Recommendations for Peptide Reconstitution:

Preparing peptide for neurotoxicity studies, to induce peptide aggregation:

The appearance of toxicity has recently been shown to correlate to the extent of beta sheet structure (S. Wang et al. [2001] *J. Biol. Chem.* **276**(45):42027-42034). Recommended preincubation is:

1. Dissolve the lyophilized peptide in 0.1% (v/v) trifluoroacetic acid in water at 10 mg/mL.
2. Dilute the peptide to 0.5–1.0 mg/mL with PBS (without Ca²⁺).
3. Incubate at 25°C for 24–48 h (24–36 h is usually sufficient).

Neurotoxic activity is usually observed at 30–100 µg/mL.

Preparing peptide for studies which require minimal peptide aggregation:

1. Dissolve the peptide at a concentration of 1 mg/mL in 100% HFIP (1,1,1,3,3,3-hexafluoro-2-propanol [Sigma-Aldrich Cat. # 32,524-4, 99.8% ACS reagent grade]).
2. Incubate at RT for 2 hours. During the incubation, vortex the peptide solution several times at moderate speed, allowing the HFIP to cover as much of the surface area as possible.
3. Dry down the HFIP/peptide solution under a gentle stream of nitrogen gas. Continue drying for an additional 10 minutes. Cap vial immediately.
4. Resuspend the peptide in 100% DMSO.
5. Incubate the peptide plus DMSO for 12 minutes at RT with periodic vortexing at moderate speed.
6. Add 50 µL of this DMSO/peptide solution dropwise to 10 mL of BSAT-DPBS (see formulation below) while vortexing at moderate speed.

Peptides prepared in this manner have been used as standards in ELISA for the detection of beta amyloid in biological samples. When using this peptide in ELISA, it is important that the assay buffer used with the peptide standards has the same composition as the samples under investigation.

Buffer Formulations:

DPBS Solution (10X Stock) (Biofluids Cat. # 316-500)

Dulbecco's PBS (DPBS w/o Mg²⁺, Ca²⁺)









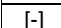
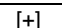



BSAT-DPBS Solution:

1X DPBS, pH 7.4

5% BSA

0.03% Tween-20

Explanation of symbols

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		In vitro diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

For Research Use Only. CAUTION: Not for human or animal therapeutic or diagnostic use.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com