



## Product Data Sheet

<b>Product Name:</b>	$\beta$ -Amyloid (1-42)-Lys(Biotin)-NH <sub>2</sub>	
<b>Catalog Number:</b>	61484-01 (0.1 mg) 61484-05 (0.5 mg)	Lot Number: See label on vial
<b>Sequence:</b>	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-Lys(Biotin)-NH <sub>2</sub> (3-letter code) DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA -K(Biotin)-NH <sub>2</sub> (1-letter code)	
<b>Molecular Weight:</b>	4868.6	
<b>% Peak Area by HPLC:</b>	≥ 95	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** Reconstitute by adding 50  $\mu$ l 1%NH<sub>4</sub>OH to 0.5 mg  $\beta$ -Amyloid 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:**  $\beta$ -Amyloid 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:** The difference between this recently released and the previous product (Cat# 23528-01) is the change at the C terminus from a carboxylic acid to an amide. The biological activity should remain the same. Elbaum, D. et al. *Biochem. Biophys. Res. Commun.* **267**, 733 (2000).

### Related Products:

Name	Cat #	Size
$\beta$ -Amyloid (1-42)-Lys(Biotin), FAM-labeled, Ex/Em=492/518 nm	23599-01	0.1 mg
FAM-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA-K(Biotin)	23598	0.5 mg
Biotin- $\beta$ -Amyloid (1-42)	23524-01	0.1 mg
Biotin-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA	23523-05	0.5 mg
Biotin-LC- $\beta$ -Amyloid (1-42)	24641-01	0.1 mg
Biotin-LC-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA	24640	0.5 mg
Biotin-LC- $\beta$ -Amyloid (1-42), mouse, rat	61718-01	0.1 mg
Biotin-LC-DAEFGHDSGFVRRHQKLVFFAEDVGSNKGAIIGLMVGGVVIA	61718-05	0.5 mg

*For Research Use Only*