



## Product Data Sheet

Product Name:	[Pro <sup>31</sup> ]- $\beta$ -Amyloid (1-42)	
Catalog Number:	62443 (0.5 mg)	Lot Number: See label on vial
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-Pro-Ile-Gly-Leu-Met-Val-Gly-Gly-Val-Val-Ile-Ala-OH (3-letter code) DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAPIGLMVGGVVIA (1-letter code)	
Molecular Weight:	4498.1	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

**Peptide Reconstitution:** Using basic buffer (1%NH<sub>4</sub>OH), reconstitute by adding 50  $\mu$ l to 0.5 mg [Pro<sup>31</sup>]- $\beta$ -Amyloid (1-42) peptide. Dilute this stock solution with working buffer, such as PBS or other buffers.

**Storage:** [Pro<sup>31</sup>]- $\beta$ -Amyloid (1-42) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and store at -20°C or lower.

**Description:** This  $\beta$ -Amyloid peptide (1-42) has Ile<sup>31</sup> substituted with Pro<sup>31</sup>. Substitution of Ile<sup>31</sup> by the helix-breaking amino acid, proline, completely abrogates the oxidative stress and neurotoxic properties of  $\beta$ -Amyloid (1-42). Ref: Butterfield, DA. *Free Rad. Res.* **36**, 1307 (2002)

### Related Products:

Name	Cat #	Size
$\beta$ -Amyloid (1-42)	24224	0.5 mg
DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAPIGLMVGGVVIA	20276	1 mg
	20276-5	5 mg
	20276-25	25 mg
	20276-100	100 mg

*For Research Use Only*