



## Product Data Sheet

---

**Product Name:** [Cys<sup>26</sup>] - $\beta$ -Amyloid (1-42), S26C  $\beta$ -Amyloid (1-42)

**Catalog Number:** AS-63672-05 (0.5 mg)                      Lot Number: See label on vial  
AS-63672-1 (1 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-Cys-Asn-Lys-Gly-Ala-Ile-Ile-Gly-  
Leu-Met-Val-Gly-Gly-Val-Val-Ile-Ala-OH (3-letter code)  
DAEFRHDSGYEVHHQKLVFFAEDVGCNKGAIIGLMVGGVVIA  
(1-letter code)

**Molecular Weight:** 4531.2

**% Peak Area by HPLC:**  $\geq 95$

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:** Using the DMSO, reconstitute by adding 50  $\mu$ l to 0.5 mg S26C  $\beta$ -Amyloid peptide. This peptide is also soluble in 1% NH<sub>4</sub>OH, but not as clear as DMSO.

**Storage:** S26C  $\beta$ -Amyloid peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Related Products:**

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
(DAEFRHDSGYEVHHQKLVFFAEDVGCNKGAIIGLMVGGVV)	AS-63671-1	1 mg
$\beta$ -Amyloid (1-42)		
(DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA)	AS-24224	0.5 mg
	AS-20276	1 mg

*For Research Use Only*