

| Product Name: | [Cys ²⁶] - β -Amyloid (1-42), S26C β -Amyloid (1-42) | | | |
|---------------------------|---|-------------------------------|--|--|
| Catalog Number: | AS-63672-05 (0.5 mg) AS-63672-1 (1 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-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: ≥ 95 | | | | |
| Appearance: | Lyophilized white powder | | | |

Peptide Reconstitution: Using the DMSO, reconstitute by adding 50 µl to 0.5 mg S26C β-Amyloid peptide. This peptide is also soluble in 1%NH₄OH, but not as clear as DMSO.

Storage: S26C β -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.

| Related Products: Name | Cat # | Size |
|--|----------------------|----------------|
| (DAEFRHDSGYEVHHQKLVFFAEDVGCNKGAIIGLMVGGVV) | AS-63671-1 | 1 mg |
| β-Amyloid (1-42) (DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA) | AS-24224 AS-20276 | 0.5 mg 1 mg |

For Research Use Only