



## Product Data Sheet

---

**Product Name:**  $\beta$ -Amyloid (17-40)

**Catalog Number:** AS-22813 (0.5 mg)                      Lot Number: See label on vial  
AS-22814 (1 mg)

**Sequence:** H-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 (3-letter code)  
LVFFAEDVGSNKGAIIGLMVGGVV (1-letter code)

**Molecular Weight:** 2392.9

**Peptide Purity:** >95%

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:**  $\beta$ -Amyloid (17-40) peptide is freely soluble in 1% NH<sub>4</sub>OH.

**Storage:**  $\beta$ -Amyloid (17-40) 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.

### Published Citations:

Morishima-Kawashima, M. et al. *Am. J. Pathol.* **157**, 2093 (2000).  
Chauhan, A. et al. *Neurochem. Res.* **25**, 423 (2004).  
Minicozzi, V. et al. *J. Biol. Chem.* 10.1074/jbc.M707109200 (2008).

### Related Products:

| Name                            | Cat #    | Size |
|---------------------------------|----------|------|
| Biotin-LVFFAEDVGSNKGAIIGLMVGGVV | AS-23541 | 1 mg |

*For Research Use Only*