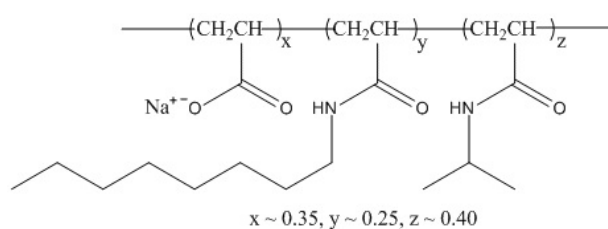




## Amphipol A8-35

Cat. No.	Amount
X-A835	50 mg



Amphipol A8-35 Structure

**For general laboratory use.****Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months**Molecular Formula:**  $(C_{6.2}H_{10.3}O_{1.35}N_{0.65}Na_{0.35})_{72}$ **Molecular Weight:** approx. 9 kDa**CAS#:** 326856-53-5**Applications:**Stabilizing agent in Cryo-EM<sup>[2-5]</sup> and X-ray crystallography<sup>[6]</sup>**Description:**

Amphipol A8-35 is a short amphipatic polymer that is specifically designed for membrane protein stabilization. The surfactant possesses a very high affinity for the transmembrane surfaces and allows to solubilize membrane proteins in a detergent-free aqueous solution<sup>[1]</sup>.

**Selected References:**

- [1] Zoonens *et al.* (2014) Amphipols for Each Season. *J Membrane Biol* **247**:759.
- [2] Chen *et al.* (2016) Structure of the STRA6 receptor for retinol uptake. *Science* **353**:887.
- [3] Zubcevic *et al.* (2016) Cryo-Electron Microscopy of the Trpv2 Ion Channel. *Nat Struct Mol Biol* **23**:180.
- [4] Bai *et al.* (2015) Sampling the conformational space of the catalytic subunit of human gamma-secretase. DOI 10.7554/eLife.11182.
- [5] Althoff *et al.* (2011) Arrangement of electron transport chain components in bovine mitochondrial supercomplex I<sub>1</sub>III<sub>2</sub>IV<sub>1</sub>. *EMBO J* **30**:4652.
- [6] Polovinkin *et al.* (2014) High-Resolution Structure of a Membrane Protein Transferred from Amphipol to a Lipidic Mesophase. *J Membrane Biol* **247**:997.