



## Product Data Sheet

---

**Product Name:**  $\beta$ -Secretase Inhibitor 1

**Catalog Number:** AS-23958 (1 mg) **Lot Number:** See label on vial

**Sequence:** H-Lys-Thr-Glu-Glu-Ile-Ser-Glu-Val-Asn-Sta-Val-Ala-Glu-Phe-OH  
(3-letter code)  
KTEEISEVN-Sta-VAEF (Sta=statine) (1-letter code)

**Molecular Weight:** 1651.9

**Peptide Purity:** >95%

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:**  $\beta$ -Secretase peptide is freely soluble in basic buffer.

**Storage:**  $\beta$ -Secretase 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.

**Description:** This statine-based substrate analog,  $\beta$ -secretase inhibitor P10–P4' statV, is used in the inhibition of  $\beta$ -secretase activity from homogenized wild-type mouse cortices and from BACE-purified from human brain. Ref: Sinha, S. et al. *Nature* **402**, 537 (1999).

*For Research Use Only*