



## Product Data Sheet

Product Name:	$\beta$ -Secretase Inhibitor 1	
Catalog Number:	23958 (1 mg) 23959 (5 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).

### Related Products:

Name	Cat #	Size
$\beta$ -Secretase Inhibitor 2	23960	1 mg
KTEEISEVN-Sta-VAEF-NH <sub>2</sub> (Sta=Statine)	23961	5 mg
$\beta$ -Secretase Substrate 1a Mca-EVKVDAEF-K(Dnp)	60268	1 mg
$\beta$ -Secretase Substrate 2 Mca-EVKMDAEFK(Dnp)	60269	1 mg

*For Research Use Only*