

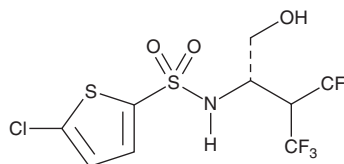
PRODUCT INFORMATION



Begacestat

Item No. 27916

CAS Registry No.: 769169-27-9
Formal Name: 5-chloro-N-[(1S)-3,3,3-trifluoro-1-(hydroxymethyl)-2-(trifluoromethyl)propyl]-2-thiophenesulfonamide
Synonym: GSI-953
MF: C₉H₈ClF₆NO₃S₂
FW: 391.7
Purity: ≥98%
UV/Vis.: λ_{max}: 262 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Begacestat is an inhibitor of γ -secretase that inhibits cleavage of amyloid precursor protein (APP) to amyloid- β (1-40) (A β 40; Item No. 21617) *in vitro* (IC₅₀ = 15 nM).¹ It is 14-fold selective for γ -secretase over Notch-1. Begacestat (5 mg/kg) reduces brain levels of A β 40 and A β 42 (Item No. 20574) in the Tg2576 transgenic mouse model of Alzheimer's disease. It also increases contextual fear conditioning freezing time, indicating reversal of contextual memory deficits, in Tg2576 mice in a dose-dependent manner.²

References

1. Mayer, S.C., Kreft, A.F., Harrison, B., *et al.* Discovery of begacestat, a Notch-1-sparing γ -secretase inhibitor for the treatment of Alzheimer's disease. *J. Med. Chem.* **51**(23), 7348-7351 (2008).
2. Martone, R.L., Zhou, H., Atchison, K., *et al.* Begacestat (GSI-953): A novel, selective thiophene sulfonamide inhibitor of amyloid precursor protein γ -secretase for the treatment of Alzheimer's disease. *J. Pharmacol. Exp. Ther.* **331**(2), 598-608 (2009).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/01/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM