

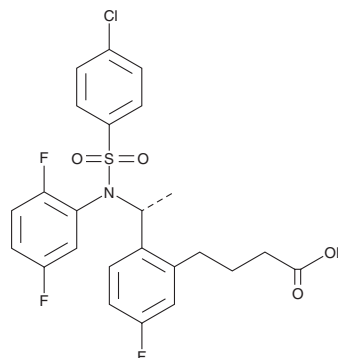
PRODUCT INFORMATION



BMS-299897

Item No. 35777

CAS Registry No.: 290315-45-6
Formal Name: 2-[(1R)-1-[[4-(4-chlorophenyl)sulfonyl](2,5-difluorophenyl)amino]ethyl]-5-fluoro-benzenebutanoic acid
MF: C₂₄H₂₁ClF₃NO₄S
FW: 511.9
Purity: ≥98%
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.

Laboratory Procedures

BMS-299897 is supplied as a solid. A stock solution may be made by dissolving the BMS-299897 in the solvent of choice, which should be purged with an inert gas. BMS-299897 is soluble in methanol.

Description

BMS-299897 is an inhibitor of γ -secretase.¹⁻³ It selectively cleaves the carboxy terminal fragment (CTF) of amyloid precursor protein (APP) over the Notch-1 CTF in HEK293 cells (IC₅₀s = 7.1 and 105.9 nM, respectively).² BMS-299897 (0.1-1 nmol/animal) reduces increases in amyloid- β (1-42) (A β 42) levels induced by A β (25-35) in the mouse hippocampus.³ It also prevents deficits induced by A β (25-35) in the novel object recognition test when tested with a one-hour, but not a 24-hour, intertrial interval in the same model.

References

1. Tian, G., Sobotka-Briner, C.D., Zysk, J., *et al.* Linear non-competitive inhibition of solubilized human gamma-secretase by pepstatin A methylester, L685458, sulfonamides, and benzodiazepines. *J. Biol. Chem.* **277**(35), 31499-31505 (2002).
2. Barten, D.M., Guss, V.L., Corsa, J.A., *et al.* Dynamics of β -amyloid reductions in brain, cerebrospinal fluid, and plasma of β -amyloid precursor protein transgenic mice treated with a γ -secretase inhibitor. *J. Pharmacol. Exp. Ther.* **312**(2), 635-643 (2005).
3. Meunier, J., Villard, V., Givalois, L., *et al.* The γ -secretase inhibitor 2-[(1R)-1-[[4-(4-chlorophenyl)sulfonyl](2,5-difluorophenyl) amino]ethyl]-5-fluorobenzenebutanoic acid (BMS-299897) alleviates A β ₁₋₄₂ seeding and short-term memory deficits in the A β ₂₅₋₃₅ mouse model of Alzheimer's disease. *Eur. J. Pharmacol.* **698**(1-3), 193-199 (2013).

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, 12/13/2022

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