

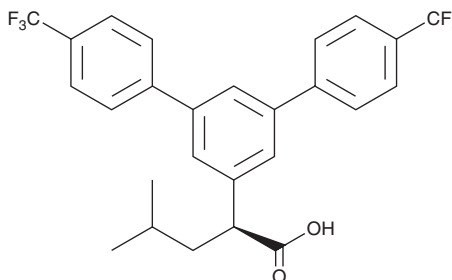
PRODUCT INFORMATION



JNJ-40418677

Item No. 21869

CAS Registry No.: 1146594-87-7
Formal Name: α S-(2-methylpropyl)-4,4''-bis(trifluoromethyl)-[1,1':3',1''-terphenyl]-5'-acetic acid
MF: C₂₆H₂₂F₆O₂
FW: 480.4
Purity: ≥98%
UV/Vis.: λ_{max} : 254 nm
Supplied as: A crystalline 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

JNJ-40418677 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-40418677 in the solvent of choice, which should be purged with an inert gas. JNJ-40418677 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JNJ-40418677 in DMF is approximately 25 mg/ml and approximately 10 mg/ml in ethanol and DMSO.

Description

JNJ-40418677 is a modulator of γ -secretase.¹ It reduces the levels of amyloid- β (1-42) (A β 42; Item No. 20574) in a cell-free γ -secretase modulation assay. It also reduces the secretion of A β 42, but not total A β , in SK-N-BE(2) human neuroblastoma cells expressing amyloid precursor protein (APP; mean IC₅₀ = 200 nM) and in primary rat cortical neurons (mean IC₅₀ = 185 nM). JNJ-40418677 (100 and 300 mg/kg) reduces the level of A β 42 in wild-type mouse brain. It decreases brain levels of A β 42, A β 40, and A β 38 in the deposited fraction, and A β 42 and A β 40 in the soluble fraction, in the Tg2576 mouse model of Alzheimer's disease when administered at doses of 60 and 120 mg/kg per day in the diet for seven months. It also reduces increases in the number of amyloid plaques in Tg2576 mouse brain compared to vehicle control animals.

Reference

1. Van Broeck, B., Chen, J.-M., Tréton, G., *et al.* Chronic treatment with a novel γ -secretase modulator, JNJ-40418677, inhibits amyloid plaque formation in a mouse model of Alzheimer's disease. *Br. J. Pharmacol.* **163**(2), 375-389 (2011).

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, 11/09/2022

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