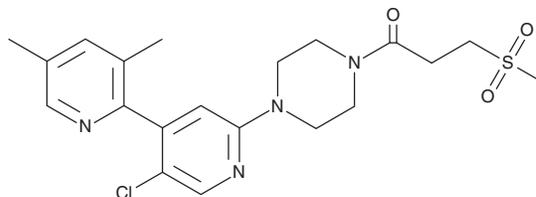


PRODUCT INFORMATION



PF-5274857
Item No. 43595

CAS Registry No.: 1373615-35-0
Formal Name: 1-[4-(5'-chloro-3,5-dimethyl[2,4'-bipyridin]-2'-yl)-1-piperazinyl]-3-(methylsulfonyl)-1-propanone
MF: C₂₀H₂₅ClN₄O₃S
FW: 437.0
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

PF-5274857 is supplied as a solid. A stock solution may be made by dissolving the PF-5274857 in the solvent of choice, which should be purged with an inert gas. PF-5274857 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

PF-5274857 is a smoothed receptor (Smo) antagonist.¹ It binds to Smo (K_i = 4.6 nM) and inhibits transcription of *Gli1* in mouse embryonic fibroblasts (MEFs; IC₅₀ = 2.7 nM). *In vivo*, PF-5274857 inhibits tumor growth and induces tumor regression in a medulloblastoma mouse allograft model when administered at doses less than 10 mg/kg and greater than 10 mg/kg, respectively.

Reference

1. Rohner, A., Spilker, M.E., Lam, J.L., *et al.* Effective targeting of hedgehog signaling in a medulloblastoma model with PF-5274857, a potent and selective smoothed antagonist that penetrates the blood-brain barrier. *Mol. Cancer Ther.* **11(1)**, 57-65 (2012).

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, 07/31/2025

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