

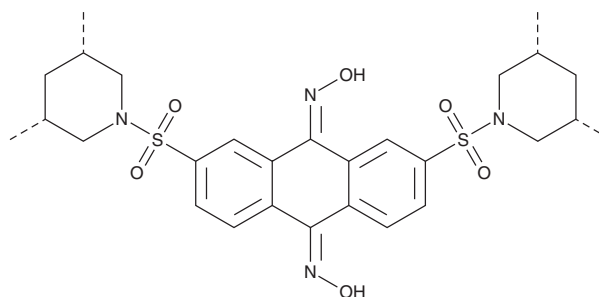
PRODUCT INFORMATION



Tegatrabetan

Item No. 37423

CAS Registry No.: 1227637-23-1
Formal Name: *rel*-2,7-bis[[[(3R,5S)-3,5-dimethyl-1-piperidinyl] sulfonyl]-9,10-anthracenedione 9,10-dioxime
Synonym: BC-2059, Tegavivint
MF: C₂₈H₃₆N₄O₆S₂
FW: 588.7
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 274, 304 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.

Laboratory Procedures

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

Description

Tegatrabetan is an inhibitor of Wnt/β-catenin signaling.¹ It reduces Wnt3a-induced transcriptional activity in a reporter activity assay using STF3a cells, which exhibit constitutive Wnt activation, when used at a concentration of 1 μM. Tegatrabetan inhibits proliferation of, and transcription in, HT-29 human colon cancer cells (IC₅₀s = 9 and 63 nM, respectively) and reduces survival in six other cancer cell lines (IC₅₀s = 14-80 nM). It induces cell cycle arrest at the G₁ phase in HCT-15 human colon cancer cells when used at a concentration of 625 nM. *In vivo*, tegatrabetan (50 mg/kg per day) reduces tumor volume and inhibits lung metastasis in a Saos-2 osteosarcoma mouse xenograft model.²

References

1. Soldi, R., Horrigan, S.K., Cholody, M.W., *et al.* Design, synthesis, and biological evaluation of a series of anthracene-9,10-dione dioxime β-catenin pathway inhibitors. *J. Med. Chem.* **58(15)**, 5854-5862 (2015).
2. Nomura, M., Rainusso, N., Lee, Y.-C., *et al.* Tegavivint and the β-catenin/ALDH axis in chemotherapy-resistant and metastatic osteosarcoma. *J. Natl. Cancer Inst.* **111(11)**, 1216-1227 (2019).

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, 08/19/2025

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