

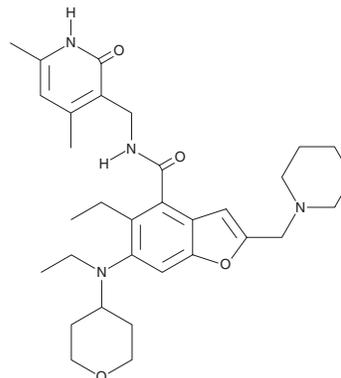
PRODUCT INFORMATION



Zeprumetostat

Item No. 44976

CAS Registry No.: 2098545-98-1
Formal Name: N-[(1,2-dihydro-4,6-dimethyl-2-oxo-3-pyridinyl)methyl]-5-ethyl-6-[ethyl(tetrahydro-2H-pyran-4-yl)amino]-2-(1-piperidinylmethyl)-4-benzofurancarboxamide
Synonym: SHR2554
MF: C₃₂H₄₄N₄O₄
FW: 548.7
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

Zeprumetostat is supplied as a solid. A stock solution may be made by dissolving the zeprumetostat in the solvent of choice, which should be purged with an inert gas. Zeprumetostat is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Zeprumetostat is an inhibitor of enhancer of zeste homolog 2 (EZH2; IC₅₀ = 0.87 nM for the wild-type enzyme).¹ It is selective for EZH2 over a panel of 24 histone and DNA methyltransferases (IC₅₀s = >10,000 nM for all) but also inhibits EZH1 (IC₅₀ = 19.1 nM). Zeprumetostat also inhibits the EZH2 mutants EZH^{2Y641C}, EZH^{2Y641F}, EZH^{2Y641N}, EZH^{2Y641S}, and EZH^{2Y641G} (IC₅₀s = 16.8, 2.68, 1.79, 5.07, and 1.13 nM, respectively). It reduces the levels of trimethylated lysine 27 on histone H3 (H3K27Me3) in Pfeiffer diffuse large B cell lymphoma (DLBCL) cells (IC₅₀ = 1.63 nM) and induces cell cycle arrest at the G₁ phase and apoptosis in the same cells. Zeprumetostat (60 and 120 mg/kg), in combination with HBI 8000 (chidamide; Item No. 13686), reduces tumor growth in patient-derived xenograft (PDX) mouse models of DLBCL.

Reference

1. Wang, X., Wang, D., Ding, N., *et al.* The synergistic anti-tumor activity of EZH2 inhibitor SHR2554 and HDAC inhibitor chidamide through ORC1 reduction of DNA replication process in diffuse large B cell lymphoma. *Cancers (Basel)* **13(17)**, 4249 (2021).

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, 02/06/2026

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