

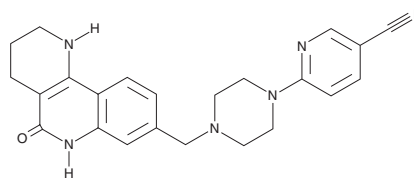
PRODUCT INFORMATION



Nesuparib

Item No. 45304

CAS Registry No.: 2055357-64-5
Formal Name: 6-[4-[(1,2,3,4,5,6-hexahydro-5-oxobenzo[h]-1,6-naphthyridin-8-yl)methyl]-1-piperazinyl]-3-pyridinecarbonitrile
Synonym: JPI-547
MF: C₂₃H₂₄N₆O
FW: 400.5
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

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

Description

Nesuparib is a dual inhibitor of the poly(ADP-ribose) polymerases PARP1/2 and tankyrases TNKS1/2.¹ It reduces the proliferation of Capan-1 pancreatic ductal adenocarcinoma (PDAC) cells with an IC₅₀ value of 0.1895 μM. Nesuparib (0.01-1 μM) selectively inhibits the proliferation of, and induces apoptosis and G₂/M phase cell cycle arrest in, HPAF-II and AsPC-1 PDAC cells expressing RNF43 loss-of-function mutations over SNU-410 PDAC cells expressing wild-type RNF43. Nesuparib (50 mg/kg) also reduces tumor growth in a Capan-1 mouse xenograft model

Reference

1. Oh, K.S., Nam, A.R., Bang, J.H., *et al.* JPI-547, a dual inhibitor of PARP/Tankyrase, shows antitumor activity against pancreatic cancers with homologous recombination repair deficiency or Wnt-addiction. *Int. J. Biol.* **21(12)**, 5460-5475 (2025).

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, 03/10/2026

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