

PRODUCT INFORMATION



YCH1899

Item No. 39737

Formal Name: 4-(3-(3-(5-bromofuran-2-yl)-5,6,7,8-tetrahydro-[1,2,4]triazolo[4,3-a]pyrazine-7-carbonyl)-4-fluorobenzyl)phthalazin-1(2H)-one

MF: C₂₅H₁₈BrFN₆O₃

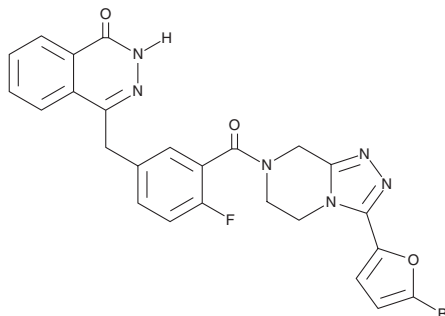
FW: 549.4

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥3 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

YCH1899 is supplied as a solid. A stock solution may be made by dissolving the YCH1899 in the solvent of choice, which should be purged with an inert gas. YCH1899 is soluble in DMSO.

Description

YCH1899 is an inhibitor of poly(ADP-ribose) polymerase 1 (PARP1) and PARP2 (IC₅₀s = <0.001 nM for both).¹ It is selective for PARP1 and PARP2 over PARP3, -4, -5A, -5B, -6, -7, -10, and -12 (IC₅₀s = 1-14.1 nM). YCH1899 inhibits the proliferation of Capan-1 cells (IC₅₀ = 0.1 nM), as well as Capan-1 cells resistant to the PARP inhibitors talazoparib (BMN 673; Item No. 19782) or olaparib (Item No. 10621; IC₅₀s = 1.13 and 0.89 nM, respectively). It inhibits the proliferation of BRCA1 mutant HCC1937 breast cancer cells (IC₅₀ = 4.54 nM). *In vivo*, YCH1899 (12.5 and 25 mg/kg) reduces tumor growth in an olaparib-resistant MDA-MB-436 breast cancer mouse xenograft model.

Reference

1. Sun, Y., Yang, H., Yuan, J., *et al.* YCH1899, a highly effective phthalazin-1(2H)-one derivative that overcomes resistance to prior PARP inhibitors. *J. Med. Chem.* **66**(17), 12284-12303 (2023).

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, 11/13/2023

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