

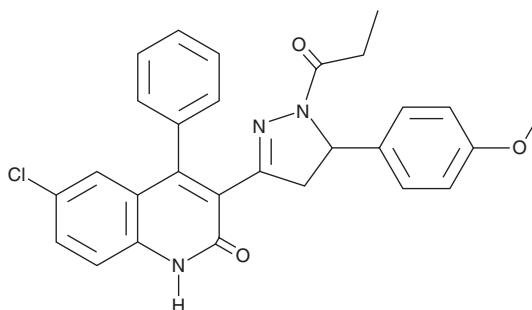
PRODUCT INFORMATION



CAY10760

Item No. 30532

CAS Registry No.: 391889-85-3
Formal Name: 6-chloro-3-[4,5-dihydro-5-(4-methoxyphenyl)-1-(1-oxopropyl)-1H-pyrazol-3-yl]-4-phenyl-2(1H)-quinolinone
MF: C₂₈H₂₄ClN₃O₃
FW: 486.0
Purity: ≥98%
UV/Vis.: λ_{max}: 234 nm
Supplied as: A crystalline 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

CAY10760 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10760 in the solvent of choice, which should be purged with an inert gas. CAY10760 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of CAY10760 in these solvents is approximately 10 and 11 mg/ml, respectively.

CAY10760 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10760 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10760 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CAY10760 is an inhibitor of the protein-protein interaction between RAD51 recombinase and BRCA2 (EC₅₀ = 19 μM in a cell-free competitive ELISA).¹ It decreases homologous recombination by 54% in wild-type BRCA2-expressing BxPC-3 pancreatic cancer cells when used at a concentration of 20 μM. CAY10760 (20 μM) decreases proliferation of BxPC-3, as well as mutant BRCA2-expressing Capan-1, cancer cells when used alone or in combination with the poly(ADP-ribose) polymerase (PARP) inhibitor olaparib (Item No. 10621).

Reference

1. Bagnolini, G., Milano, D., Manerba, M., *et al.* Synthetic lethality in pancreatic cancer: Discovery of a new RAD51-BRCA2 small molecule disruptor that inhibits homologous recombination and synergizes with olaparib. *J. Med. Chem.* **63**(5), 2588-2619 (2020).

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, 03/21/2024

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