

PRODUCT INFORMATION



PDD00017273

Item No. 19511

CAS Registry No.: 1945950-21-9
Formal Name: 1-[(1,3-dimethyl-1H-pyrazol-5-yl)methyl]-1,2,3,4-tetrahydro-N-(1-methylcyclopropyl)-3-[(2-methyl-5-thiazolyl)methyl]-2,4-dioxo-6-quinazolinesulfonamide

MF: C₂₃H₂₆N₆O₄S₂

FW: 514.6

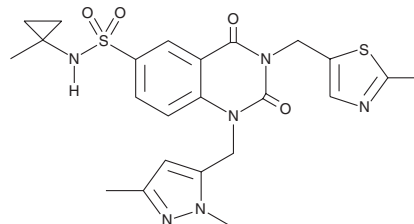
Purity: ≥98%

UV/Vis.: λ_{max}: 227 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

PDD00017273 is supplied as a solid. A stock solution may be made by dissolving the PDD00017273 in the solvent of choice, which should be purged with an inert gas. PDD00017273 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

PDD00017273 is a poly(ADP-ribose) glycohydrolase (PARG) inhibitor (IC₅₀ = 26 nM in an enzyme assay).¹ It maintains persistence of nuclear PAR chains in HeLa cells induced by the DNA damaging agent methylmethanesulfonate (MMS; IC₅₀ = 37 nM). PDD00017273, in combination with MMS, increases nuclear phosphorylated H2AX (γH2AX) intensity, a marker of DNA double strand breaks, in MCF-7 cells in a concentration-dependent manner. It reduces clonogenic growth of ZR-75-1 and MDA-MB-436 cells but not HCC1937 cells (IC₅₀s = 0.2, 0.8, and >10 μM, respectively).

Reference

1. James, D.I., Smith, K.M., Jordan, A.M., *et al.* First-in-class chemical probes against poly(ADP-ribose) glycohydrolase (PARG) inhibit DNA repair with differential pharmacology to olaparib. *ACS Chem Biol.* **11**(11), 3179-3190 (2016).

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, 12/22/2022

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