

PRODUCT INFORMATION



YU238259

Item No. 25452

CAS Registry No.: 1943733-16-1

Formal Name: N-[2-(5-chloro-2-pyridinyl)ethyl]-4-[[[(4-methoxyphenyl)sulfonyl]amino]methyl]-benzamide

MF: C₂₂H₂₂ClN₃O₄S

FW: 459.9

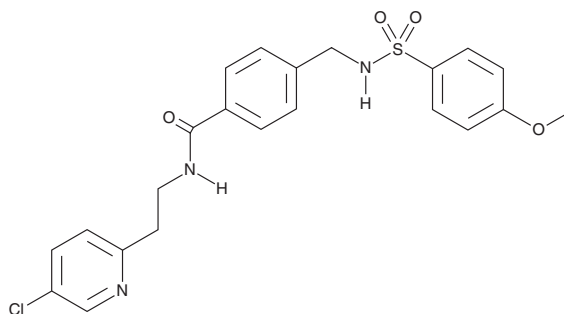
Purity: ≥95%

UV/Vis.: λ_{max}: 239 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

YU238259 is supplied as a crystalline solid. A stock solution may be made by dissolving the YU238259 in the solvent of choice, which should be purged with an inert gas. YU238259 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of YU238259 in these solvents is approximately 30 mg/ml.

YU238259 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, YU238259 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. YU238259 has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

YU238259 is an inhibitor of homology-dependent DNA repair (HDR) that is lethal to *BRCA2*^{-/-} but not *BRCA2*^{+/+} PEO1/4 cells (LD₅₀s = 8.7 and >100 μM, respectively).¹ YU238259 selectively inhibits HDR over non-homologous end joining (NHEJ), decreasing the number of U2OS cells with ionizing radiation-induced *BRCA1* foci but not 53BP1 or pDNA-PK foci. It also increases the number of ionizing radiation-induced double strand breaks (DSBs) in DLD-1 cells. YU238259 (1-20 μM) is cytotoxic to *BRCA2*^{-/-} DLD-1, *ATM*^{-/-} GM05849, and *PTEN*^{-/-} U251 cells but has no effect on wild-type cells. *In vivo*, YU238259 (3 mg/kg) delays tumor growth and increases survival in a *BRCA2*^{-/-}, but not a *BRCA*^{+/+}, DLD-1 mouse xenograft model.

Reference

1. Stachelek, G.C., Peterson-Roth, E., Liu, Y., *et al.* YU238259 is a novel inhibitor of homology-dependent DNA repair that exhibits synthetic lethality and radiosensitization in repair-deficient tumors. *Mol. Cancer Res.* **13**(10), 1389-1397 (2015).

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, 11/23/2022

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