

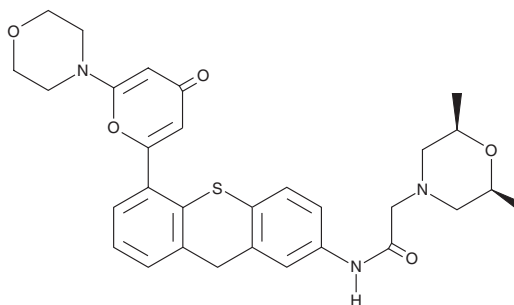
PRODUCT INFORMATION



Ku-60019

Item No. 17502

CAS Registry No.: 925701-46-8
Formal Name: *rel*-2*R*,6*S*-dimethyl-N-[5-[6-(4-morpholinyl)-4-oxo-4*H*-pyran-2-yl]-9*H*-thioxanthen-2-yl]-4-morpholineacetamide
MF: 547.7
FW: C₃₀H₃₃N₃O₅S
Purity: ≥98%
UV/Vis.: λ_{max}: 244, 272 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

Ku-60019 is supplied as a crystalline solid. A stock solution may be made by dissolving the Ku-60019 in the solvent of choice, which should be purged with an inert gas. Ku-60019 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Ku-60019 in ethanol and DMSO is approximately 50 mg/ml and approximately 20 mg/ml in ethanol and DMF.

Ku-60019 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Ku-60019 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Ku-60019 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ataxia-telangiectasia mutated (ATM) is a serine/threonine kinase that activates checkpoint signaling following double strand DNA breaks and genotoxic stress. Ku-60019 is a potent, reversible inhibitor of ATM kinase (IC₅₀ = 6.3 nM), blocking the phosphorylation of ATM substrate proteins.^{1,2} It is much less effective or without effect against a panel of 229 other kinases.¹ Ku-60019 sensitizes glioma cells to radiation and inhibits migration and invasion of glioma cells *in vitro*.^{1,2} It produces radiosensitization and increases survival *in vivo* when administered intra-tumorally in orthotopic xenograft models of glioblastoma multiforme.³ Ku-60019 is particularly effective in producing lethality in cells with mutant p53 or that are deficient in PTEN.^{3,4}

References

1. Golding, S.E., Rosenberg, E., Valerie, N., *et al.* Improved ATM kinase inhibitor KU-60019 radiosensitizes glioma cells, compromises insulin, AKT and ERK prosurvival signaling, and inhibits migration and invasion. *Mol. Cancer Ther.* **8**(10), 2894-2902 (2009).
2. Golding, S.E., Rosenberg, E., Adams, B.R., *et al.* Dynamic inhibition of ATM kinase provides a strategy for glioblastoma multiforme radiosensitization and growth control. *Cell Cycle* **11**(6), 1167-1173 (2012).
3. Biddlestone-Thorpe, L., Sajjad, M., Rosenberg, E., *et al.* ATM kinase inhibition preferentially sensitizes p53-mutant glioma to ionizing radiation. *Clin. Cancer Res.* **19**(12), 3189-3200 (2013).
4. McCabe, N., Hanna, C.D., Walker, S.M., *et al.* Mechanistic rationale to target PTEN-deficient tumour cells with inhibitors of the DNA damage response kinase ATM. *Cancer Res.* **75**(11), 2159-2165 (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, 10/24/2022

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