

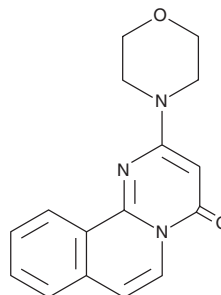
PRODUCT INFORMATION



Compound 401

Item No. 21769

CAS Registry No.: 168425-64-7
Formal Name: 2-(4-morpholinyl)-4H-pyrimido[2,1-a]isoquinolin-4-one
MF: C₁₆H₁₅N₃O₂
FW: 281.3
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 252, 276 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

Compound 401 is supplied as a crystalline solid. A stock solution may be made by dissolving the compound 401 in the solvent of choice, which should be purged with an inert gas. Compound 401 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of compound 401 in ethanol is approximately 0.5 mg/ml and approximately 1 mg/ml in DMSO and DMF.

Description

Compound 401 is an inhibitor of DNA-dependent protein kinase (DNA-PK) and mammalian target of rapamycin (mTOR; IC₅₀s = 0.28 and 5.3 μM, respectively).¹ It is selective for DNA-PK and mTOR over PI3K, ATM, and ATR (IC₅₀s = >100 μM for all). Compound 401 (10 μM) inhibits phosphorylation of the mTOR targets S6K1 and Akt in Rat-1 fibroblasts and in M059J glioma cells that lack DNA-PK.² It inhibits proliferation of mouse embryonic fibroblasts (MEFs) lacking tuberous sclerosis complex 1 (TSC1^{-/-}; IC₅₀ = 2 μM), a complex associated with hamartomas that display hyperactive mTOR signaling, but not TSC1^{+/+} MEFs. Compound 401 also induces apoptosis in TSC1^{-/-} MEFs.

References

1. Griffin, R.J., Fontana, G., Golding, B.T., *et al.* Selective benzopyranone and pyrimido[2,1-a]isoquinolin-4-one inhibitors of DNA-dependent protein kinase: Synthesis, structure-activity studies, and radiosensitization of a human tumor cell line in vitro. *J. Med. Chem.* **48**(2), 569-585 (2005).
2. Ballou, L.M., Selinger, E.S., Choi, J.Y., *et al.* Inhibition of mammalian target of rapamycin signaling by 2-(morpholin-1-yl)pyrimido[2,1-a]isoquinolin-4-one. *J. Biol. Chem.* **282**(33), 24463-24470 (2007).

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, 11/28/2022

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