

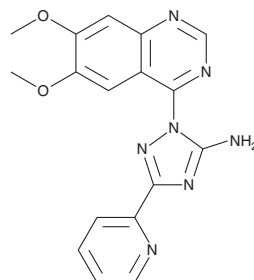
PRODUCT INFORMATION



CP 466,722

Item No. 25417

CAS Registry No.: 1080622-86-1
Formal Name: 1-(6,7-dimethoxy-4-quinazoliny)-3-(2-pyridiny)-1H-1,2,4-triazol-5-amine
MF: C₁₇H₁₅N₇O₂
FW: 349.4
Purity: ≥98%
UV/Vis.: λ_{max}: 243, 364 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

CP 466,722 is supplied as a crystalline solid. A stock solution may be made by dissolving the CP 466,722 in the solvent of choice, which should be purged with an inert gas. CP 466,722 is soluble in the organic solvent chloroform at a concentration of approximately 2 mg/ml.

Description

CP 466,722 is an inhibitor of ataxia-telangiectasia mutated (ATM) kinase that inhibits ionizing radiation-induced ATM autophosphorylation and phosphorylation of the ATM targets SMC1, Chk2, and p53 in HeLa cells when used at a concentration of 10 μM.¹ It is selective for ATM over PI3K and PIKK in HFF(hTERT) and A-T(hTERT) fibroblasts, respectively. CP 466,722 decreases cell survival in response to ionizing radiation in HeLa cells. It is also cytotoxic to MCF-7 and SKBr-3 breast cancer cells (IC₅₀s = 16.92 and 12.78 μM, respectively).¹

Reference

1. Rainey, M.D., Charlton, M.E., Stanton, R.V., *et al.* Transient inhibition of ATM kinase is sufficient to enhance cellular sensitivity to ionizing radiation. *Cancer Res.* **68(18)**, 7466-7474 (2008).

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, 11/11/2022

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