

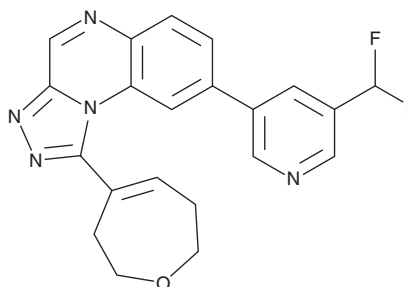
PRODUCT INFORMATION



BAY-8400

Item No. 43707

CAS Registry No.: 2763602-59-9
Formal Name: 8-[5-(difluoromethyl)-3-pyridinyl]-1-(2,3,6,7-tetrahydro-4-oxepinyl)-[1,2,4]triazolo[4,3-a]quinoxaline
MF: C₂₁H₁₇F₂N₅O
FW: 393.4
Purity: ≥98%
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

BAY-8400 is supplied as a solid. A stock solution may be made by dissolving the BAY-8400 in the solvent of choice, which should be purged with an inert gas. BAY-8400 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

BAY-8400 is an inhibitor of DNA-dependent protein kinase (DNA-PK; IC₅₀ = 0.081 μM).¹ It is selective for DNA-PK over ataxia-telangiectasia mutated kinase (ATM) and mTOR (IC₅₀s = 19.3 and 1.91 μM, respectively) and a panel of 365 additional kinases at 1 μM but does inhibit ataxia-telangiectasia and Rad3-related protein/kinase (ATR) and PI3Kβ (IC₅₀s = 0.394 and 0.117 μM, respectively). BAY-8400 inhibits the formation of γH2AX in ATM-negative HT-144 melanoma cells (IC₅₀ = 0.069 nM) and reduces the proliferation of HT-29 colon cancer cells when used alone or in combination with bleomycin. *In vivo*, BAY-8400 (150 mg/kg, p.o.) reduces tumor area in an LNCaP prostate cancer mouse xenograft model when administered in combination with BAY 2315497.

Reference

- Berger, M., Wortmann, L., Buchgraber, P., *et al.* BAY-8400: A novel potent and selective DNA-PK inhibitor which shows synergistic efficacy in combination with targeted alpha therapies. *J. Med. Chem.* **64**(17), 12723-12737 (2021).

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, 09/04/2025

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