

PRODUCT INFORMATION

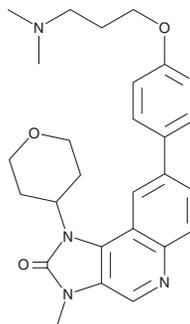


AZD 0156

Item No. 25648

CAS Registry No.: 1821428-35-6
Formal Name: 8-[6-[3-(dimethylamino)propoxy]-3-pyridinyl]-1,3-dihydro-3-methyl-1-(tetrahydro-2H-pyran-4-yl)-2H-imidazo[4,5-c]quinolin-2-one

MF: C₂₆H₃₁N₅O₃
FW: 461.6
Purity: ≥95%
UV/Vis.: λ_{max}: 210, 273 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

AZD 0156 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 0156 in the solvent of choice, which should be purged with an inert gas. AZD 0156 is soluble in the organic solvent DMSO.

Description

AZD 0156 is an inhibitor of the ataxia-telangiectasia mutated (ATM) kinase (IC₅₀s = 0.04 and 0.57 nM in enzyme and cell-based assays, respectively).¹ It is selective for ATM over other PI3K-related kinase (PIKK) family kinases including DNA-PK, mTOR, PI3Kα, PI3Kβ, PI3Kγ, and PI3Kδ (IC₅₀s = 0.14, 0.20, 0.32, 1.8, 1.1, and 0.27 μM, respectively, in enzyme assays). AZD 0156 (20 mg/kg) enhances the antitumor activity of irinotecan (Item No. 14180) in an SW620 colon cancer mouse xenograft model and olaparib (Item No. 10621) in an HBCx-10 breast cancer patient-derived xenograft (PDX) mouse model when administered at a dose of 5 mg/kg, leading to tumor regression in both models.

Reference

1. Pike, K.G., Barlaam, B., Cadogan, E., *et al.* The identification of potent, selective, and orally available inhibitors of ataxia telangiectasia mutated (ATM) kinase: The discovery of AZD0156 (8-{6-[3-(dimethylamino)propoxy]pyridin-3-yl}-3-methyl-1-(tetrahydro-2 H-pyran-4-yl)-1,3-dihydro-2 H-imidazo[4,5-c]quinolin-2-one). *J. Med. Chem.* **61**(9), 3823-3841 (2018).

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, 12/12/2022

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