

PRODUCT INFORMATION

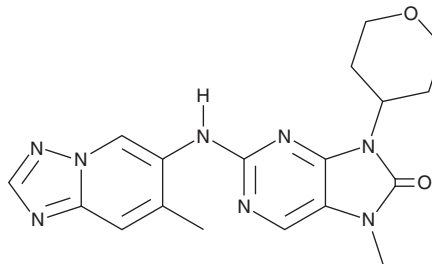


AZD 7648

Item No. 28598

CAS Registry No.: 2230820-11-6
Formal Name: 7,9-dihydro-7-methyl-2-[(7-methyl[1,2,4]triazolo[1,5-a]pyridin-6-yl)amino]-9-(tetrahydro-2H-pyran-4-yl)-8H-purin-8-one

MF: C₁₈H₂₀N₈O₂
FW: 380.4
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 263, 319 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.

Description

AZD 7648 is an inhibitor of DNA-dependent protein kinase (DNA-PK; IC₅₀ = 0.63 nM in an enzyme assay).¹ It is selective for DNA-PK over TTK (IC₅₀ = 5.012 μM), as well as JAK1, JAK2, JAK3, Aurora A, and Aurora B (IC₅₀s = >10 μM for all). AZD 7648 reduces tumor growth in a FaDu ATM KO pharynx cancer mouse xenograft model when administered at doses of 37.5 or 75 mg/kg twice per day and induces tumor regression at these doses when administered in combination with the PARP inhibitor olaparib (Item No. 10621).

Reference

1. Finlay, M. and Verschoyle, R. Amino-triazolopyridine compounds and their use in treating cancer. *Astrazeneca AB and Cancer Research Technology Limited. WO2018/114999AI* (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, 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