

PRODUCT INFORMATION

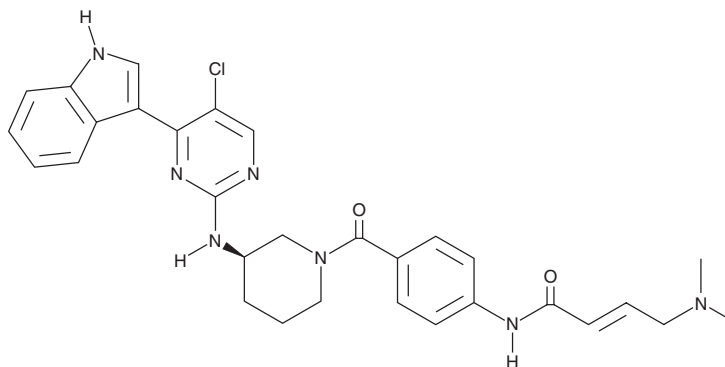


THZ531

Item No. 26386

CAS Registry No.: 1702809-17-3
Formal Name: (2E)-N-[4-[[[(3R)-3-[[5-chloro-4-(1H-indol-3-yl)-2-pyrimidinyl]amino]-1-piperidinyl]carbonyl]phenyl]-4-(dimethylamino)-2-butenamide

MF: C₃₀H₃₂ClN₇O₂
FW: 558.1
Purity: ≥95%
UV/Vis.: λ_{max}: 214, 251, 352 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

THZ531 is supplied as a crystalline solid. A stock solution may be made by dissolving the THZ531 in the solvent of choice, which should be purged with an inert gas. THZ531 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of THZ531 in these solvents is approximately 25, 20, and 30 mg/ml, respectively.

Description

THZ531 is an inhibitor of cyclin-dependent kinase 12 (Cdk12) and Cdk13 (IC₅₀s = 158 and 69 nM for Cdk12/cyclin K and Cdk13/cyclin K, respectively).¹ It is selective for Cdk12/cyclin K and Cdk13/cyclin K over Cdk9/cyclin T1 and Cdk7/cyclin H/MAT1 (IC₅₀s = 10.5 and 8.5 μM, respectively). THZ531 inhibits proliferation of Jurkat cells with an IC₅₀ value of 50 nM and induces apoptosis in a concentration-dependent manner. It also reduces expression of genes associated with the DNA damage response, including *BRCA1*, *FANCF*, and *ERCC4*, in Jurkat cells when used at a concentration of 50 nM.

Reference

1. Zhang, T., Kwiatkowski, N., Olson, C.M., *et al.* Covalent targeting of remote cysteine residues to develop CDK12 and CDK13 inhibitors. *Nat. Chem. Biol.* **12**(10), 876-884 (2016).

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, 09/13/2019

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