

# PRODUCT INFORMATION

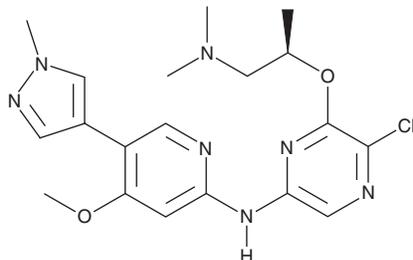


**CCT244747**

Item No. 25666

**CAS Registry No.:** 1404095-34-6  
**Formal Name:** 3-[(1R)-2-(dimethylamino)-1-methylethoxy]-5-[[4-methoxy-5-(1-methyl-1H-pyrazol-4-yl)-2-pyridinyl]amino]-2-pyrazinecarbonitrile

**MF:** C<sub>20</sub>H<sub>24</sub>N<sub>8</sub>O<sub>2</sub>  
**FW:** 408.5  
**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

CCT244747 is supplied as a solid. A stock solution may be made by dissolving the CCT244747 in the solvent of choice, which should be purged with an inert gas. CCT244747 is soluble in DMSO.

## Description

CCT244747 is a potent inhibitor of checkpoint kinase 1 (Chk1; IC<sub>50</sub> = 8 nM).<sup>1</sup> It is selective for Chk1 over FLT3, Chk2, and CDK1 (IC<sub>50</sub>s = 600, >10,000, and >10,000 nM, respectively) as well as a panel of 121 additional kinases at a concentration of 10 μM. CCT244747 reverses cell cycle arrest at the G<sub>2</sub>/M phase induced by etoposide (Item No. 12092) and SN-38 (Item No. 15632) in HT-29 cells and gemcitabine (Item Nos. 11690 | 9003096) in SW620 cells, indicating cellular G<sub>2</sub> checkpoint abrogation. It potentiates the genotoxicity of SN-38, gemcitabine, and etoposide in HT-29, SW620, MiaPaCa-2, and Calu-6 cells with potentiation index (PI) values ranging from 1.9-8.5, 2.6-12.2, 5-16, and 1.4-5.6, respectively. *In vivo*, CCT244747 (75 mg/kg), in combination with irinotecan (Item No. 14180) or gemcitabine, increases the delay of tumor growth in SW620 and Calu-6 mouse xenograft models, respectively.

## Reference

- Walton, M.I., Eve, P.D., Hayes, A., *et al.* CCT244747 is a novel potent and selective CHK1 inhibitor with oral efficacy alone and in combination with genotoxic anticancer drugs. *Clin. Cancer Res.* **18**(20), 5650-5661 (2012).

### 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, 10/19/2022

## 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