

PRODUCT INFORMATION



CCT245737

Item No. 28775

CAS Registry No.: 1489389-18-5
Formal Name: 5-[[4-[[[(2R)-2-morpholinylmethyl]amino]-5-(trifluoromethyl)-2-pyridinyl]amino]-2-pyrazinecarbonitrile

MF: C₁₆H₁₆F₃N₇O

FW: 379.3

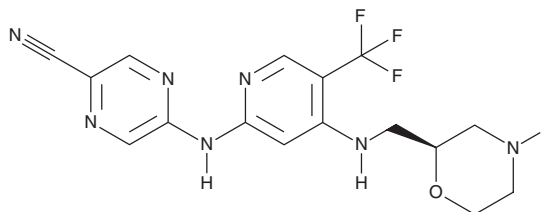
Purity: ≥98%

UV/Vis.: λ_{max}: 224, 293, 339 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

CCT245737 is supplied as a crystalline solid. A stock solution may be made by dissolving the CCT245737 in the solvent of choice, which should be purged with an inert gas. CCT245737 is soluble in the organic solvent DMSO.

Description

CCT245737 is a potent inhibitor of checkpoint kinase 1 (Chk1; IC₅₀ = 1.3 nM).¹ It is selective for Chk1 over Chk2 (IC₅₀ = 2,440 nM). CCT245737 inhibits G₂ checkpoint arrest induced by etoposide (Item No. 12092) in HT-29 colon cancer cells (IC₅₀ = 30 nM). *In vivo*, CCT245737 (300 mg/kg) inhibits Chk1 autophosphorylation induced by gemcitabine (Item No. 11690) in an SW620 colon cancer mouse xenograft model. It reduces tumor volume when administered alone or in combination with gemcitabine in an HT-29 mouse xenograft model. CCT245737 (150 mg/kg) reduces inguinal, brachial/axillary, and mesenteric lymph node weights in the Eμ-Myc driven transgenic mouse transplant model of lymph gland-infiltrating B cell lymphoma.²

References

- Osborne, J.D., Matthews, T.P., McHardy, T., *et al.* Multiparameter lead optimization to give an oral checkpoint kinase 1 (CHK1) inhibitor clinical candidate: (R)-5-((4-((morpholin-2-ylmethyl)amino)-5-(trifluoromethyl)pyridin-2-yl)amino)pyrazine-2-carbonitrile (CCT245737). *J. Med. Chem.* **59(11)**, 5221-5237 (2016).
- Walton, M.I., Eve, P.D., Hayes, A., *et al.* The clinical development candidate CCT245737 is an orally active CHK1 inhibitor with preclinical activity in RAS mutant NSCLC and Eμ-MYC driven B-cell lymphoma. *Oncotarget* **7(3)**, 2329-2342 (2015).

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, 12/05/2022

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