

PRODUCT INFORMATION



CCT241161

Item No. 25536

CAS Registry No.: 1163719-91-2
Formal Name: N-[4-[(3,4-dihydro-3-oxopyrido[2,3-b]pyrazin-8-yl)oxy]-2-(methylthio)phenyl]-N'-(3-(1,1-dimethylethyl)-1-phenyl-1H-pyrazol-5-yl)-urea

MF: C₂₈H₂₇N₇O₃S

FW: 541.6

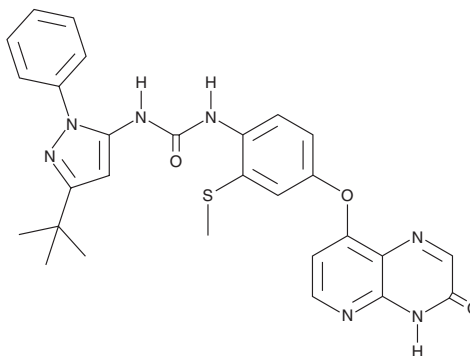
Purity: ≥98%

UV/Vis.: λ_{max}: 235 nm

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

CCT241161 is supplied as a solid. A stock solution may be made by dissolving the CCT241161 in the solvent of choice. CCT241161 is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

CCT241161 is a multi-kinase inhibitor that inhibits B-RAF, B-RAF^{V600E}, C-RAF, Src, and LCK (IC₅₀s = 252, 15, 6, 15, and 3 nM, respectively).¹ It is selective for RAFs, Src, LCK, and MAPKs in a panel of 63 kinases when used at a concentration of 1 μM. CCT241161 inhibits MEK and ERK signaling in B-RAF mutant WM266.4 cells, but not B-RAF wild-type D35 cells, and inhibits growth of B-RAF mutant melanoma cells in a concentration-dependent manner. *In vivo*, CCT241161 (20 mg/kg per day) induces tumor regression in a B-RAF mutant A375 mouse xenograft model. CCT241161 also inhibits ERK and Src signaling and induces tumor regression in B-RAF inhibitor-resistant patient-derived xenograft (PDX) mouse models including those resistant to both dabrafenib (Item No. 16989) and trametinib (Item No. 16292).

Reference

1. Girotti, M.R., Lopes, F.J.P., Preece, N., *et al.* Paradox-breaking RAF inhibitors that also target SRC are effective in drug-resistant BRAF mutant melanoma. *Cancer Cell* **31**(3), 85-96 (2017).

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, 10/20/2022

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