

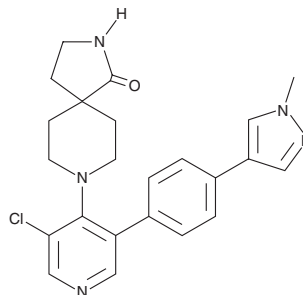
PRODUCT INFORMATION



CCT251545

Item No. 18761

CAS Registry No.: 1661839-45-7
Formal Name: 8-[3-chloro-5-[4-(1-methyl-1H-pyrazol-4-yl)phenyl]-4-pyridinyl]-2,8-diazaspiro[4.5]decan-1-one
MF: C₂₃H₂₄ClN₅O
FW: 421.9
Purity: ≥95%
UV/Vis.: λ_{max}: 274 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

CCT251545 is supplied as a crystalline solid. A stock solution may be made by dissolving the CCT251545 in the solvent of choice, which should be purged with an inert gas. CCT251545 is soluble in the organic solvent DMSO at a concentration of approximately 50 mg/ml.

Description

CCT251545 is an orally bioavailable inhibitor of Wnt signaling (IC₅₀ = 5 nM).¹ It is a potent and selective inhibitor of the human mediator complex-associated protein kinases Cdk8 and Cdk19.² Through its effects on Cdk8 and Cdk19, CCT251545 alters Wnt pathway-regulated gene expression.² It reduces tumor growth in COLO 205 human colon cancer xenografts after oral dosing (70 mg/kg bid) and in Wnt-dependent tumors.^{1,2}

References

1. Mallinger, A., Crumpler, S., Pichowicz, M., *et al.* Discovery of potent, orally bioavailable, small-molecule inhibitors of WNT signaling from a cell-based pathway screen. *J. Med. Chem.* **58**(4), 1717-1735 (2015).
2. Dale, T., Clarke, P.A., Esdar, C., *et al.* A selective chemical probe for exploring the role of CDK8 and CDK19 in human disease. *Nat. Chem. Biol.* **11**(12), 973-980 (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, 11/07/2022

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