

PRODUCT INFORMATION



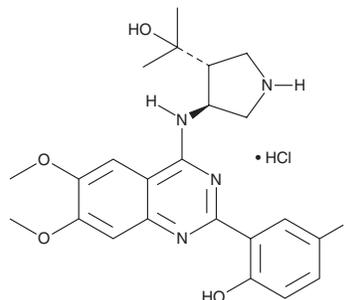
CCT241533 (hydrochloride)

Item No. 19178

Sold under license from Cancer Research Technology, LTD.

CAS Registry No.: 1431697-96-9
Formal Name: (3R,4S)-rel-4-[[2-(5-fluoro-2-hydroxyphenyl)-6,7-dimethoxy-4-quinazoliny]amino]- α,α -dimethyl-3-pyrrolidinemethanol, monohydrochloride

MF: C₂₃H₂₇FN₄O₄ • HCl
FW: 479.0
Purity: ≥98%
UV/Vis.: λ_{max} : 222, 255, 334 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

CCT241533 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the CCT241533 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. CCT241533 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CCT241533 (hydrochloride) in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of CCT241533 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of CCT241533 (hydrochloride) in PBS, pH 7.2, is approximately 500 $\mu\text{g/ml}$. We do not recommend storing the aqueous solution for more than one day.

Description

CCT241533 is a selective inhibitor of checkpoint kinase 2 (Chk2; IC₅₀ = 3 nM).¹ In human cancer cell lines, CCT241533 blocks Chk2 activity and potentiates the cytotoxicity of the PARP inhibitors, rucaparib (Item No. 15643) and olaparib (Item No. 10621).²

References

1. Caldwell, J.J., Welsh, E.J., Matijssen, C., *et al.* Structure-based design of potent and selective 2-(quinazolin-2-yl)phenol inhibitors of checkpoint kinase 2. *J. Med. Chem.* **54**(2), 580-590 (2011).
2. Anderson, V.E., Walton, M.I., Eve, P.D., *et al.* CCT241533 is a potent and selective inhibitor of CHK2 that potentiates the cytotoxicity of PARP inhibitors. *Cancer Res.* **71**(2), 463-472 (2011).

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, 11/29/2022

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