

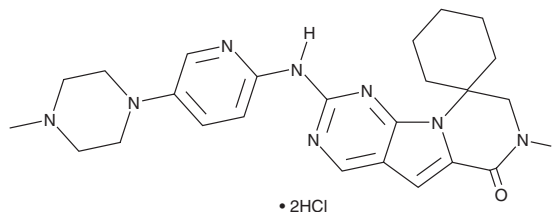
PRODUCT INFORMATION



Trilaciclib (hydrochloride)

Item No. 37297

CAS Registry No.: 1977495-97-8
Formal Name: 7',8'-dihydro-2'-[[5-(4-methyl-1-piperazinyl)-2-pyridinyl]amino]-spiro[cyclohexane-1,9'-(6'H)-pyrazino[1',2':1,5]pyrrolo[2,3-d]pyrimidin]-6'-one, dihydrochloride
Synonym: G1T28
MF: C₂₄H₃₀N₈O • 2HCl
FW: 519.5
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 277, 366 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

Trilaciclib (hydrochloride) is supplied as a solid. Aqueous solutions of trilaciclib (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. Trilaciclib (hydrochloride) is slightly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day. It is also slightly soluble in ethanol.

Description

Trilaciclib is an inhibitor of cyclin-dependent kinase 4 (Cdk4)/cyclin D1 and Cdk6/cyclin D3 (IC₅₀s = 1 and 4 nM, respectively).¹ It is selective for Cdk4/cyclin D1 and Cdk6/cyclin D3 over Cdk2/cyclin A, Cdk2/cyclin E, Cdk5/p25, Cdk5/p35, Cdk7/cyclin H/Mat1, and Cdk9/cyclin T (IC₅₀s = 1,290, 2,510, 1,710, 1,240, 4,640, and 50 nM, respectively). It halts the cell cycle at the G₁ phase in Cdk4/6-dependent Hs68 cells (EC₅₀ = 30 nM) but not Cdk4/6-independent A2058 cells. Trilaciclib (10-1,000 nM) prevents DNA damage and apoptosis induced by the DNA enzyme topoisomerase I inhibitor camptothecin (Item No. 11694) in Hs68 cells used as a model of Cdk4/6-dependent hematopoietic stem and progenitor cells (HSPCs). It induces reversible cell cycle arrest of HSPCs in mice when administered at doses of 50, 100, or 150 mg/kg and prevents apoptosis of bone marrow cells induced by etoposide (Item No. 12092) in mice at 100 mg/kg. Trilaciclib (150 mg/kg) also prevents myelosuppression induced by 5-fluorouracil (5-FU; Item No. 14416) in mice. Formulations containing trilaciclib have been used in the treatment of late-stage small cell lung cancer to decrease the incidence of chemotherapy-induced myelosuppression.

Reference

1. Bisi, J.E., Sorrentino, J.A., Roberts, P.J., et al. Preclinical characterization of G1T28: A novel CDK4/6 inhibitor for reduction of chemotherapy-induced myelosuppression. *Mol. Cancer Ther.* **15**(5), 783-793 (2016).

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, 04/18/2023

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