

PRODUCT INFORMATION

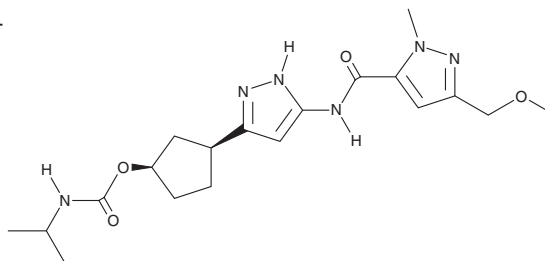


PF-07104091

Item No. 41356

CAS Registry No.: 2460249-19-6
Formal Name: carbamic acid N-(1-methylethyl)-(1R,3S)-3-[5-[[[3-(methoxymethyl)-1-methyl-1H-pyrazol-5-yl]carbonyl]amino]-1H-pyrazol-3-yl]cyclopentyl ester

Synonym: Tagtociclib
MF: C₁₉H₂₈N₆O₄
FW: 404.5
Purity: ≥95%
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

PF-07104091 is supplied as a solid. A stock solution may be made by dissolving the PF-07104091 in the solvent of choice, which should be purged with an inert gas. PF-07104091 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

PF-07104091 is an inhibitor of cyclin-dependent kinase 2 (Cdk2; $K_i = 1.16$ nM).¹ It is selective for Cdk2 over glycogen synthase kinase 3 β (GSK3 β ; $K_i = 537.81$ nM). PF-07104091 decreases proliferation of TOV-21G and OVCAR-3 ovarian and HCT116 colorectal cancer cells (IC_{50} s = 4.8, 0.59, and 0.88 μ M, respectively).² *In vivo*, PF-07104091 (50 mg/kg per day) reduces tumor volume and weight in an HCT116 mouse xenograft model.

References

- Behenna, D.C., Freeman-Cook, K.D., Hoffman, R.L., *et al.* CDK2 Inhibitors. *Pfizer, Inc.* **WO 2020/157652 A2** (2020).
- Liu, Z., Yang, Y., Sun, X., *et al.* Discovery of novel antitumor small-molecule agent with dual action of CDK2/p-RB and MDM2/p53. *Molecules* **29**(3), 725 (2024).

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, 08/16/2024

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