

PRODUCT INFORMATION



PF-06726304

Item No. 44260

CAS Registry No.: 1616287-82-1
Formal Name: 5,8-dichloro-2-[(1,2-dihydro-4,6-dimethyl-2-oxo-3-pyridinyl)methyl]-7-(3,5-dimethyl-4-isoxazolyl)-3,4-dihydro-1(2H)-isoquinolinone

MF: C₂₂H₂₁Cl₂N₃O₃

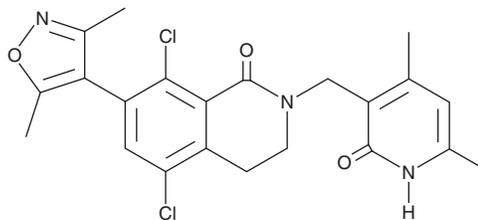
FW: 446.3

Purity: ≥98%

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-06726304 is supplied as a solid. A stock solution may be made by dissolving the PF-06726304 in the solvent of choice, which should be purged with an inert gas. PF-06726304 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

PF-06726304 is an inhibitor of enhancer of zeste homolog 2 (EZH2; $K_i = 0.7$ nM).¹ *In vivo*, PF-06726304 (200 and 300 mg/kg) reduces tumor volume and intratumor histone methylation in a Karpas-422 mouse xenograft model.

Reference

1. Kung, P.P., Rui, E., Bergqvist, S., *et al.* Design and synthesis of pyridone-containing 3,4-dihydroisoquinoline-1(2H)-ones as a novel class of enhancer of zeste homolog 2 (EZH2) inhibitors. *J. Med. Chem.* **59**(18), 8306-8325 (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, 09/15/2025

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