

PRODUCT INFORMATION

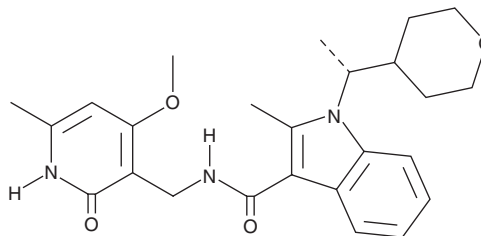


CPI-360

Item No. 19125

CAS Registry No.: 1802175-06-9
Formal Name: N-[(1,2-dihydro-4-methoxy-6-methyl-2-oxo-3-pyridinyl)methyl]-2-methyl-1-[(1R)-1-(tetrahydro-2H-pyran-4-yl)ethyl]-1H-indole-3-carboxamide

MF: $C_{25}H_{31}N_3O_4$
FW: 437.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 212, 292 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

CPI-360 is supplied as a crystalline solid. A stock solution may be made by dissolving the CPI-360 in the solvent of choice, which should be purged with an inert gas. CPI-360 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CPI-360 in these solvents is approximately 30, 25, and 10 mg/ml, respectively.

CPI-360 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CPI-360 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. CPI-360 has a solubility of approximately 0.12 mg/ml in a 1:7 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CPI-360 is a selective EZH2 inhibitor with IC_{50} values of 0.5 and 2.5 nM for wild-type EZH2 and Y641N mutant EZH2, respectively.¹ It decreases cellular levels of H3K27me3 and H3K27me2 (EC_{50} s = 56 and 65 nM, respectively), triggering cell cycle arrest and ultimately resulting in apoptosis in a large panel of non-Hodgkin's lymphoma (NHL) cell lines.¹ Twice daily, subcutaneous administration of 200 mg/kg of CPI-360 reduced tumor growth of KARPAS-422 xenografts in mice.¹

Reference

1. Bradley, W.D., Arora, S., Busby, J., *et al.* EZH2 inhibitor efficacy in non-Hodgkin's lymphoma does not require suppression of H3K27 monomethylation. *Chem. Biol.* **21**, 1-13 (2014).

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, 10/03/2022

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