

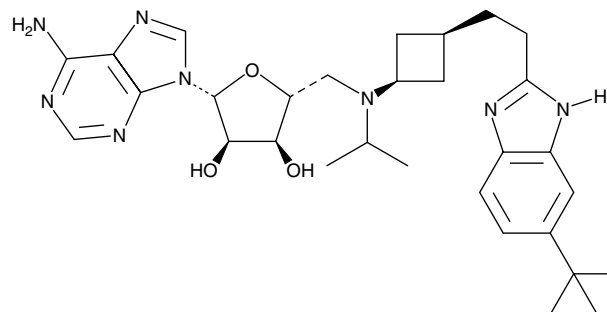
Product Information



EPZ5676

Item No. 16175

CAS Registry No.: 1380288-87-8
Formal Name: 9-[5-deoxy-5-[[*cis*-3-[2-[6-(1,1-dimethylethyl)-1H-benzimidazol-2-yl]ethyl]cyclobutyl](1-methylethyl)amino]-β-D-ribofuranosyl]-9H-purin-6-amine
MF: C₃₀H₄₂N₈O₃
FW: 562.7
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 255 nm



Laboratory Procedures

For long term storage, we suggest that EPZ5676 be stored as supplied at -20°C. It should be stable for at least two years. EPZ5676 is supplied as a crystalline solid. A stock solution may be made by dissolving the EPZ5676 in the solvent of choice. EPZ5676 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of EPZ5676 in ethanol and DMF is approximately 30 mg/ml and approximately 25 mg/ml in DMSO.

EPZ5676 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, EPZ5676 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. EPZ5676 has a solubility of approximately 0.1 mg/ml in a 1:8 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

DOT1L is a non-SET domain containing methyltransferase whose function is important for the transcriptional activation of certain genes, DNA damage repair, and cell cycle regulation. DOT1L is known to play an essential role in MLL-rearranged leukemias, and thus has emerged as a drug target for these leukemias.^{1,2} EPZ5676 is a highly potent aminonucleoside inhibitor of DOT1L histone methyltransferase activity (K_i = 80 pM) that demonstrates 37,000-fold selectivity over other methyltransferases.^{3,4} In various human leukemia cell lines, EPZ5676 has been shown to inhibit the methylation of lysine 79 on histone 3 (IC₅₀s = 3-5 nM), decreasing MLL-fusion target gene expression (IC₅₀s = 53-67 nM).³ It can selectively inhibit the proliferation of acute leukemia cell lines bearing MLL translocations (IC₅₀s range from 3.5 nM - 1.3 μM).³ Continuous intravenous infusion of EPZ5676 at a dose of 35 mg/kg has been reported to cause complete and sustained tumor regression in a rat xenograft model of MLL-rearranged leukemia.³

Reference

1. Nguyen, A.T., He, J., Taranova, O., *et al.* Essential role of DOT1L in maintaining normal adult hematopoiesis. *Cell Res.* **21(9)**, 1370-1373 (2011).
2. Bernt, K.M. and Armstrong, S.A. A role for DOT1L in MLL-rearranged leukemias. *Epigenomics* **3(6)**, 667-670 (2011).
3. Daigle, S.R., Olhava, E.J., Therkelsen, C.A., *et al.* Potent inhibition of DOT1L as treatment of MLL-fusion leukemia. *Blood* **122(6)**, 1017-1025 (2013).
4. Basavapathruni, A., Olhava, E.J., Daigle, S.R., *et al.* Nonclinical pharmacokinetics and metabolism of EPZ5676, a novel DOT1L histone methyltransferase inhibitor. *Biopharm. Drug Dispos.* **35(4)**, 237-252 (2014).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/16175

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 08/01/2014

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com