

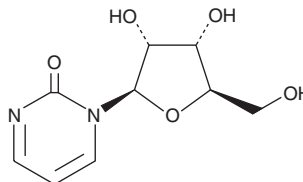
PRODUCT INFORMATION



Zebularine

Item No. 10975

CAS Registry No.: 3690-10-6
Formal Name: 1H-β-D-ribofuranosyl-2-pyrimidinone
MF: C₉H₁₂N₂O₅
FW: 228.2
Purity: 98%
UV/Vis.: λ_{max}: 310 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

Zebularine is supplied as a crystalline solid. A stock solution may be made by dissolving the zebularine in the solvent of choice, which should be purged with an inert gas. Zebularine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of zebularine in these solvents is approximately 0.25, 14, and 5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of zebularine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of zebularine in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Zebularine, a cytidine analog, is a DNA methylation inhibitor that acts by forming a covalent complex with DNA methyltransferases (DNMTs).¹ By stabilizing the binding of DNMTs to DNA, zebularine hinders methylation and decreases dissociation, trapping DNMT and preventing its turnover.² Zebularine preferentially targets cancer cells in various human carcinoma cell lines demonstrating IC₂₀ values of 20, 10, 20, and 5 μM for reducing proliferation of TK6, Jurkat, KG-1, and HCT116 cells, respectively.³ Zebularine exhibits low toxicity in mice even after prolonged administration (0.2 mg/ml drinking water for 113 days in a mouse intestinal cancer model).⁴

References

1. Zhou, L., Cheng, X., Connolly, B.A., *et al.* Zebularine: A novel DNA methylation inhibitor that forms a covalent complex with DNA methyltransferases. *J. Mol. Biol.* **321**(4), 591-599 (2002).
2. Champion, C., Guianvarc'h, D., Sénamaud-Beaufort, C., *et al.* Mechanistic insights on the inhibition of c5 DNA methyltransferases by zebularine. *PLoS One* **5**(8), e12388 (2010).
3. Stresemann, C., Brueckner, B., Musch, T., *et al.* Functional diversity of DNA methyltransferase inhibitors in human cancer cell lines. *Cancer Res.* **66**(5), 2794-2800 (2006).
4. Yoo, C.B., Chuang, J.C., Byun, H.M., *et al.* Long-term epigenetic therapy with oral zebularine has minimal side effects and prevents intestinal tumors in mice. *Cancer Prev. Res.* **1**(4), 233-240 (2008).

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, 02/21/2024

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