

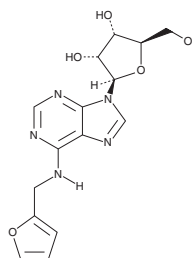
PRODUCT INFORMATION



Kinetin Riboside

Item No. 36570

CAS Registry No.: 4338-47-0
Formal Name: N-(2-furanylmethyl)-adenosine
Synonyms: N⁶-Furfuryladosine, NSC 120958
MF: C₁₅H₁₇N₅O₅
FW: 347.3
Purity: ≥98%
UV/Vis.: λ_{max}: 267 nm
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

Kinetin riboside is supplied as a solid. A stock solution may be made by dissolving the kinetin riboside in the solvent of choice, which should be purged with an inert gas. Kinetin riboside is soluble in the organic solvent dimethyl formamide at a concentration of approximately 3 mg/ml.

Description

Kinetin riboside is a purine derivative and nucleoside with anticancer activity.¹ It induces the production of reactive oxygen species (ROS), cellular ATP depletion, and apoptosis in HepG2 cells. Kinetin riboside (10 μM) reduces cyclin D1 and cyclin D2 levels, as well as induces cell cycle arrest at the G₁ phase and apoptosis in patient-derived multiple myeloma cells.² *In vivo*, kinetin riboside (85 mg/kg) reduces tumor growth in OCI-My5 and RPMI-8226 myeloma mouse xenograft models.

References

- Orlicka-Płocka, M., Gurda-Wozna, D., Fedoruk-Wyzomirska, A., *et al.* Circumventing the Crabtree effect: Forcing oxidative phosphorylation (OXPHOS) via galactose medium increases sensitivity of HepG2 cells to the purine derivative kinetin riboside. *Apoptosis* **25**(11-12), 835-852 (2020).
- Tiedemann, R.E., Mao, X., Shi, C.-X., *et al.* Identification of kinetin riboside as a repressor of CCND1 and CCND2 with preclinical antimyeloma activity. *J. Clin. Invest.* **118**(5), 1750-1764 (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, 12/20/2022

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