

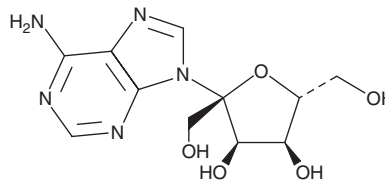
PRODUCT INFORMATION



Psicofuranine

Item No. 19574

CAS Registry No.: 1874-54-0
Formal Name: 1'-C-(hydroxymethyl)-adenosine
Synonyms: Angustmycin C, NSC 53104
MF: C₁₁H₁₅N₅O₅
FW: 297.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Special Conditions: Avoid freeze/thaw cycles; After reconstitution protect from light

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Psicofuranine is supplied as a solid. A stock solution may be made by dissolving the psicofuranine in the solvent of choice. Psicofuranine is soluble in the organic solvent DMSO.

Description

Psicofuranine is an antibiotic and antitumor compound that acts as an inhibitor of xanthosine monophosphate (XMP) aminase (IC₅₀ = 67 μM), causing guanine deficiency in enteric bacteria.¹⁻³

References

1. Hanka, L.J. Mechanism of action of psicofuranine. *J. Bacteriol.* **80**, 30-36 (1960).
2. Udaka, S. and Moyed, H.S. Inhibition of parental and mutant xanthosine 5'-phosphate aminases by psicofuranine. *J. Biol. Chem.* **238**, 2797-2803 (1963).
3. Fukuyama, T.T. Formation of an adenylyl xanthosine monophosphate intermediate by xanthosine 5'-phosphate aminase and its inhibition by psicofuranine. *J. Biol. Chem.* **241**(20), 4745-4749 (1966).

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, 12/22/2022

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