

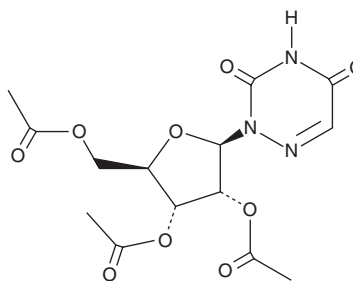
PRODUCT INFORMATION



Azaribine

Item No. 40131

CAS Registry No.: 2169-64-4
Formal Name: 2-(2,3,5-tri-O-acetyl-β-D-ribofuranosyl)-1,2,4-triazine-3,5(2H,4H)-dione
Synonyms: 6-Azauridine Triacetate, NSC 67239
MF: C₁₄H₁₇N₃O₉
FW: 371.3
Purity: ≥95%
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

Azaribine is supplied as a solid. A stock solution may be made by dissolving the azaribine in the solvent of choice, which should be purged with an inert gas. Azaribine is soluble in chloroform and acetonitrile.

Description

Azaribine is a prodrug form of the nucleoside analog and immunosuppressant 6-azauridine.^{1,2} It inhibits viral replication in influenza A- or influenza B virus-infected MDCK cells (EC₅₀s = 0.29-0.8 μM).³ Formulations containing azaribine have previously been used in the treatment of psoriasis.

References

1. Merkel, P.A., Letourneau, E.N., and Polisson, R.P. Investigational agents for rheumatoid arthritis. *Rheum. Dis. Clin. North Am.* **21(3)**, 779-796 (1995).
2. Creasey, W.A., Fink, M.E., Handschumacher, R.E., *et al.* Clinical and pharmacological studies with 2',3',5'-triacetyl-6-azauridine. *Cancer Res.* **23**, 444-453 (1963).
3. Park, J.-G., Ávila-Pérez, G., Nogales, A., *et al.* Identification and characterization of novel compounds with broad-spectrum antiviral activity against influenza A and B viruses. *J. Virol.* **94(7)**, e02149-19 (2020).

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, 12/06/2023

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