

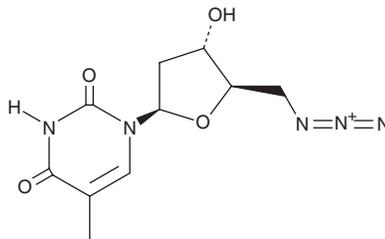
# PRODUCT INFORMATION



## 5'-Azido-5'-deoxythymidine

Item No. 44620

CAS Registry No.: 19316-85-9  
Synonyms: 5'-AZT, NSC 254064  
MF:  $C_{10}H_{13}N_5O_4$   
FW: 267.2  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

5'-Azido-5'-deoxythymidine is supplied as a solid. A stock solution may be made by dissolving the 5'-azido-5'-deoxythymidine in the solvent of choice, which should be purged with an inert gas. 5'-Azido-5'-deoxythymidine is soluble ( $\geq 10$  mg/ml) in DMSO.

### Description

5'-Azido-5'-deoxythymidine is a nucleoside derivative and an inhibitor of poly(ADP-ribose) polymerase (PARP;  $K_i = 12 \mu M$ ).<sup>1</sup> It is also a synthetic intermediate that has been used in the synthesis of antiviral prodrugs.<sup>2</sup>

### References

1. Pivazyan, A.D., Birks, E.M., Wood, T.G., *et al.* Inhibition of poly(ADP-ribose)polymerase activity by nucleoside analogs of thymidine. *Biochem. Pharmacol.* **44(5)**, 947-953 (1992).
2. Lin, T.-S. Synthesis and *in vitro* antiviral activity of 3'-O-acyl derivatives of 5'-amino-5'-deoxythymidine: Potential prodrugs for topical application. *J. Pharm. Sci.* **73(11)**, 1568-1571 (1984).

#### 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, 12/17/2025

#### 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