

PRODUCT INFORMATION



5'-Deoxy-5'-iodothymidine

Item No. 44622

CAS Registry No.: 25953-14-4

Synonym: NSC 338954

MF: $C_{10}H_{13}IN_2O_4$

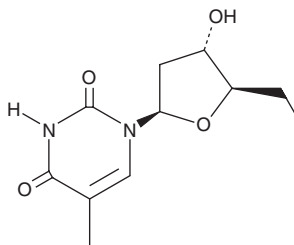
FW: 352.1

Purity: $\geq 95\%$

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

5'-Deoxy-5'-iodothymidine is supplied as a solid. A stock solution may be made by dissolving the 5'-deoxy-5'-iodothymidine in the solvent of choice, which should be purged with an inert gas. 5'-Deoxy-5'-iodothymidine is soluble (≥ 10 mg/ml) in DMSO.

Description

5'-Deoxy-5'-iodothymidine is an iodinated derivative of thymidine (Item No. 20519).¹ It has been used as an intermediate in the synthesis of thymidylate kinase (TMPK) inhibitors and as a starting material for anticancer agents.^{1,2}

References

1. Suthagar, K., Jiao, W., Munier-Lehmann, H., *et al.* Synthesis of sulfamide analogues of deoxythymidine monophosphate as potential inhibitors of mycobacterial cell wall biosynthesis. *Carbohydr. Res.* **457**, 32-40 (2018).
2. Chen, X., Bastow, K., Goz, B., *et al.* Synthesis and evaluation of novel thymidine analogs as antitumor and antiviral agents. *J. Med. Chem.* **39**(17), 3412-3417 (1996).

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, 12/02/2025

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