

PRODUCT INFORMATION



2',3'-Dideoxythymidine-5'-O-triphosphate (sodium salt)

Item No. 38376

Formal Name: ((2S,5R)-5-(5-methyl-2,4-dioxo-3,4-dihydropyrimidin-1(2H)-yl)tetrahydrofuran-2-yl)methyl triphosphate

Synonym: ddTTP

MF: C₁₀H₁₃N₂O₁₃P₃ • 4Na

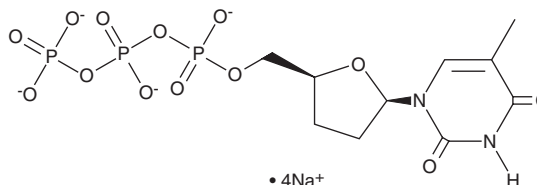
FW: 554.1

Purity: ≥95%

Supplied as: A solution in water

Storage: -80°C

Stability: ≥2 years



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

Description

2',3'-Dideoxythymidine-5'-O-triphosphate (ddTTP) is a reverse transcriptase inhibitor ($K_{i,s} = 40$ and 25 nM for the visna virus and HIV-1 enzymes, respectively).^{1,2} It also inhibits DNA polymerase γ ($K_i = 400$ nM for the bovine enzyme).³ ddTTP has been used to terminate chain extension produced by the *Taq* polymerase in PCR assays.⁴

References

1. Frank, K.B., McKernan, P.A., Smith, R.A., *et al.* Visna virus as an in vitro model for human immunodeficiency virus and inhibition by ribavirin, phosphonoformate, and 2',3'-dideoxynucleosides. *Antimicrob. Agents Chemother.* **31(9)**, 1369-1374 (1987).
2. Ueno, T., Shirasaka, T., and Mitsuya, H. Enzymatic characterization of human immunodeficiency virus type 1 reverse transcriptase resistant to multiple 2',3'-dideoxynucleoside 5'-triphosphates. *J. Biol. Chem.* **270(40)**, 23605-23611 (1995).
3. Li, Y., Mitaxov, V., and Waksman, G. Structure-based design of *Taq* DNA polymerases with improved properties of dideoxynucleotide incorporation. *Proc. Natl. Acad. Sci. USA* **96(17)**, 9491-9496 (1999).
4. Izuta, S., Saneyoshi, T., Sakurai, T., *et al.* The 5'-triphosphates of 3'-azido-3'-deoxythymidine and 2', 3'-dideoxynucleosides inhibit DNA polymerase γ by different mechanisms. *Biochim. Biophys. Res. Commun.* **179(2)**, 776-783 (1991).

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, 04/06/2023

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