

PRODUCT INFORMATION

2',3'-Dideoxycytidine-5'-O-(1-thiotriphosphate) (sodium salt)

Item No. 38461

Formal Name: 2',3'-dideoxy-cytidine, 5'→P''-ester with thiotriphosphoric acid ((HO)2P(O)OP(O)(OH)OP(O)(OH)(SH)), tetrasodium salt

Synonyms: α[S]ddCTP, ddCTP-α-S

MF: C₉H₁₂N₃O₁₁P₃S • 4Na

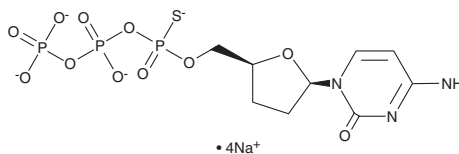
FW: 555.2

Purity: ≥95% (a mixture of Rp and Sp isomers)

Supplied as: A solution in water

Storage: -80°C

Stability: ≥4 years



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

Description

2',3'-Dideoxycytidine-5'-O-(1-thiotriphosphate) (ddCTP-α-S) is a sulfur-containing nucleoside triphosphate derivative. Thiol-containing dideoxynucleoside triphosphates (α-S-ddNTPs) incorporated into allele-specific primers prevent complete digestion by 5'-phosphodiesterase, which allows for alkylation and identification of SNP-containing alleles.¹ ddCTP-α-S has been used for chain extension in PCR assays.² ddCTP-α-S is a mixture of Rp and Sp isomers.

References

1. Sauer, S., Lechner, D., Berlin, K.D., et al. A novel procedure for efficient genotyping of single nucleotide polymorphisms. *Nucleic Acids Res.* **28(5)**, E13 (2000).
2. Sauer, S. and Gut, I.G. Extension of the GOOD assay for genotyping single nucleotide polymorphisms by matrix-assisted laser desorption/ionization mass spectrometry. *Rapid Commun. Mass Spectrom.* **17(12)**, 1265-1272 (2003).

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, 04/15/2026

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