

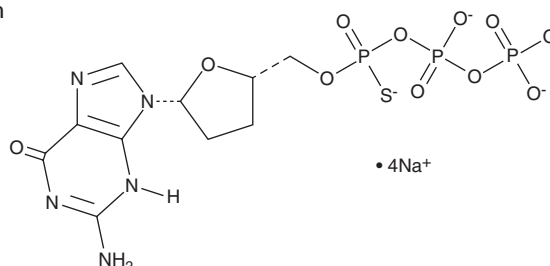
PRODUCT INFORMATION



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

Item No. 38462

Formal Name: 2',3'-dideoxy-guanosine, 5'→P''-ester with thiotriphosphoric acid ((HO)₂P(O)OP(O)(OH)OP(O)(OH)(SH)), tetrasodium salt
Synonyms: ddGTP-α-S, α-S-ddGTP
MF: C₁₀H₁₂N₅O₁₁P₃S • 4Na
FW: 595.2
Purity: ≥95% (a mixture of Rp and Sp isomers)
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'-Dideoxyguanosine-5'-O-(1-thiotriphosphate) (ddGTP-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 ddGTP-α-S has been used for chain extension using a mutated polymerase selective for dideoxynucleotides and chain termination in allele-specific PCR assays.^{1,2} ddGTP-α-S is a mixture of Rp and Sp isomers.

References

1. Tost, J., Schatz, P.F., Schuster, M., *et al.* Analysis and accurate quantification of CpG methylation by MALDI mass spectrometry. *Nucleic Acids Res.* **31(9)**, e50 (2003).
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, 06/08/2023

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