

# PRODUCT INFORMATION



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

Item No. 38375

**Formal Name:** 2',3'-dideoxy-guanosine  
5'-(tetrahydrogen triphosphate),  
tetrasodium salt

**Synonym:** ddGTP

**MF:** C<sub>10</sub>H<sub>12</sub>N<sub>5</sub>O<sub>12</sub>P<sub>3</sub> • 4Na

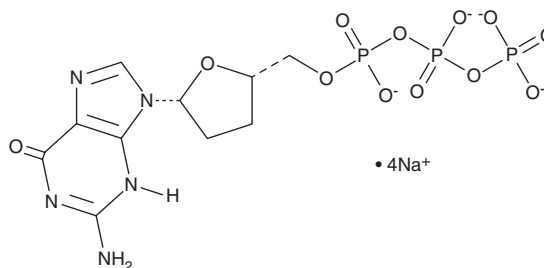
**FW:** 579.1

**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'-Dideoxyguanosine-5'-O-triphosphate (ddGTP) is a reverse transcriptase inhibitor ( $K_i$ s = 13 and 6.7 nM for the visna virus and HIV-1 enzymes, respectively).<sup>1,2</sup> ddGTP has been used in the synthesis of oligonucleotides containing a terminal ddG that are resistant to cleavage by severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) exonuclease.<sup>3</sup> ddGTP has also been used to terminate chain extension produced by the *Taq* polymerase in PCR assays.<sup>4</sup>

### 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. Wang, X., Tao, C., Morozova, I., *et al.* Identifying structural features of nucleotide analogues to overcome SARS-CoV-2 exonuclease activity. *Viruses* **14(7)**, 1413 (2022).
4. 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).

#### 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, 04/03/2023

#### 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