

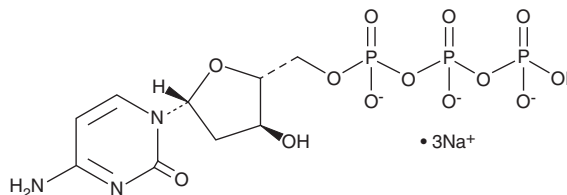
PRODUCT INFORMATION



2'-Deoxycytidine-5'-O-triphosphate (sodium salt)

Item No. 33484

CAS Registry No.: 109909-44-6
Formal Name: 2'-deoxy-cytidine 5'-(tetrahydrogen triphosphate), trisodium salt
Synonym: dCTP
MF: $C_9H_{13}N_3O_{13}P_3 \cdot 3Na$
FW: 533.1
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 273 nm
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

2'-Deoxycytidine-5'-O-triphosphate (dCTP) (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the dCTP (sodium salt) in water. The solubility of dCTP (sodium salt) in water is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

dCTP is a nucleoside triphosphate.¹ It has commonly been used in PCR reaction mixtures as a substrate for DNA synthesis by DNA polymerases.

Reference

1. Mohsen, M.G., Ji, D., and Kool, E.T. Polymerase synthesis of four-base DNA from two stable dimeric nucleotides. *Nucleic Acids Res.* **47(18)**, 9495-9501 (2019).

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, 08/14/2023

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