

PRODUCT INFORMATION

5'-DMTr-dC (Ac)-Methyl Phosphonamidite

Item No. 43384

CAS Registry No.: 178094-78-5

Formal Name: N-acetyl-5'-O-[bis(4-methoxyphenyl)phenylmethyl]-2'-deoxy-cytidine, 3'-[P-methyl-N-bis(1-methylethyl)phosphonamidite]

Synonyms: 5'-DMT-N⁴-Acetyldeoxycytidine
3'-Methylphosphonamidite, 5'-DMT dC^{Ac} 3'-Methylphosphonamidite

MF: C₃₉H₄₉N₄O₇P

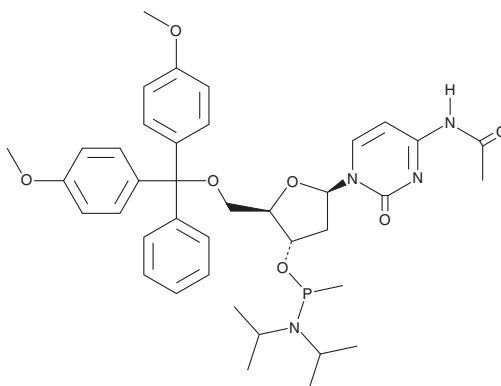
FW: 716.8

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

5'-DMTr-dC (Ac)-methyl phosphonamidite is supplied as a solid. A stock solution may be made by dissolving the 5'-DMTr-dC (Ac)-methyl phosphonamidite in the solvent of choice, which should be purged with an inert gas. 5'-DMTr-dC (Ac)-methyl phosphonamidite is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

5'-DMTr-dC (Ac)-methyl phosphonamidite is a nucleoside building block that has been used in the synthesis of oligonucleoside methylphosphonates.¹

Reference

1. Reddy, M.P., Farooqui, F., and Hanna, N.B. Elimination of transamination side product by the use of dC^{Ac} methylphosphonamidite in the synthesis of oligonucleoside methylphosphonates. *Tetrahedron Lett.* **37**(48), 8691-8694 (1996).

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, 02/02/2026

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