

PRODUCT INFORMATION



5-Me-3'-dU-2'-phosphoramidite

Item No. 43328

CAS Registry No.: 142103-12-6

Formal Name: 5'-O-[bis(4-methoxyphenyl)phenylmethyl]-3'-deoxy-5-methyluridine, 2'-[2-cyanoethyl N,N-bis(1-methylethyl)phosphoramidite]
Synonym: 3'-Deoxy-5'-O-DMT-5-methyluridine
2'-CE Phosphoramidite

MF: $C_{40}H_{49}N_4O_8P$

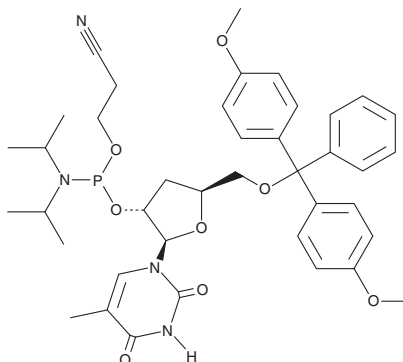
FW: 744.8

Purity: $\geq 98\%$

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-Me-3'-dU-2'-phosphoramidite is supplied as a solid. A stock solution may be made by dissolving the 5-Me-3'-dU-2'-phosphoramidite in the solvent of choice, which should be purged with an inert gas. 5-Me-3'-dU-2'-phosphoramidite is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

5-Me-3'-dU-2'-phosphoramidite is a nucleoside building block.^{1,2} It has been used in the synthesis of 2'5'-linked oligonucleotides.

References

1. Osawa, T., Hitomi, Y., Wakita, S., *et al.* Synthesis and hybridizing properties of isoDNAs including 3'-O,4'-C-ethyleneoxy-bridged 5-methyluridine derivatives. *Bioorg. Med. Chem.* **26(14)**, 3875-3881 (2018).
2. Rizzo, C.J., Dougherty, J.P., and Breslow, R. 3'-Deoxy-2'-phosphoramidites of adenosine and 5-methyluridine used for the solid phase synthesis of unnatural 3'-deoxy-2'-5''-oligonucleotides. *Tetrahedron Lett.* **33(29)**, 4129-4132 (1992).

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, 04/28/2025

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