

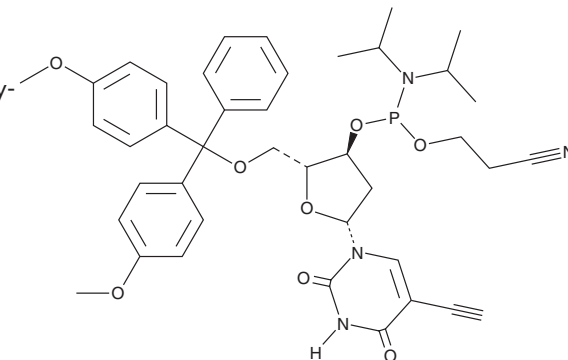
PRODUCT INFORMATION



5-Ethynyl DU CEP

Item No. 43315

CAS Registry No.: 615288-66-9
Formal Name: 5'-O-[bis(4-methoxyphenyl)phenylmethyl]-2'-deoxy-5-ethynyl-3'-[2-cyanoethyl N,N-bis(1-methylethyl)phosphoramidite] uridine
Synonyms: EdU CEP, 5-Ethynyl-2'-deoxyuridine 2-Cyanoethyl-N,N-diisopropylphosphoramidite
MF: C₄₁H₄₇N₄O₈P
FW: 754.8
Purity: ≥95%
Supplied as: A crystalline 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-Ethynyl DU CEP is supplied as a solid. A stock solution may be made by dissolving the 5-ethynyl DU CEP in the solvent of choice, which should be purged with an inert gas. 5-Ethynyl DU CEP is soluble (≥10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

5-Ethynyl DU CEP is a clickable nucleoside building block.¹ It has been used in the synthesis of spin-labeled oligonucleotides.

Reference

1. Jakobsen, U., Shelke, S.A., Vogel, S., *et al.* Site-directed spin-labeling of nucleic acids by click chemistry: Detection of abasic sites in duplex DNA by EPR spectroscopy. *J. Am. Chem. Soc.* **132**(30), 10424-10428 (2010).

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, 07/03/2025

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