

PRODUCT INFORMATION



2'-O-Propargyl A(Bz)-3'-phosphoramidite

Item No. 43628

CAS Registry No.: 171486-59-2
Formal Name: 3'-[2-cyanoethyl N,N-bis(1-methylethyl)phosphoramidite], N-benzoyl-5'-O-[bis(4-methoxyphenyl)phenylmethyl]-2'-O-2-propyn-1-yl-adenosine

MF: C₅₀H₅₄N₇O₈P

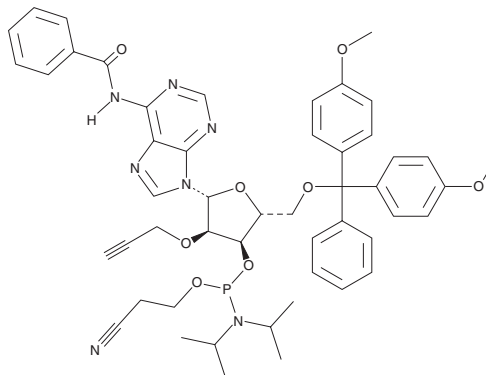
FW: 912.0

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

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

Description

2'-O-Propargyl A(Bz)-3'-phosphoramidite is a nucleoside building block that has been used in the synthesis of propargylated ribonucleosides for DNA crosslinking.¹

Reference

1. Pujari, S.S. and Seela, F. Cross-linked DNA: Propargylated ribonucleosides as "click" ligation sites for bifunctional azides. *J. Org. Chem.* **77**(9), (2012).

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, 10/09/2025

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