

PRODUCT INFORMATION

N⁶-Bz-5'-O-DMTr-2'-O-hexadecanyl Adenosine 3'-CED Phosphoramidite

Item No. 43602

CAS Registry No.: 2382942-35-8

Formal Name: 3'-[2-cyanoethyl N,N-bis(1-methylethyl)
phosphoramidite] N-benzoyl-5'-O-
[bis(4-methoxyphenyl)phenylmethyl]-2'
O-hexadecyl-adenosine

Synonyms: N⁶-Bz-5'-O-DMTr-2'-O-palmityl Adenosine
3'-CED Phosphoramidite

MF: C₆₃H₈₄N₇O₈P

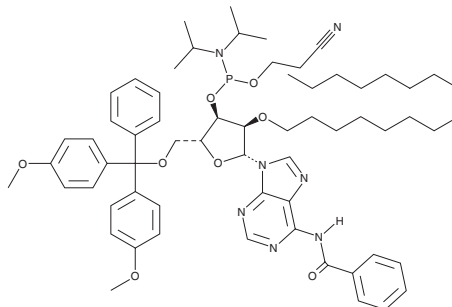
FW: 1,098.4

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

N⁶-Bz-5'-O-DMTr-2'-O-hexadecanyl adenosine 3'-CED phosphoramidite is supplied as a solid. A stock solution may be made by dissolving the N⁶-Bz-5'-O-DMTr-2'-O-hexadecanyl adenosine 3'-CED phosphoramidite in the solvent of choice, which should be purged with an inert gas. N⁶-Bz-5'-O-DMTr-2'-O-hexadecanyl adenosine 3'-CED phosphoramidite is sparingly soluble (1-10 mg/ml) in acetonitrile and DMSO.

Description

N⁶-Bz-5'-O-DMTr-2'-O-hexadecanyl adenosine 3'-CED phosphoramidite is a nucleoside building block containing a C16 alkyl chain at the 2'-position.

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, 06/10/2026

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