

PRODUCT INFORMATION



N⁶-Methyl-rA Phosphoramidite

Item No. 43600

CAS Registry No.: 588698-79-7

Formal Name: 3'-[2-cyanoethylN,N-bis(1-methylethyl)phosphoramidite]5'-O-[bis(4-methoxyphenyl)phenylmethyl]-2'-O-[(1,1-dimethylethyl)dimethylsilyl]-N-methyl-adenosine

MF: C₄₇H₆₄N₇O₇PSi

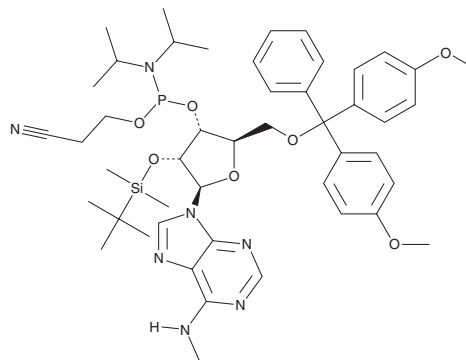
FW: 898.1

Purity: ≥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

N⁶-Methyl-rA phosphoramidite is supplied as a solid. A stock solution may be made by dissolving the N⁶-methyl-rA phosphoramidite in the solvent of choice, which should be purged with an inert gas. N⁶-Methyl-rA phosphoramidite is soluble (≥10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

N⁶-Methyl-rA phosphoramidite is a nucleoside building block that has been used in the synthesis of methylated oligonucleotides.¹

Reference

1. Dai, Q., Fong, R., Saikia, M., *et al.* Identification of recognition residues for ligation-based detection and quantitation of pseudouridine and N⁶-methyladenosine. *Nucleic Acids Res.* **35(18)**, 6322-6329 (2007).

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, 08/04/2025

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