

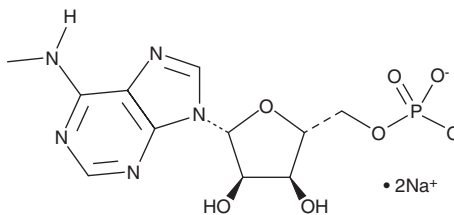
PRODUCT INFORMATION



N⁶-Methyladenosine 5'-monophosphate (sodium salt)

Item No. 23382

CAS Registry No.: 81921-35-9
Formal Name: N-methyl-5'-adenylic acid, disodium salt
Synonym: N⁶-Methyl-AMP
MF: C₁₁H₁₄N₅O₇P • 2Na
FW: 405.2
Purity: ≥95%
UV/Vis.: λ_{max}: 210, 266 nm
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

N⁶-Methyladenosine 5'-monophosphate (sodium salt) is supplied as a crystalline solid. Aqueous solutions of N⁶-methyladenosine 5'-monophosphate (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of N⁶-methyladenosine 5'-monophosphate (sodium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N⁶-Methyladenosine 5'-monophosphate is an activator of glycogen phosphorylase *b* ($K_a = 22 \mu\text{M}$).¹ It is also a non-competitive inhibitor of rat adenylate kinase II ($K_i = 4.2 \text{ mM}$).² N⁶-Methyladenosine 5'-monophosphate (sodium salt) has been used for immunoprecipitation of N⁶-methyladenosine.³

References

1. Morange, M., García Blanco, F., Vandebunder, B., *et al.* AMP Analogs : Their function in the activation of glycogen phosphorylase *b*. *Eur. J. Biochem.* **65**(2), 553-563 (1976).
2. Hai, T.T., Picker, D., Abo, M., *et al.* Species- or isozyme-specific enzyme inhibitors. 7. Selective effects in inhibitions of rat adenylate kinase isozymes by adenosine 5'-phosphate derivatives. *J. Med. Chem.* **25**(7), 806-812 (1982).
3. Fry, N.J., Law, B.A., Ilkayeva, O.R., *et al.* N⁶-methyladenosine is required for the hypoxic stabilization of specific mRNAs. *RNA* **23**(9), 1444-1455 (2017).

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, 12/19/2022

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