

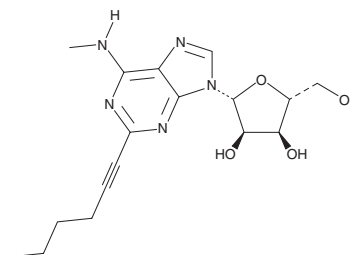
PRODUCT INFORMATION



HEMADO

Item No. 21015

CAS Registry No.: 403842-38-6
Formal Name: 2-(1-hexyn-1-yl)-N-methyl-adenosine
MF: C₁₇H₂₃N₅O₄
FW: 361.4
Purity: ≥98%
UV/Vis.: λ_{max}: 236, 274, 298 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

HEMADO is supplied as a crystalline solid. A stock solution may be made by dissolving the HEMADO in the solvent of choice. HEMADO is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of HEMADO in these solvents is approximately 20 and 5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of HEMADO can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of HEMADO in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

HEMADO is a selective agonist of the adenosine A₃ receptor (K_is = 1.1, 327, and 1,230 nM for A₃, A₁, and A_{2A} receptors, respectively).¹ HEMADO reduces release of lactate dehydrogenase (LDH) and creatine kinase-MB (CK-MB) from myocardium in isolated rat hearts subjected to ischemia-reperfusion injury.²

References

1. Volpini, R., Costanzi, S., Lambertucci, C., *et al.* N⁶-Alkyl-2-alkynyl derivatives of adenosine as potent and selective agonists at the human adenosine A₃ receptor and a starting point for searching A_{2B} ligands. *J. Med. Chem.* **45**(15), 3271-3279 (2002).
2. Amani, M., Jeddi, S., Ahmadiasl, N., *et al.* Effect of HEMADO on level of CK-MB and LDH enzymes after ischemia/reperfusion injury in isolated rat heart. *Bioimpacts* **3**(2), 101-104 (2013).

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, 09/30/2022

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