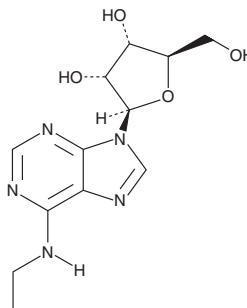


PRODUCT INFORMATION

N⁶-Ethyladenosine

Item No. 28428

CAS Registry No.: 14357-08-5
Formal Name: N-ethyl-adenosine
Synonym: NSC 516603
MF: C₁₂H₁₇N₅O₄
FW: 295.3
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 267 nm
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-Ethyladenosine is supplied as a solid. A stock solution may be made by dissolving the N⁶-ethyladenosine in the solvent of choice, which should be purged with an inert gas. N⁶-Ethyladenosine is soluble in the organic solvent DMSO at a concentration of approximately 75 mg/ml.

Description

N⁶-Ethyladenosine is an adenosine A₁ and A₃ receptor agonist.¹ It binds selectively to rat adenosine A₁ and human adenosine A₃ receptors over rat adenosine A_{2A} and A₃ receptors (K_is = 4.9, 4.7, 8,900 and 1,050 nM, respectively, for the recombinant receptors in CHO cells). N⁶-Ethyladenosine completely inhibits forskolin-induced cAMP accumulation in CHO cells expressing the human recombinant adenosine A₃ receptor when used at a concentration of 10 μM.

Reference

1. Gao, Z.-G., Blaustein, J.B., Gross, A.S., *et al.* N⁶-Substituted adenosine derivatives: Selectivity, efficacy, and species differences at A₃ adenosine receptors. *Biochem. Pharmacol.* **65**(10), 1675-1684 (2003).

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, 10/13/2022

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