

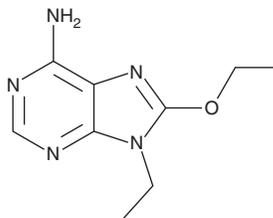
PRODUCT INFORMATION



ANR 94

Item No. 43476

CAS Registry No.: 634924-89-3
Formal Name: 8-ethoxy-9-ethyl-9H-purin-6-amine
Synonym: 8-ethoxy-9-Ethyladenine
MF: C₉H₁₃N₅O
FW: 207.2
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

ANR 94 is supplied as a solid. A stock solution may be made by dissolving the ANR 94 in the solvent of choice, which should be purged with an inert gas. ANR 94 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

ANR 94 is an adenosine A_{2A} receptor antagonist (K_i = 0.046 μM).¹ It is selective for the adenosine A_{2A} receptor over adenosine A₁, -A_{2B}, and -A₃ receptors (K_is = 2.4, >30, and 21 μM, respectively). ANR 94 (5 mg/kg) reduces the number of parkinsonian-like tremors induced by tacrine (Item No. 70240) in rats.² It decreases neurodegeneration and neuroinflammation in the substantia nigra pars compacta in a mouse model of MPTP-induced Parkinson's disease when administered at a dose of 0.5 mg/kg. ANR 94 (5 mg/kg) increases alcohol intake but not water or food intake in alcohol-preferring rats.³

References

1. Klotz, K.-N., Kachler, S., Lambertucci, C., *et al.* 9-Ethyladenine derivatives as adenosine receptor antagonists: 2- and 8-substitution results in distinct selectivities. *Naunyn Schmiedebergs Arch Pharmacol.* **367(6)**, 629-634 (2003).
2. Pinna, A., Tronci, E., Schintu, N., *et al.* A new ethyladenine antagonist of adenosine A_{2A} receptors: Behavioral and biochemical characterization as an antiparkinsonian drug. *Neuropharmacology* **58(3)**, 613-623 (2009).
3. Micioni Di Bonaventura, M.V., Cifani, C., Lambertucci, C., *et al.* Effects of A_{2A} adenosine receptor blockade or stimulation on alcohol intake in alcohol-preferring rats. *Psychopharmacol. (Berl)* **219(4)**, 945-957 (2011).

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, 07/02/2025

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