

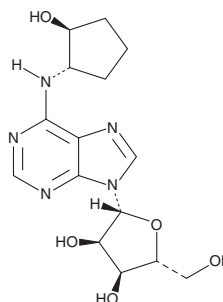
PRODUCT INFORMATION



GR79236

Item No. 28430

CAS Registry No.: 124555-18-6
Formal Name: N-[(1S,2S)-2-hydroxycyclopentyl]-adenosine
MF: $C_{15}H_{21}N_5O_5$
FW: 351.4
Purity: $\geq 98\%$
UV/Vis.: $\lambda_{\max}$: 270 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

GR79236 is supplied as a solid. A stock solution may be made by dissolving the GR79236 in the solvent of choice, which should be purged with an inert gas. GR79236 is soluble in organic solvents such as ethanol and DMSO. It is also soluble in water. The solubility of GR79236 in ethanol is approximately 15 mg/ml and approximately 30 mg/ml in DMSO and water. We do not recommend storing the aqueous solution for more than one day.

Description

GR79236 is a selective agonist of adenosine A_1 receptors (K_i s = 3.1 and 1,300 nM for adenosine A_1 and A_2 receptors, respectively, in rat brain).¹ GR79236 inhibits cAMP accumulation induced by isoprenaline (isoproterenol; Item No. 15592) in DDT1MF-2 hamster smooth muscle cells (IC_{50} = 2.6 nM) but only decreases contractility in isolated guinea pig atria with an EC_{50} value of 1,500 nM. It also inhibits lipolysis induced by norepinephrine (Item No. 16673) in rat adipocytes (EC_{50} = 0.83 nM).² GR79236 decreases locomotor activity and inhibits DMCM-induced seizures in mice (ED_{50} s = 0.13 and 0.3 mg/kg, respectively).¹ It also decreases mean blood pressure by 74% in anesthetized rats when administered at a dose of 0.1 mg/kg.

References

1. Knutsen, L.J.S., Lau, J., Petersen, H., *et al.* N-Substituted adenosines as novel neuroprotective A_1 agonists with diminished hypotensive effects. *J. Med. Chem.* **42**, 3463-3477 (1999).
2. Merkel, L.A., Hawkins, E.D., Colussi, D.J., *et al.* Cardiovascular and antilipolytic effects of the adenosine agonist GR79236. *Pharmacology* **51(4)**, 224-236 (1995).

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, 12/05/2022

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