

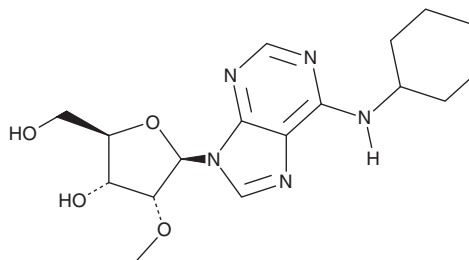
PRODUCT INFORMATION



SDZ-WAG 994

Item No. 43521

CAS Registry No.: 130714-47-5
Formal Name: N-cyclohexyl-2'-O-methyl-adenosine
MF: C₁₇H₂₅N₅O₄
FW: 363.4
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

SDZ-WAG 994 is supplied as a solid. A stock solution may be made by dissolving the SDZ-WAG 994 in the solvent of choice, which should be purged with an inert gas. SDZ-WAG 994 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

SDZ-WAG 994 is an adenosine A₁ receptor agonist.¹ It selectively binds to adenosine A₁ receptors over adenosine A₂ receptors (K_ds = 19.95 and 25,109 nM, respectively) and induces adenylate cyclase activity in rat adipocyte membranes (EC₅₀ = 398 nM). *In vivo*, SDZ-WAG 994 (0.14-1.4 μmol/kg) induces hypotension and bradycardia in spontaneously hypertensive rats.

Reference

1. Wagner, H., Milavec-Krizman, M., Gadiant, F., *et al.* General pharmacology of SDZ WAG 994, a potent selective and orally active adenosine A₁ receptor agonist. *Drug Dev. Res.* **34(3)**, 276-288 (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, 06/27/2025

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