

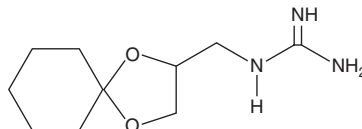
PRODUCT INFORMATION



Guanadrel

Item No. 33927

CAS Registry No.: 40580-59-4
Formal Name: N-(1,4-dioxaspiro[4.5]dec-2-ylmethyl)-guanidine
MF: $C_{10}H_{19}N_3O_2$
FW: 213.3
Purity: $\geq 95\%$
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

Guanadrel is supplied as a solid. A stock solution may be made by dissolving the guanadrel in the solvent of choice, which should be purged with an inert gas. Guanadrel is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of guanadrel in these solvents is approximately 30 mg/ml.

Description

Guanadrel is an antihypertensive agent.¹ *In vivo*, guanadrel (5 mg/kg) inhibits reflex increases in mean arterial and femoral perfusion blood pressures induced by bilateral carotid occlusion in dogs. Formulations containing guanadrel have been used in the treatment of hypertension.

Reference

1. Finnerty, F.A., Jr. and Brogden, R.N. Guanadrel. A review of its pharmacodynamic and pharmacokinetic properties and therapeutic use in hypertension. *Drugs* **30(1)**, 22-31 (1985).

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, 02/08/2023

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