PRODUCT INFORMATION



Guanoclor

Item No. 27659

CAS Registry No.: 5001-32-1

Formal Name: 2-[2-(2,6-dichlorophenoxy)ethyl]-

hydrazinecarboximidamide

Synonym: Guanochlor MF: $C_9H_{12}CI_2N_4O$

263.1 FW: ≥95% **Purity:**

Supplied as: A crystalline solid

Storage: -20°C Stability: ≥4 years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Guanoclor is an antihypertensive agent that binds to non-adrenergic sites in pig kidney membranes (IC $_{50}$ = 10 nM). 1,2 It inhibits the Na $^+$ /H $^+$ exchanger in pig kidney membranes (IC $_{50}$ = 1 μ M) and cell-free bovine dopamine β -hydroxylase (IC₅₀ = 90 μ M). Guanoclor inhibits electrically stimulated norepinephrine release from spleen.¹ It decreases norepinephrine levels in dog heart, spleen, hypothalamus, and adrenal gland when administered at a dose of 10 mg/kg per day for six weeks. It decreases adrenal catecholamine levels in rats when administered at a dose of 20 mg/kg per day for six weeks. Guanoclor (20 mg/kg) decreases experimentally induced hypertension and the tyramine-induced pressor response and increases the norepinephrine-induced pressor response in non-anesthetized dogs.

References

- 1. Lawrie, T.D., Lorimer, A.R., McAlpine, S.G., et al. Clinical trial and pharmacological study of compound 1029. ("Vatensol"). Br. Med. J. 1(5380), 402-406 (1964).
- 2. Vigne, P., Lazdunski, M., and Frelin, C. Guanabenz, guanochlor, guanoxan and idazoxan bind with high affinity to non-adrenergic sites in pig kidney membranes. Eur. J. Pharmacol. 160(2), 295-298 (1989).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information Buyer agrees to purchase the mater can be found on our website.

Copyright Cayman Chemical Company, 10/17/2022

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