

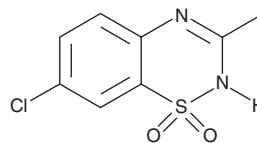
PRODUCT INFORMATION



Diazoxide

Item No. 14576

CAS Registry No.: 364-98-7
Formal Name: 7-chloro-3-methyl-1,1-dioxide-2H-1,2,4-benzothiadiazine
Synonyms: NSC 64198, NSC 76130, SRG 95213
MF: C₈H₇ClN₂O₂S
FW: 230.7
Purity: ≥99%
UV/Vis.: λ_{max}: 267 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

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

Description

Diazoxide is an activator of sulfonylurea receptor 1 (SUR1) linked to ATP-sensitive potassium channel K_{ir}6.2 (EC₅₀ = 14.1 μM in a FLIPR assay using HEK293 cells).¹ It also activates SUR2A/K_{ir}6.2 and SUR2B/K_{ir}6.2 channels in HEK293T cells in a patch-clamp assay when used at concentrations of 30 and 300 μM.² Diazoxide inhibits glucose-induced insulin release from isolated rat pancreatic β cells and induces relaxation of isolated rat aortic rings precontracted with potassium chloride (IC₅₀s = 22.6 and 22.4 μM, respectively).³ It reduces mean arterial pressure and cerebral blood flow in spontaneously hypertensive rats when administered intravenously as a 5 mg/kg bolus dose.⁴ Diazoxide (50 mg/kg, i.p.) increases blood glucose levels in mice.⁵ Formulations containing diazoxide have been used in the treatment of hypoglycemia.

References

1. Gopalakrishnan, M., Molinari, E.J., Char-Change, S., *et al.* Pharmacology of human sulphonylurea receptor SUR1 and inward rectifier K⁺ channel Kir6.2 combination expressed in HEK-293 cells. *Br. J. Pharmacol.* **129(7)**, 1323-1332 (2000).
2. Matsuoka, T., Matsushita, K., Katayama, Y., *et al.* C-Terminal tails of sulfonylurea receptors control ADP-induced activation and diazoxide modulation of ATP-sensitive K⁺ channels. *Circ. Res.* **87(10)**, 873-880 (2000).
3. de Tullio, P., Becker, B., Boverie, S., *et al.* Toward tissue-selective pancreatic B-cells K_{ATP} channel openers belonging to 3-alkylamino-7-halo-4H-1,2,4-benzothiadiazine 1,1-dioxides. *J. Med. Chem.* **46(15)**, 3342-3353 (2003).
4. Barry, D.I., Strandgaard, S., Graham, D.I., *et al.* Effect of diazoxide-induced hypotension on cerebral blood flow in hypertensive rats. *Eur. J. Clin. Invest.* **13(3)**, 201-207 (1983).
5. Foy, J.M. and Furman, B.L. Effect of single dose administration of diuretics on the blood sugar of alloxan-diabetic mice or mice made hyperglycaemic by the acute administration of diazoxide. *Br. J. Pharmacol.* **47(1)**, 124-132 (1973).

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, 01/23/2020

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