

Diazoxide

K⁺ channel activator

ATP-sensitive K⁺ channel activator. Effective concentrations in patch clamp experiments: 100-600µM. Blocks desensitization of AMPA receptors. Acts in pancreatic β-cells, where it is more potent than pinacidil (Prod. No. ALX-550-285). Antihypertensive and vasodilatory agent.

Citations: 8

[View Online »](#)

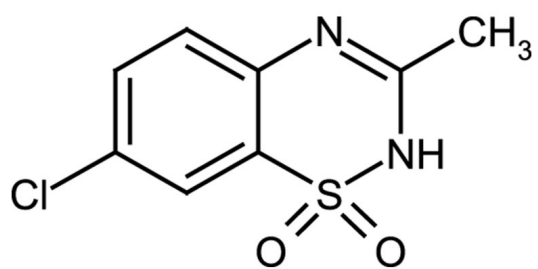
Ordering Information

[Order Online »](#)

ALX-550-260-M500	500mg
ALX-550-260-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 7-Chloro-3-methyl-2H-1,2,4-benzothiadiazine 1,1-dioxide

Appearance White to off-white powder.

CAS 364-98-7

Couple Target Potassium channel

Couple Type Activator

Formula $C_8H_7ClN_2O_2S$

MI 14: 3004

MW 230.7

Purity ≥98% (HPLC)

RTECS DK8185000

Solubility Soluble in DMSO (25mg/ml); insoluble in water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)