

# Diazoxide

K<sup>+</sup> channel activator

ATP-sensitive K<sup>+</sup> 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 »](#)

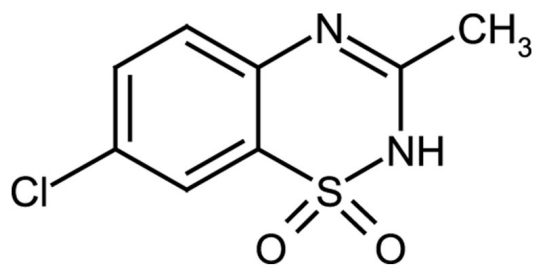
## Ordering Information

[Order Online »](#)

ALX-550-260-M500	500mg
ALX-550-260-46190	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)