

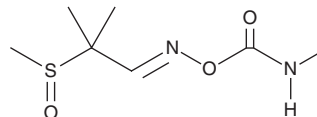
PRODUCT INFORMATION



Aldicarb Sulfoxide

Item No. 24211

CAS Registry No.: 1646-87-3
Formal Name: O-[(methylamino)carbonyl]oxime 2-methyl-2-(methylsulfinyl)-propanal
MF: C₇H₁₄N₂O₃S
FW: 206.3
Purity: ≥98%
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

Aldicarb sulfoxide is supplied as a solid. A stock solution may be made by dissolving the aldicarb sulfoxide in water. We do not recommend storing the aqueous solution for more than one day.

Description

Aldicarb sulfoxide is a metabolite and degradation product of the carbamate pesticide aldicarb (Item No. 18466).¹ Aldicarb sulfoxide (10 µg/ml) reduces the number of male *T. semipenetrans* second-stage larvae that reach the third, fourth, and adult stages of development. It inhibits carboxylesterase and cholinesterase in zebrafish (IC₅₀ = 10 µM for both).²

References

1. Huang, S.P. and Van Gundy, S.D. Effects of aldicarb and its sulfoxide and sulfone on the biology of *Tylenchulus semipenetrans*. *J. Nematol.* **10(1)**, 100-106 (1978).
2. Küster, E. and Altenburger, R. Suborganismic and organismic effects of aldicarb and its metabolite aldicarb-sulfoxide to the zebrafish embryo (*Danio rerio*). *Chemosphere* **68(4)**, 751-760 (2007).

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, 12/06/2022

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