

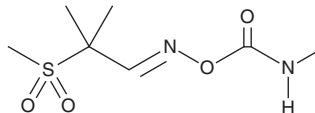
# PRODUCT INFORMATION



## Aldicarb sulfone

Item No. 24210

CAS Registry No.: 1646-88-4  
Formal Name: O-[(methylamino)carbonyl]oxime  
2-methyl-2-(methylsulfonyl)-propanal  
MF: C<sub>7</sub>H<sub>14</sub>N<sub>2</sub>O<sub>4</sub>S  
FW: 222.3  
Purity: ≥95%  
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 sulfone is supplied as a solid. A stock solution may be made by dissolving the aldicarb sulfone in the solvent of choice, which should be purged with an inert gas. Aldicarb sulfone is soluble in the organic solvent DMSO at a concentration of approximately 34 mg/ml.

### Description

Aldicarb sulfone is a metabolite and degradation product of the carbamate pesticide aldicarb (Item No. 18466).<sup>1</sup> Aldicarb sulfone (100 µg/ml) reduces the number of male *T. semipenetrans* second-stage larvae that reach the third, fourth, and adult stages of development.

### Reference

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).

#### 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, 11/16/2022

#### 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