

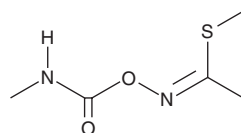
PRODUCT INFORMATION



Methomyl

Item No. 25610

CAS Registry No.: 16752-77-5
Formal Name: N-[[[(methylamino)carbonyl]oxy]-ethanimidothioic acid, methyl ester
MF: C₅H₁₀N₂O₂S
FW: 162.2
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

Methomyl is supplied as a solid. A stock solution may be made by dissolving the methomyl in the solvent of choice, which should be purged with an inert gas. Methomyl is slightly soluble in chloroform and methanol.

Description

Methomyl is a carbamate insecticide and molluscicide.^{1,2} It induces mortality of tobacco budworm larvae when applied at a concentration of 0.56 kg/hectare and southern armyworm larvae at a concentration of 0.01 and 100 ppm.³ Methomyl is toxic to the land snail *M. obstructa* (LD₅₀ = 11.9 µg/snail).² It is also toxic to rats *via* oral but not dermal administration (LD₅₀s = 30 and >2,000 mg/kg).¹ Formulations containing methomyl have been used in the control of insects in agriculture.

References

1. European Food Safety Authority. Conclusion regarding the peer review of the pesticide risk assessment of the active substance methomyl. *EFSA Scientific Report* **222**, 1-99 (2008).
2. Hussein, H.I., Al-Rajhy, D., El-Shahawi, F.I., *et al.* Molluscicidal activity of *Pergularia tomentosa* (L.), methomyl and methiocarb, against land snails. *Int. J. Pest Mgt.* **45(3)**, 211-213 (2010).
3. Wolfenbarger, D.A. and Redfern, R.E. Toxicity of five carbamate insecticides to the two-spotted spider mite and larvae of the southern armyworm and the tobacco budworm. *J. Econ. Entomol.* **61(2)**, 580-581 (1968).

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, 09/12/2023

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