

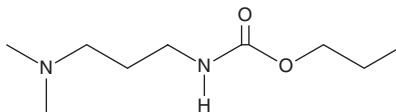
PRODUCT INFORMATION



Propamocarb

Item No. 39502

CAS Registry No.: 24579-73-5
Formal Name: N-[3-(dimethylamino)propyl]-
carbamic acid, propyl ester
Synonyms: Nor-Am 39744, SN 39744
MF: $C_9H_{20}N_2O_2$
FW: 188.3
Purity: $\geq 98\%$
Supplied as: A neat oil
Storage: -20°C
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Propamocarb is supplied as a neat oil. A stock solution may be made by dissolving the propamocarb in the solvent of choice, which should be purged with an inert gas. Propamocarb is soluble in the organic solvent methanol.

Description

Propamocarb is a carbamate fungicide.¹ It is active against the plant pathogenic fungi *P. aphanidermatum*, *P. dissotocum*, *P. irregulare*, and *P. ultimum* when used at a concentration of 400 $\mu\text{g/ml}$.² Propamocarb potentiates estrogen receptor-mediated transactivation induced by 17 β -estradiol (Item No. 10006315) in MCF-7 cells in a reporter assay but has no effect when used alone.¹ It has been found in spring water.³

References

1. Andersen, H.R., Vinggaard, A.M., Rasmussen, T.H., *et al.* Effects of currently used pesticides in assays for estrogenicity, androgenicity, and aromatase activity *in vitro*. *Toxicol. Appl. Pharmacol.* **179**(1), 1-12 (2002).
2. Moorman, G.W., and Kim, S.H. Species of *Pythium* from greenhouses in Pennsylvania exhibit resistance to propamocarb and mefenoxam. *Plant Dis.* **88**(6), 630-632 (2004).
3. Nantia, E.A., Moreno-González, D., Manfo, F.P.T., *et al.* Characterization of carbamate pesticides in natural water from Cameroon. *Anal. Lett.* **50**(9), 1397-1409 (2017).

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, 07/09/2025

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