

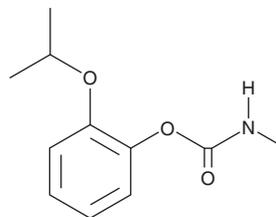
PRODUCT INFORMATION



Propoxur

Item No. 25628

CAS Registry No.: 114-26-1
Formal Name: 2-(1-methylethoxy)-phenol 1-(N-methylcarbamate)
Synonyms: ENT 25,671, NSC 379584
MF: C₁₁H₁₅NO₃
FW: 209.2
Purity: ≥95%
UV/Vis.: λ_{max}: 215, 272 nm
Supplied as: A crystalline 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

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

Description

Propoxur is a carbamate insecticide, an inhibitor of acetylcholinesterase (IC₅₀ = 4.3 μM), and an antagonist of α4β4 subunit-containing nicotinic acetylcholine receptors (IC₅₀ = 165 μM for the rat receptor).^{1,2} It is lethal to German cockroaches (*B. germanica*) with LD₅₀ values of 2.45, 4.73, and 1.06 μg per insect for newly emerged males, newly emerged females, and 4-week-old nymphs, respectively. Propoxur induces lactate dehydrogenase (LDH) release from and inhibits the growth of flounder gill cells (IC₅₀s = 86.59 and 89.96 μg/ml, respectively) via induction of necrosis.³ It also induces yolk sac and pericardial edema in zebrafish embryos when used at concentrations of 100 and 200 μg/ml. Propoxur (8.3 mg/kg) inhibits rat blood and brain cholinesterase, decreases ambulation and rearing in an open field test, and increases latency to escape foot shock after an audio stimulus in the two-way active avoidance box in rats.⁴ Formulations containing propoxur have been used in the control of insects in agriculture, as well as residential areas.

References

1. Qian, K., Wei, X., Zeng, X., *et al.* Stage-dependent tolerance of the German cockroach, *Blattella germanica* for dichlorvos and propoxur. *J. Insect Sci.* **10**(1), 201 (2010).
2. Smulders, C.J., Bueters, T.J., Van Kleef, R.G., *et al.* Selective effects of carbamate pesticides on rat neuronal nicotinic acetylcholine receptors and rat brain acetylcholinesterase. *Toxicol. Appl. Pharmacol.* **193**(2), 139-146 (2003).
3. Pandey, M.R. and Guo, H. Evaluation of cytotoxicity, genotoxicity and embryotoxicity of insecticide propoxur using flounder gill (FG) cells and zebrafish embryos. *Toxicol. In Vitro* **28**(3), 340-353 (2014).
4. Thiesen, F.V., Barros, H.M.T., Tannhauser, M., *et al.* Behavioral changes and cholinesterase activity of rats acutely treated with propoxur. *Jpn. J. Pharmacol.* **79**(1), 25-31 (1999).

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, 09/18/2023

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