

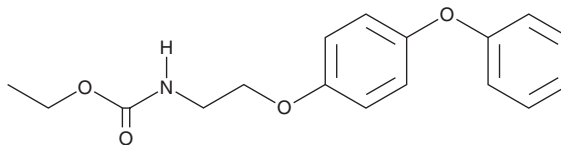
PRODUCT INFORMATION



Fenoxycarb

Item No. 25607

CAS Registry No.: 72490-01-8
Formal Name: N-[2-(4-phenoxyphenoxy)ethyl]-
carbamic acid, ethyl ester
Synonyms: ABG 6215, Ro 13-5223
MF: C₁₇H₁₉NO₄
FW: 301.3
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

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

Description

Fenoxycarb is a non-neurotoxic carbamate insecticide that acts as an insect growth regulator via juvenile hormone-like activity.¹ It inhibits terminal development of first instar and newly transformed second instar nymphs of Florida red scale (*C. aonidum*) when used at a concentration of 0.0125% AI.² Fenoxycarb (5 and 10 mg AI/colony) reduces the colony size index of laboratory colonies of red imported fire ants (*S. invicta*) by 93.6 to 95.9% at 8 weeks post-treatment.³ It is toxic to *D. magna* (LC₅₀ = 0.5 mg a.s./L) and fish including *O. mykiss*, *L. macrochirus*, *C. carpio*, *I. punctatus*, and *C. variegatus* (LC₅₀s = 0.66-1.5 mg a.s./L), but is not toxic to rats (LD₅₀ = >10,000 mg/kg).⁴ Fenoxycarb is also an antagonist of α4β4-, α4β2-, α3β4-, and α3β2-containing rat nicotinic acetylcholine receptors (nAChRs; IC₅₀s = 3, 2.4, 1.8, and 7.6 μM, respectively) but not rat brain acetylcholinesterase (AChE; IC₅₀ = >1,000 μM).⁵

References

1. Grenier, S. and Grenier, A.M. Fenoxycarb, a fairly new insect growth regulator: A review of its effects on insects. *Ann. Appl. Biol.* **122**(2), 369-403 (1993).
2. Peleg, B.A. Effect of a new insect growth regulator, RO 13-5223, on hymenopterous parasites of scale insects. *Phytoparasitica* **10**(1), 27-31 (1983).
3. Banks, W.A., Williams, D.F., and Lofgren, C.S. Effectiveness of fenoxycarb for control of red imported fire ants (Hymenoptera: Formicidae). *J. Econ. Entomol.* **81**(1), 83-87 (1988).
4. European Food Safety Authority. Conclusion on the peer review of the pesticide risk assessment of the active substance fenoxycarb. *EFSA J.* **8**(12), 1779 (2010).
5. 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).

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, 10/06/2022\

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