

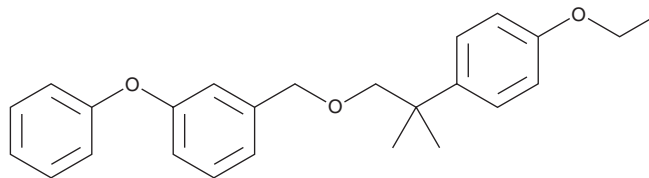
PRODUCT INFORMATION



Etofenprox

Item No. 25785

CAS Registry No.: 80844-07-1
Formal Name: 1-[[2-(4-ethoxyphenyl)-2-methylpropoxy]methyl]-3-phenoxy-benzene
Synonym: SA 130301
MF: C₂₅H₂₈O₃
FW: 376.5
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

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

Description

Etofenprox is a non-ester pyrethroid insecticide.¹ It induces toxicity in the housefly *M. domestica* (LD₅₀ = 23 ng/fly). Etofenprox (0.005-0.02%) induces 100% mortality in green leafhopper adults and brown planthopper and white-backed planthopper nymphs on rice plants within 24 hours of application.² It is toxic to the aquatic invertebrate *D. magna* (EC₅₀ = 1.2 µg a.s./L) and the fish *O. mykiss* (LC₅₀ = 2.7 µg a.s./L) but not rats (LD₅₀ = >2,000 mg/kg).³ Formulations containing etofenprox have been used in the control of agricultural pests.

References

1. Pedersen, L.-E.K. The potency of cyclopropane pyrethroid ethers against susceptible and resistant strains of the house fly *Musca domestica*. *Experientia* **42(9)**, 1057-1058 (1986).
2. Krishnaiah, N.V. and Kalode, M.B. Efficacy of ethofenprox against leafhoppers and planthoppers, with emphasis on preventing resurgence of brown planthopper, *Nilaparvata lugens* (Stål), in rice. *Crop Prot.* **12(7)**, 532-538 (1993).
3. European Food Safety Authority. Conclusion regarding the peer review of the pesticide risk assessment of the active substance etofenprox. *EFSA Scientific Report* **213**, 1-131 (2008).

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, 12/22/2022

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