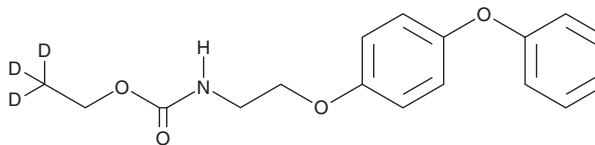


PRODUCT INFORMATION



Fenoxycarb-d₃ Item No. 39598

CAS Registry No.: 2468639-22-5
Formal Name: N-[2-(4-phenoxyphenoxy)ethyl]-
carbamic acid, ethyl-2,2,2-d₃ ester
MF: C₁₇H₁₆D₃NO₄
FW: 304.4
Chemical Purity: ≥95% (Fenoxycarb)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
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-d₃ is intended for use as an internal standard for the quantification of fenoxycarb (Item No. 25607) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Fenoxycarb-d₃ is supplied as a solid. A stock solution may be made by dissolving the fenoxycarb-d₃ in the solvent of choice, which should be purged with an inert gas. Fenoxycarb-d₃ is slightly soluble in chloroform and methanol.

Description

Fenoxycarb-d₃ is intended for use as an internal standard for the quantification of fenoxycarb (Item No. 25607) by GC- or LC-MS. Fenoxycarb is a carbamate insect growth regulator.^{1,2} It inhibits embryonic and larval development in cat flea (*C. felis*) eggs and larvae, respectively, when applied to filter paper discs at a concentration of 1 µg/cm².¹ Fenoxycarb (1% w/w) decreases the number of worker red imported fire ants (*S. invicta*) by 95.9% after eight weeks when provided as bait.² It is toxic to *D. magna* (LC₅₀ = 0.5 mg/L) and fish, including *O. mykiss*, *L. macrochirus*, *C. carpio*, *I. punctatus*, and *C. variegatus* (LC₅₀s = 0.66-1.5 mg/L), but is not toxic to rats (LD₅₀ = >10,000 mg/kg).³ Fenoxycarb is also an antagonist of α4β4-, α4β2-, α3β4-, and α3β2 subunit-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).¹

References

1. Marchiondo, A.A., Riner, J.L., Sonenshine, D.E., *et al.* Ovicidal and larvicidal modes of action of fenoxycarb against the cat flea (Siphonaptera: Pulicidae). *J. Med. Entomol.* **27**(5), 913-921 (1990).
2. 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).
3. 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).
4. 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, 07/14/2025

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