

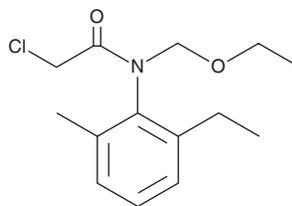
PRODUCT INFORMATION



Acetochlor

Item No. 24131

CAS Registry No.: 34256-82-1
Formal Name: 2-chloro-N-(ethoxymethyl)-N-(2-ethyl-6-methylphenyl)-acetamide
MF: C₁₄H₂₀ClNO₂
FW: 269.8
Purity: ≥95%
Supplied as: An 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

Acetochlor is supplied as an oil. A stock solution may be made by dissolving the acetochlor in the solvent of choice, which should be purged with an inert gas. Acetochlor is slightly soluble in chloroform, DMSO, and water.

Description

Acetochlor is an herbicide that completely inhibits shoot growth of oat (*A. sativa*) and mustard (*S. alba*) plants when used at a concentration of 2 kg/hectare.¹ It accelerates thyroid hormone-induced metamorphosis of and alters thyroid hormone-responsive gene expression in *X. laevis*. *In vivo*, acetochlor induces formation of thyroid, bone, stomach, and nasal tumors in rats as well as liver and lung tumors in mice.³ Acetochlor also induces pericardial edema, thrombosis, circulation defects, and a reduction in the number of cardiomyocytes in zebrafish larvae when administered in tank water at the maximum non-lethal concentration (MNL) of 9.6 µg/ml.⁴ Formulations containing acetochlor have been used as herbicides in agricultural settings.

References

1. Jablonkai, I. Alkylating reactivity and herbicidal activity of chloroacetamides. *Pest. Manag. Sci.* **59(4)**, 443-450 (2003).
2. Crump, D., Werry, K., Veldhoen, N., et al. Exposure to the herbicide acetochlor alters thyroid hormone-dependent gene expression and metamorphosis in *Xenopus Laevis*. *Environ. Health Perspect.* **110(12)**, 1199-1205 (2002).
3. Hurley, P.M. Mode of carcinogenic action of pesticides inducing thyroid follicular cell tumors in rodents. *Environ. Health Perspect.* **106(8)**, 437-445 (1998).
4. Liu, H., Chu, T., Chen, L., et al. *In vivo* cardiovascular toxicity induced by acetochlor in zebrafish larvae. *Chemosphere* **181**, 600-608 (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, 07/25/2023

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