

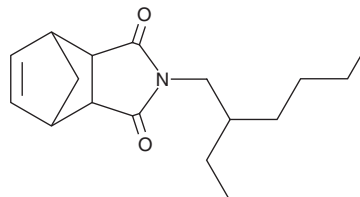
PRODUCT INFORMATION



MGK-264

Item No. 25819

CAS Registry No.: 113-48-4
Formal Name: 2-(2-ethylhexyl)-3a,4,7,7a-tetrahydro-4,7-methano-1H-isoindole-1,3(2H)-dione
Synonyms: ENT 8184, NSC 36678, NSC 406879, Octyl Bicycloheptene Dicarboximide
MF: C₁₇H₂₅NO₂
FW: 275.4
Purity: ≥95%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

MGK-264 is supplied as a liquid. A stock solution may be made by dissolving the MGK-264 in the solvent of choice, which should be purged with an inert gas. MGK-264 is slightly soluble in chloroform, methanol, and DMSO.

Description

MGK-264 is a dicarboxamide synergist used to enhance the effects of pesticides.¹ It is not lethal to house flies when used at a concentration of 5,000 ppm and induces only 26.4% mortality of two-spotted spider mites when used at a concentration of 1,000 ppm. MGK-264 has synergistic effects on toxicity when used in combination with the insecticide cyhexatin in susceptible strains of *S. littoralis* (LD₅₀S = 0.98 and 40 µg/larva with or without MGK-264, respectively) and in field strains (LD₅₀S = 0.15 and 1.1 µg/larva with or without MGK-264, respectively).² It also decreases the fecundity of adult *L. acuminata* and reduces survival of the offspring after hatching when used in combination with sublethal doses of a variety of plant-derived molluscicides.³ Formulations containing MGK-264 have been used in the control of insects in agricultural, aquatic, commercial, industrial, and residential areas.

References

1. Kilsheimer, J.A. and Kaufman, H.A. Pesticidal composition comprising a benzothienyl carbamate and synergist. *Rhone-Poulenc, Inc. US3689658A* (1972).
2. Radwan, H.S.A., Riskallah, M.R., and El-Keie, I.A. Synergistic effects on the toxicity of organotins on cotton leafworms. *Toxicology* **14(3)**, 193-198 (1979).
3. Singh, K. and Singh, D.K. Effect of different combinations of MGK-264 or piperonyl butoxide with plant-derived molluscicides on snail reproduction. *Arch. Environ. Contam. Toxicol.* **38(2)**, 182-190 (2000).

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, 07/26/2023

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