

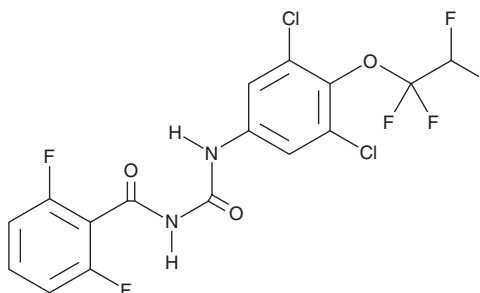
PRODUCT INFORMATION



Hexaflumuron

Item No. 18598

CAS Registry No.: 86479-06-3
Formal Name: N-[[[3,5-dichloro-4-(1,1,2,2-tetrafluoroethoxy)phenyl]amino]carbonyl]-2,6-difluorobenzamide
Synonyms: AI-3-29832, OMS 3031, XRD-473
MF: C₁₆H₈Cl₂F₆N₂O₃
FW: 461.1
Purity: ≥95%
UV/Vis.: λ_{max}: 209, 254 nm
Supplied as: A crystalline 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

Hexaflumuron is supplied as a crystalline solid. A stock solution may be made by dissolving the hexaflumuron in the solvent of choice, which should be purged with an inert gas. Hexaflumuron is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of hexaflumuron in these solvents is approximately 10 mg/ml.

Hexaflumuron is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, hexaflumuron should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Hexaflumuron has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Hexaflumuron is a chitin synthesis inhibitor used to bait and eliminate termite colonies.¹ Hexaflumuron spreads efficiently through entire populations through mutual exchange of liquids *via* regurgitation.² Termites are unable to metabolize hexaflumuron and clearance is slow, resulting in up to 100% elimination. Hexaflumuron has also been tested for use with the raisin moth, cowpea weevil, and Asiatic rice borer with positive results.^{3,4} Formulations containing hexaflumuron have been used in the control of insects in residential areas.

References

1. Evans, T.A. and Iqbal, N. Termite (order Blattodea, infraorder Isoptera) baiting 20 years after commercial release. *Pest. Manag. Sci.* **71**(7), 897-906 (2015).
2. Sheets, J.J., Karr, L.L., and Dripps, J.E. Kinetics of uptake, clearance, transfer, and metabolism of hexaflumuron by eastern subterranean termites (Isoptera: Rhinotermitidae). *J. Econ. Entomol.* **93**(3), 871-877 (2000).
3. Khajepour, S., Izadi, H., and Asari, M.J. Evaluation of two formulated chitin synthesis inhibitors, hexaflumuron and lufenuron against the raisin moth, *Ephestia figulilella*. *J. Insect. Sci.* **12**, 102 (2012).
4. Abo-Elghar, G.E., El-Sheikh, A.E., El-Sayed, F.M., et al. Persistence and residual activity of an organophosphate, pirimiphos-methyl, and three IGRs, hexaflumuron, teflubenzuron and pyriproxyfen, against the cowpea weevil, *Callosobruchus maculatus* (Coleoptera: Bruchidae). *Pest. Manag. Sci.* **60**(1), 95-102 (2004).

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