

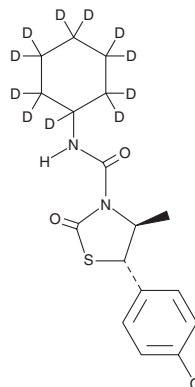
PRODUCT INFORMATION



Hexythiazox-d₁₁

Item No. 39602

Formal Name: (4R,5R)-rel-5-(4-chlorophenyl)-N-(cyclohexyl-d₁₁)-4-methyl-2-oxo-3-thiazolidinecarboxamide
MF: C₁₇H₁₀ClD₁₁N₂O₂S
FW: 363.9
Chemical Purity: ≥95% (Hexythiazox)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₁₁); ≤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

Hexythiazox-d₁₁ is intended for use as an internal standard for the quantification of hexythiazox (Item No. 25626) 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).

Hexythiazox-d₁₁ is supplied as a solid. A stock solution may be made by dissolving the hexythiazox-d₁₁ in the solvent of choice, which should be purged with an inert gas. Hexythiazox-d₁₁ is soluble in DMSO and methanol.

Description

Hexythiazox is an acaricide.^{1,2} It induces toxicity in the summer and winter eggs of the European spider mite (*P. ulmi*; LC₅₀s = 2.2 and 20 ppm, respectively).² Hexythiazox is toxic to bluegill (*L. macrochirus*; LC₅₀ = 3.2 mg/L) and *D. magna* (EC₅₀ = 0.36 mg/L) but not rats (LD₅₀ = >5,000 mg/kg).² Formulations containing hexythiazox have been used in the control of mites in agriculture.

References

1. Nauen, R., Stumpf, N., Elbert, A., *et al.* Acaricide toxicity and resistance in larvae of different strains of *Tetranychus urticae* and *Panonychus ulmi* (Acari: Tetranychidae). *Pest Manag. Sci.* **57**(3), (2001).
2. European Food Safety Authority. Conclusion on the peer review of the pesticide risk assessment of the active substance hexythiazox. *EFSA J.* **8**(10), 1722 (2010).

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, 07/17/2025

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