

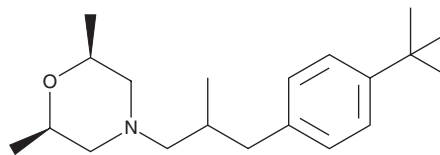
PRODUCT INFORMATION



Fenpropimorph

Item No. 20875

CAS Registry No.: 67564-91-4
Formal Name: (2R,6S)-rel-4-[3-[4-(1,1-dimethylethyl)phenyl]-2-methylpropyl]-2,6-dimethylmorpholine
Synonym: BAS 42100F
MF: C₂₀H₃₃NO
FW: 303.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 263, 271 nm
Supplied as: A neat 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

Fenpropimorph is supplied as a neat oil. A stock solution may be made by dissolving the fenpropimorph in the solvent of choice, which should be purged with an inert gas. Fenpropimorph is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fenpropimorph in ethanol and DMF is approximately 15 mg/ml and approximately 5 mg/ml in DMSO.

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

Description

Fenpropimorph is a fungicide that inhibits the sterol pathway.¹ It inhibits Δ^8 - Δ^7 -sterol isomerase in yeast at low concentrations, with Δ^{14} -sterol reductase being blocked at higher levels, preventing the biosynthesis of ergosterol (Item No. 19850).² Fenpropimorph is also able to inhibit sterol synthesis in certain plants and mammalian cells.^{1,3} Formulations containing fenpropimorph have been used in the control of fungi in agriculture.

References

1. Costet, M.F., Achouri, M.E., Charlet, M., *et al.* Ecdysteroid biosynthesis and embryonic development are disturbed in insects (*Locusta migratoria*) reared on plant diet (*Triticum sativum*) with a selectively modified sterol profile. *Proc. Natl. Acad. Sci. USA* **84**(3), 643-647 (1987).
2. Marcireau, C., Guilloton, M., and Karst, F. *In vivo* effects of fenpropimorph on the yeast *Saccharomyces cerevisiae* and determination of the molecular basis of the antifungal property. *Antimicrob. Agents Chemother.* **34**(6), 989-993 (1990).
3. Corio-Costet, M.F., Gerst, N., Benveniste, P., *et al.* Inhibition by the fungicide fenpropimorph of cholesterol biosynthesis in 3T3 fibroblasts. *Biochem J.* **256**(3), 829-834 (1988).

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