

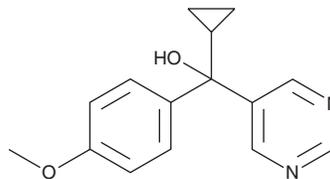
PRODUCT INFORMATION



Ancymidol

Item No. 33311

CAS Registry No.: 12771-68-5
Formal Name: α -cyclopropyl- α -(4-methoxyphenyl)-5-pyrimidinemethanol
Synonym: EL 531
MF: C₁₅H₁₆N₂O₂
FW: 256.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 224 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Ancymidol is supplied as a crystalline solid. A stock solution may be made by dissolving the ancymidol in the solvent of choice, which should be purged with an inert gas. Ancymidol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ancymidol in these solvents is approximately 1 mg/ml in ethanol and approximately 30 mg/ml in DMSO and DMF.

Ancymidol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ancymidol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ancymidol has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ancymidol is a plant growth regulator.¹ It inhibits oxidation of the gibberellin biosynthetic intermediates *ent*-kaur-16-ene, *ent*-kaur-16-en-19-ol, and *ent*-kaur-16-en-19-al, but not *ent*-kaur-16-en-19-oic acid, in immature *M. macrocarpus* seed endosperm preparations.² Ancymidol slows shoot and root elongation of *Phaseolus vulgaris* when used at concentrations ranging from 0.316 to 316 $\mu\text{g}/\text{ml}$.¹ Formulations containing ancymidol have been used as plant growth regulators in aquatic areas.

References

1. Shive, J.B. and Sisler, H.D. Effects of ancymidol (a growth retardant) and triarimol (a fungicide) on the growth, sterols, and gibberellins of *Phaseolus vulgaris* (L.). *Plant Physiol.* **57(4)**, 640-644 (1976).
2. Coolbaugh, R.C., Hirano, S.S., and West, C.A. Studies on the specificity and site of action of α -cyclopropyl- α -[*p*-methoxyphenyl]-5-pyrimidine methyl alcohol (ancymidol), a plant growth regulator. *Plant Physiol.* **62(4)**, 571-576 (1978).

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