

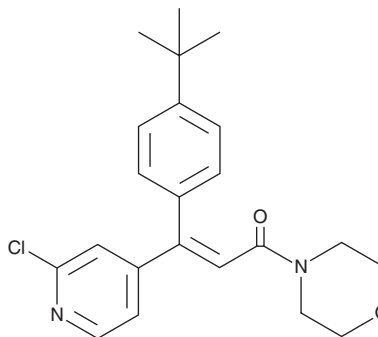
PRODUCT INFORMATION



Pyrimorph

Item No. 37286

CAS Registry No.: 868390-90-3
Formal Name: 3-(2-chloro-4-pyridinyl)-3-[4-(1,1-dimethylethyl)phenyl]-1-(4-morpholinyl)-2-propen-1-one
MF: C₂₂H₂₅ClN₂O₂
FW: 384.9
Purity: ≥90%
UV/Vis.: λ_{max}: 244 nm
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

Pyrimorph is supplied as a solid. A stock solution may be made by dissolving the pyrimorph in the solvent of choice, which should be purged with an inert gas. Pyrimorph is soluble in the organic solvent dimethyl formamide at a concentration of approximately 1 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of pyrimorph can be prepared by directly dissolving the solid in aqueous buffers. The solubility of pyrimorph in PBS (pH 7.2) is approximately 0.20 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Pyrimorph is a fungicide.¹ It is active against the plant pathogenic fungi *P. infestans*, *P. capsica*, and *R. solani* (EC₅₀s = 1.36, 0.72, and 4.44 µg/ml, respectively). Pyrimorph is toxic to zebrafish (LC₅₀ = 19.8 mg/L).²

References

1. Chen, X., Yuan, H., Qin, Z., *et al.* Preliminary studies on antifungal activity of pyrimorph. *Nongyaoxue Xuebao* **9**(3), 229-234 (2007).
2. Zhao, C., Liu, B., Wang, J., *et al.* Acute toxicity and bioconcentration of pyrimorph in zebrafish, *Brachydanio rerio*. *Pest. Manag. Sci.* **67**(9), 1178-1183 (2011).

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, 07/12/2023

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