

PRODUCT INFORMATION



Picoxystrobin

Item No. 39545

CAS Registry No.: 117428-22-5

Formal Name: α E-(methoxymethylene)-2-[[[6-(trifluoromethyl)-2-pyridinyl]oxy]methyl]-benzeneacetic acid, methyl ester

MF: $C_{18}H_{16}F_3NO_4$

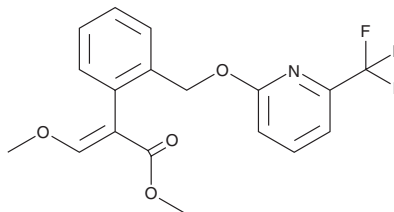
FW: 367.3

Purity: $\geq 98\%$

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

Picoxystrobin is supplied as a solid. A stock solution may be made by dissolving the picoxystrobin in the solvent of choice, which should be purged with an inert gas. Picoxystrobin is slightly soluble (0.1-1 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

Picoxystrobin is a strobilurin fungicide.¹ It is active against the plant pathogenic fungus *R. herpotrichoides* when used at a concentration of 0.75 mg/L.² Picoxystrobin (100 $\mu\text{g/L}$) induces tail deformities, pericardial edema, and spine curvature in zebrafish embryos and is toxic to zebrafish embryos ($\text{LC}_{50} = 213.8 \mu\text{g/L}$).³ Formulations containing picoxystrobin have been used in the control of fungal diseases in agriculture.

References

1. McCreary, C.M., Depuydt, D., Vyn, R.J., et al. Fungicide efficacy of dry bean white mold [*Sclerotinia sclerotiorum* (Lib.) de Bary, causal organism] and economic analysis at moderate to high disease pressure. *Crop Prot.* **82**, 75-81 (2016).
2. Maliński, Z.T. Studies on *in vitro* activity of picoxystrobin and its mixtures with other fungicides against *Ramulispora herpotrichoides* and *Ramulispora acuformis*. *Pestycydy* 75-86 (2008).
3. Jia, W., Mao, L., Zhang, L., et al. Effects of two strobilurins (azoxystrobin and picoxystrobin) on embryonic development and enzyme activities in juveniles and adult fish livers of zebrafish (*Danio rerio*). *Chemosphere* **207**, 573-580 (2018).

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, 07/21/2025

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