

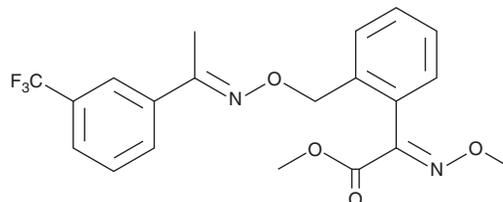
PRODUCT INFORMATION



Trifloxystrobin

Item No. 25825

CAS Registry No.: 141517-21-7
Formal Name: αE-(methoxyimino)-2-[[[(E)-[1-[3-(trifluoromethyl)phenyl]ethylidene]amino]oxy)methyl]-benzeneacetic acid, methyl ester
Synonym: CGA 279202
MF: C₂₀H₁₉F₃N₂O₄
FW: 408.4
Purity: ≥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

Trifloxystrobin is supplied as a solid. A stock solution may be made by dissolving the trifloxystrobin in the solvent of choice, which should be purged with an inert gas. Trifloxystrobin is slightly soluble in chloroform and methanol.

Description

Trifloxystrobin is a fungicide.^{1,2} It inhibits the growth of *R. solani* isolates from sugar beet crops ($EC_{50}s = 0.14-823.54 \mu\text{g/ml}$).¹ Trifloxystrobin (100-1,000 ppm) inhibits mycelial growth of *C. gloeosporioides* when used in combination with tebuconazole (Item No. 24052).² It is toxic to various aquatic species, including *R. arenarum*, *P. santafecinus*, *E. bicolor*, and *L. latrans* tadpoles ($LC_{50}s = 0.22, 0.14, 0.1, \text{ and } 0.26 \text{ mg/L}$, respectively).³ It also reduces mobility of tadpoles and decreases predation of trifloxystrobin-exposed, but not non-exposed, tadpoles by eels. Formulations containing trifloxystrobin have been used in the control of fungi in agriculture and aquatic areas.

References

1. Arabiat, S. and Khan, M. Sensitivity of *Rhizoctonia solani* AG-2-2 from sugar beet to fungicides. *Plant Dis.* **100(12)**, 2427-2433 (2016).
2. Dev, D. and Narendrappa, T. *In vitro* evaluation of fungicides against *Colletotrichum gloeosporioides* (Penz.) Penz and Sacc. causing anthracnose of pomegranate (*Punica granatum* L.). *J. Appl. Nat. Sci.* **8(4)**, 2268-2272 (2017).
3. Junges, C.M., Peltzer, P.M., Lajmanovich, R.C., et al. Toxicity of the fungicide trifloxystrobin on tadpoles and its effect on fish-tadpole interaction. *Chemosphere* **87(11)**, 1348-1354 (2012).

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, 07/26/2023

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