

# PRODUCT INFORMATION



## Trifloxystrobin-d<sub>6</sub>

Item No. 39591

**Formal Name:** methyl (E)-2-(methoxyimino)-2-(2-(((E)-1-(3-(trifluoromethyl)phenyl)ethylidene)amino)oxy)methyl-d<sub>2</sub>)phenyl-3,4,5,6-d<sub>4</sub>)acetate

**MF:** C<sub>20</sub>H<sub>13</sub>D<sub>6</sub>F<sub>3</sub>N<sub>2</sub>O<sub>4</sub>

**FW:** 414.4

**Chemical Purity:** ≥95% (Trifloxystrobin)

**Deuterium**

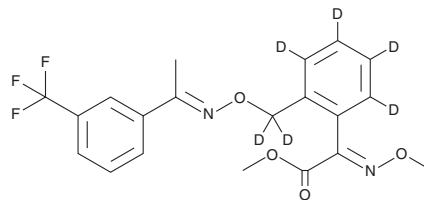
**Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>6</sub>); ≤1% d<sub>0</sub>

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Special Conditions:** Hygroscopic



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

### Laboratory Procedures

Trifloxystrobin-d<sub>6</sub> is intended for use as an internal standard for the quantification of trifloxystrobin (Item No. 25825) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Trifloxystrobin-d<sub>6</sub> is supplied as a solid. A stock solution may be made by dissolving the trifloxystrobin-d<sub>6</sub> in the solvent of choice, which should be purged with an inert gas. Trifloxystrobin-d<sub>6</sub> is slightly soluble in chloroform and methanol.

### Description

Trifloxystrobin is a fungicide.<sup>1,2</sup> It inhibits the growth of *R. solani* isolates from sugar beet crops (EC<sub>50</sub>s = 0.14-823.54 µg/ml).<sup>1</sup> Trifloxystrobin (100-1,000 ppm) inhibits mycelial growth of *C. gloeosporioides* when used in combination with tebuconazole (Item No. 24052).<sup>2</sup> It is toxic to various aquatic species, including *R. arenarum*, *P. santafecinus*, *E. bicolor*, and *L. latrans* tadpoles (LC<sub>50</sub>s = 0.22, 0.14, 0.1, and 0.26 mg/L, respectively).<sup>3</sup> 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, 04/28/2026

#### 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