

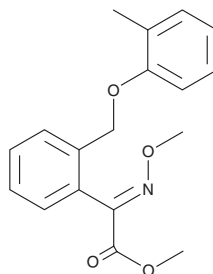
PRODUCT INFORMATION



Kresoxim-methyl

Item No. 25816

CAS Registry No.: 143390-89-0
Formal Name: α E-(methoxyimino)-2-[(2-methylphenoxy)methyl]-benzeneacetic acid, methyl ester
MF: $C_{18}H_{19}NO_4$
FW: 313.4
Purity: $\geq 95\%$
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

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

Description

Kresoxim-methyl is a strobilurin fungicide.¹ It inhibits conidial germination of *V. inaequalis* isolates from apple orchards ($EC_{50}s = 0.00033\text{--}0.0078$ mg/L). Kresoxim-methyl also inhibits mycelial growth ($EC_{50} = 0.240$ mg/L) and is fungicidal against *Saprolegnia* (MIC = 1 mg/L).² It increases intracellular calcium levels and disrupts the mitochondrial membrane potential in mouse cortical cultures in a concentration-dependent manner.³ Kresoxim-methyl is toxic to goldfish (*C. auratus*; $LC_{50} = 0.807$ mg/L).² Formulations containing kresoxim-methyl have been used in the control of fungi in agriculture.

References

1. Fiaccadori, R., Cicognani, E., Alberoni, G., et al. Sensitivity to strobilurin fungicides of Italian *Venturia inaequalis* populations with different origin and scab control. *Pest Manag. Sci.* **67**(5), 535-540 (2011).
2. Hu, X.-G., Liu, L., Hu, K., et al. *In vitro* screening of fungicidal chemicals for antifungal activity against *Saprolegnia*. *J. World Aquac. Soc.* **44**(4), 528-535 (2013).
3. Regueiro, J., Olguín, N., Simal-Gándara, J., et al. Toxicity evaluation of new agricultural fungicides in primary cultured cortical neurons. *Environ. Res.* **140**, 37-44 (2015).

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