

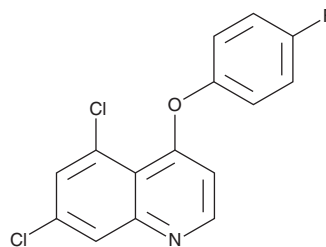
PRODUCT INFORMATION



Quinoxifen

Item No. 42938

CAS Registry No.: 124495-18-7
Formal Name: 5,7-dichloro-4-(4-fluorophenoxy)-quinoline
Synonym: DE-795
MF: $C_{15}H_8Cl_2FNO$
FW: 308.1
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

Quinoxifen is a quinoline fungicide.¹ It reduces the incidence of grape powdery mildew and the mean proportion of infected berries when applied at a concentration of 292 ml/ha to grape vines at the end of bloom stage.² Quinoxifen reduces the feeding rate of *G. fossarum* ($EC_{50} = 23.8 \mu g/L$).¹ It also reduces androgen-mediated transcriptional activity in a reporter assay using MDA-kb2 cells ($IC_{20} = 4.79 \mu M$) and is cytotoxic to MDA-kb2 cells ($EC_{20} = 10.1 \mu M$).³

References

1. Zubrod, J.P., Baudy, P., Schulz, R., *et al.* Effects of current-use fungicides and their mixtures on the feeding and survival of the key shredder *Gammarus fossarum*. **150**, 133-143 (2014).
2. Warneke, B., Thiessen, L.D., and Mahaffee, W.F. Effect of fungicide mobility and application timing on the management of grape powdery mildew. *Plant Dis.* **104(4)**, 1167-1174 (2020).
3. Orton, F., Rosivatz, E., Scholze, M., *et al.* Widely used pesticides with previously unknown endocrine activity revealed as *in vitro* antiandrogens. *Environ. Health Perspect.* **119(6)**, 794-800 (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, 05/07/2025

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