

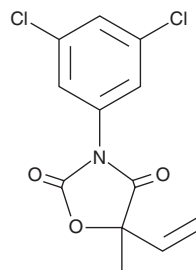
PRODUCT INFORMATION



Vinclozolin

Item No. 23939

CAS Registry No.: 50471-44-8
Formal Name: 3-(3,5-dichlorophenyl)-5-ethenyl-5-methyl-2,4-oxazolidinedione
Synonym: (±)-Vinclozolin
MF: C₁₂H₉Cl₂NO₃
FW: 286.1
Purity: ≥98%
Supplied as: A crystalline 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

Vinclozolin is supplied as a crystalline solid. A stock solution may be made by dissolving the vinclozolin in the solvent of choice, which should be purged with an inert gas. Vinclozolin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of vinclozolin in these solvents is approximately 30 mg/ml.

Vinclozolin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, vinclozolin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Vinclozolin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Vinclozolin is a dicarboximide fungicide.¹ It is active against the plant pathogenic fungi *S. sclerotiorum* and *N. crassa* (EC₅₀s = 0.18 and 1.4 µg/ml, respectively).^{1,2} Vinclozolin decreases the incidence of leaf drop in lettuce fields experimentally infected with *S. sclerotiorum* when applied at a concentration of 1,121 g of active ingredient/hectare (AI/ha).¹

References

1. Matheron, M.E. and Matejka, J.C. In vitro and field comparison of six new fungicides with iprodione and vinclozolin for control of leaf drop of lettuce caused by *Sclerotinia sclerotiorum*. *Plant Dis.* **73**(9), 727-730 (1989).
2. Grindle, M. Isolation and characterization of vinclozolin resistant mutants of *Neurospora crassa*. *Trans. Br. Mycol. Soc.* **82**(4), 635-643 (1984).

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