

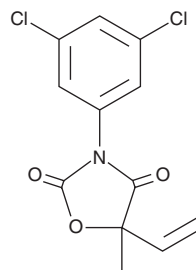
# PRODUCT INFORMATION



## 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:** C<sub>12</sub>H<sub>9</sub>Cl<sub>2</sub>NO<sub>3</sub>  
**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.<sup>1</sup> It is active against the plant pathogenic fungi *S. sclerotiorum* and *N. crassa* (EC<sub>50</sub>s = 0.18 and 1.4 µg/ml, respectively).<sup>1,2</sup> Vinclozolin (0.56 or 1.12 kg/ha) decreases the incidence of leaf drop in lettuce fields experimentally infected with *S. sclerotiorum*.<sup>3</sup>

### 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).
3. Patterson, C.L. and Grogan, R.G. Differences in epidemiology and control of lettuce drop caused by *Sclerotinia minor* and *S. sclerotiorum*. *Plant Dis.* **69(9)**, 766-770 (1985).

#### 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, 07/09/2025

#### 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](http://WWW.CAYMANCHEM.COM)