

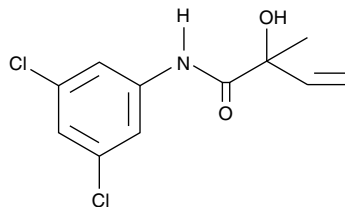
Product Information



Vinclozolin M2

Item No. 10007452

CAS Registry No.: 83792-61-4
Formal Name: N-(3,5-dichlorophenyl)-2-hydroxy-2-methyl-3-butenamide
Synonym: M2
MF: C₁₁H₁₁Cl₂NO₂
FW: 260.1
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 216, 249 nm



Laboratory Procedures

For long term storage, we suggest that vinclozolin M2 (M2) be stored as supplied at -20°C. It should be stable for at least two years.

M2 is supplied as a crystalline solid. A stock solution may be made by dissolving the M2 in an organic solvent purged with an inert gas. M2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of M2 in these solvents is at least 30 mg/ml.

M2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, M2 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. M2 has a solubility of 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.

Vinclozolin is a dicarboximide fungicide widely used in Europe and the United States for control of diseases caused by *B. cinerea*, *S. sclerotiorum*, and *Monilinia* species in grapes, fruits, vegetables, ornamental plants, and turfgrass.¹ Metabolites of vinclozolin, M1 and M2, are effective antagonists of the androgen receptor in rats exhibiting K_i values of 92 and 9.7 μM respectively.² In fathead minnow, M1 and M2 failed to compete for high-affinity, low capacity testosterone binding sites in brain and ovary cytosolic fractions suggesting that they are not anti-androgens in this species.³

References

1. Pothuluri, J.V., Freeman, J.P., Heinze, T.M., *et al.* Biotransformation of vinclozolin by the fungus *Cunninghamella elegans*. *J. Agric. Food Chem.* **48**, 6138-6148 (2000).
2. Kelce, W.R., Monosson, E., Gamcsik, M.P., *et al.* Environmental hormone disruptors: Evidence that vinclozolin developmental toxicity is mediated by antiandrogenic metabolites. *Toxicol. Appl. Pharmacol.* **126**, 276-285 (1994).
3. Makynen, E.A., Kahl, M.D., Jensen, K.M., *et al.* Effects of the mammalian antiandrogen vinclozolin on development and reproduction of the fathead minnow. *Aquatic Toxicology* **48**, 461-475 (2000).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10007452

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 11/29/2012

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com