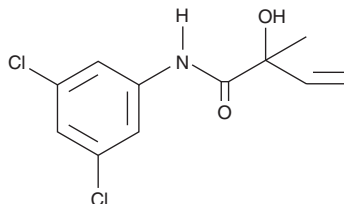


PRODUCT INFORMATION



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%
UV/Vis.: λ_{max}: 216, 249 nm
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 M₂ is supplied as a crystalline solid. A stock solution may be made by dissolving the vinclozolin M₂ in the solvent of choice, which should be purged with an inert gas. Vinclozolin M₂ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of vinclozolin M₂ in these solvents is approximately 30 mg/ml.

Vinclozolin M₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, vinclozolin M₂ should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Vinclozolin M₂ 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 M₂ is an active metabolite of vinclozolin (Item No. 23939).¹ It is formed from vinclozolin by successive esterase activity and decarboxylation of vinclozolin in *C. elegans* and by decarboxylation in human liver microsomes.^{1,2} Vinclozolin M₂ is an antagonist of the mineralocorticoid receptor (IC₅₀ = 1,400 nM) and androgen receptor (IC₅₀ = 0.17 nM) in reporter assays using MCF-7 cells.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**(12), 6138-6148 (2000).
2. Cruz-Hurtado, M., López-González, M.d.L., Mondragón, V., *et al.* *In vitro* phase I metabolism of vinclozolin by human liver microsomes. *Xenobiotica* **49**(8), 895-904 (2019).
3. Molina-Molina, J.-M., Hillenweck, A., Jouanin, I., *et al.* Steroid receptor profiling of vinclozolin and its primary metabolites. *Toxicol. Appl. Pharmacol.* **216**(1), 44-54 (2006).

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, 01/18/2024

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