

# PRODUCT INFORMATION

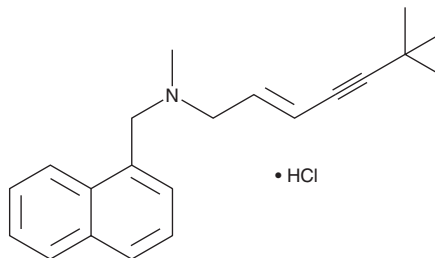


## Terbinafine (hydrochloride)

Item No. 10011619

**CAS Registry No.:** 78628-80-5  
**Formal Name:** N-[(2E)-6,6-dimethyl-2-hepten-4-yn-1-yl]-N-methyl-1-naphthalenemethanamine, monohydrochloride

**MF:** C<sub>21</sub>H<sub>25</sub>N • HCl  
**FW:** 327.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 223, 284 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

Terbinafine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the terbinafine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Terbinafine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of terbinafine (hydrochloride) in these solvents is approximately 30, 12.5, and 14 mg/ml, respectively.

Terbinafine (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, terbinafine (hydrochloride) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Terbinafine (hydrochloride) has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Terbinafine is a broad-spectrum antifungal agent that has activity against *T. rubrum*, *T. mentagrophytes*, *T. verrucosum*, *E. floccosum*, *M. canis*, *A. fumigatus*, and *S. schenckii* (MIC<sub>50</sub>s = 0.003-0.8 µg/ml).<sup>1</sup> It selectively inhibits *C. albicans* squalene epoxidase over rat liver epoxidase (IC<sub>50</sub>s = 0.03 and 77 µM, respectively).<sup>2</sup> Terbinafine (90-120 µM) induces cell cycle arrest at the G<sub>0</sub>/G<sub>1</sub> phase in COLO 205 tumor cells and human umbilical vein endothelial cells (HUVECs).<sup>3,4</sup> Formulations containing terbinafine have been used in the treatment of nail and skin fungal infections.

### References

1. Petranyi, G., Meingassner, J.G., and Mieth, H. Antifungal activity of the allylamine derivative terbinafine *in vitro*. *Antimicrob. Agents Chemother.* **31(9)**, 1365-1368 (1987).
2. Ryder, N.S. and Dupont, M.-C. Inhibition of squalene epoxidase by allylamine antimycotic compounds. A comparative study of the fungal and mammalian enzymes. *Biochem. J.* **230(3)**, 765-770 (1985).
3. Lee, W.-S., Chen, R.-J., Wang, Y.-J., *et al.* *In vitro* and *in vivo* studies of the anticancer action of terbinafine in human cancer cell lines: G<sub>0</sub>/G<sub>1</sub> p53-associated cell cycle arrest. *Int. J. Cancer* **106**, 125-137 (2003).
4. Ho, P.-Y., Liang, Y.-C., Ho, Y.-S., *et al.* Inhibition of human vascular endothelial cells proliferation by terbinafine. *Int. J. Cancer* **111(1)**, 51-59 (2004).

#### 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, 12/19/2022

#### 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