

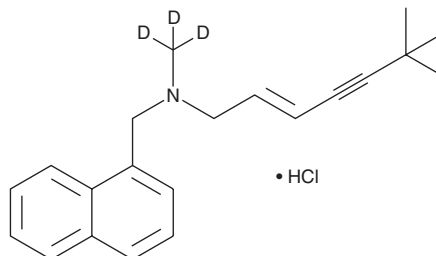
PRODUCT INFORMATION



Terbinafine-d₃ (hydrochloride)

Item No. 26392

CAS Registry No.: 1310012-15-7
Formal Name: N-[-6,6-dimethyl-2-hepten-4-yn-1-yl]-N-(methyl-d₃)-1-naphthalenemethanamine, monohydrochloride
MF: C₂₁H₂₂D₃N • HCl
FW: 330.9
Chemical Purity: ≥98% (Terbinafine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A 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-d₃ (hydrochloride) is intended for use as an internal standard for the quantification of terbinafine (Item No. 10011619) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Terbinafine-d₃ (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the terbinafine-d₃ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Terbinafine-d₃ (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of terbinafine-d₃ (hydrochloride) in these solvents is approximately 30, 10, and 14 mg/ml, respectively.

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* (MIC₅₀s = 0.003-0.8 µg/ml).¹ It selectively inhibits *C. albicans* squalene epoxidase over rat liver epoxidase (IC₅₀s = 0.03 and 77 µM, respectively).² Terbinafine (90-120 µM) induces cell cycle arrest at the G₀/G₁ phase in COLO 205 tumor cells and human umbilical vein endothelial cells (HUVECs).^{3,4} 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: G₀/G₁ p53-associated cell cycle arrest. *Int. J. Cancer* **106(1)**, 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, 10/06/2022

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