

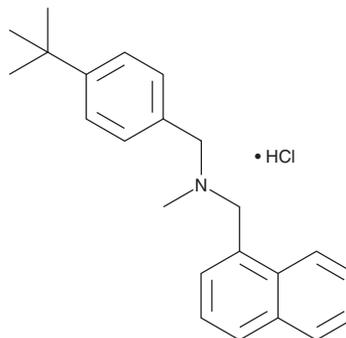
# PRODUCT INFORMATION



## Butenafine (hydrochloride)

Item No. 23797

**CAS Registry No.:** 101827-46-7  
**Formal Name:** N-[[4-(1,1-dimethylethyl)phenyl]methyl]-N-methyl-1-naphthalenemethanamine, monohydrochloride  
**MF:** C<sub>23</sub>H<sub>27</sub>N • HCl  
**FW:** 353.9  
**Purity:** ≥98%  
**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

Butenafine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the butenafine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Butenafine (hydrochloride) is slightly soluble in methanol and DMSO.

### Description

Butenafine is a benzylamine antifungal agent that has broad-spectrum activity against 87 strains of dermatophytes (MICs = 0.0015-0.05 µg/ml), 15 strains of *Aspergillus* (MICs = 0.025-0.78 µg/ml), 4 strains of *Cryptococcus* (MICs = 0.78-1.56 µg/ml), and 67 strains of *Candida* (MICs = 3.13 to >100 µg/ml).<sup>1</sup> It acts by inhibiting squalene epoxidase.<sup>2</sup> Butenafine (50 µM) inhibits the synthesis of ergosterol (Item No. 19850), a sterol in the cell membranes of fungi. Formulations containing butenafine have been used in the treatment of tinea pedis, tinea corporis, and tinea cruris.

### References

1. Maeda, T., Takase, M., Ishibashi, A., *et al.* Synthesis and antifungal activity of butenafine hydrochloride (KP-363), a new benzylamine antifungal agent. *Yakugaku Zasshi* **111**(2), 126-137 (1991).
2. Hiratani, T., Asagi, Y., and Yamaguchi, H. Studies on antifungal mechanism of action of butenafine hydrochloride. II. Comparison in the response to drug treatment between a wild-type strain and tolclate-resistant mutant strains of *Sporothrix schenckii*. *Nihon Ishinkin Gakkai Zasshi* **32**(2), 139-149 (1991).

#### 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