

PRODUCT INFORMATION



Naftifine (hydrochloride)

Item No. 19234

CAS Registry No.: 65473-14-5

Formal Name: N-methyl-N-[(2E)-3-phenyl-2-propen-1-yl]-1-naphthalenemethanamine, monohydrochloride

Synonym: AW 105-843

MF: $C_{21}H_{21}N \cdot HCl$

FW: 323.9

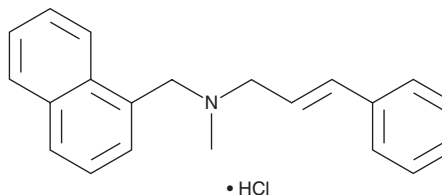
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 222, 255, 282 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Naftifine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the naftifine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Naftifine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of naftifine (hydrochloride) in ethanol and DMSO is approximately 5 mg/ml and approximately 10 mg/ml in DMSO.

Naftifine (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, naftifine (hydrochloride) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Naftifine (hydrochloride) has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Naftifine is an allylamine compound that blocks ergosterol biosynthesis in fungi by inhibiting squalene epoxidase ($K_i = 1.1 \mu M$ for *C. albicans* epoxidase).¹

Reference

1. 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**, 765-770 (1985).

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, 11/16/2022

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