

Terbinafine HCl

Squalene epoxidase inhibitor

Inhibits ergosterol biosynthesis at the level of squalene epoxidase. Potent non-competitive inhibitor of squalene epoxidase from *Candida* ($K_i=30$ nM), inhibiting mammalian enzymes at higher concentrations (4-77 μ M). It inhibits HUVEC proliferation and angiogenesis and displays interesting anti-inflammatory and free radical scavenging activities. Clinically useful antifungal agent.

Citations: 5

[View Online »](#)

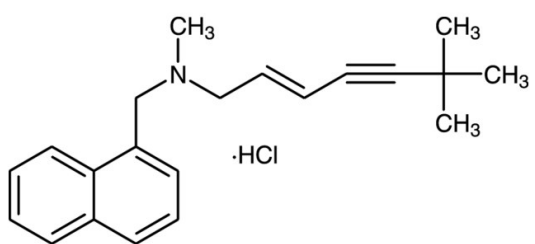
Ordering Information

[Order Online »](#)

BML-EI318-0100	100mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name trans-N-(6,6-Dimethyl-2-hepten-4-ynyl)-N-methyl-1-naphthylmethylamine . hydrochloride

Appearance White to off-white solid.

CAS 78628-80-5

Couple Target Squalene epoxidase

Couple Type Inhibitor

Formula $C_{21}H_{25}N \cdot HCl$

MW 327.9

Purity $\geq 98\%$ (UHPLC)

Solubility Soluble in ethanol (45 mg/ml), DMSO (30 mg/ml) or water (3 mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)