

PRODUCT DATA SHEET



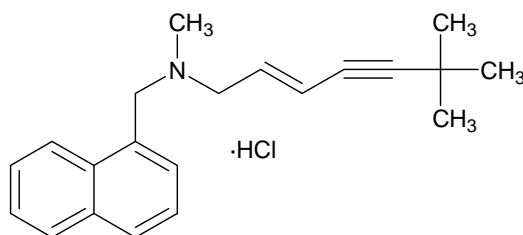
PRODUCT: Terbinafine
trans-N-(6,6-Dimethyl-2-hepten-4-ynyl)-N-methyl-1-naphthylmethylamine-hydrochloride

CAS NO: 78628-80-5

CATALOG NO.: EI-318

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: $C_{21}H_{25}N \cdot HCl$

MOLECULAR WEIGHT: 327.9

PURITY: 98% (TLC)

SOLUBILITY: Soluble in ethanol (45 mg/ml), DMSO (30 mg/ml), Water (3 mg/ml)

STORAGE: Store, as supplied, at room temperature for up to 1 year. Store solutions at $-20^{\circ}C$ for up to 3 months.

APPLICATION NOTES: Inhibits ergosterol biosynthesis at the level of squalene epoxidase.¹ Potent non-competitive inhibitor of squalene epoxidase from *Candida* ($K_i = 30$ nM), inhibiting mammalian enzymes at higher concentrations (4-77 μM).^{1,2} It inhibits HUVEC proliferation and angiogenesis³ and displays interesting anti-inflammatory⁴ and free radical scavenging⁵ activities. Clinically useful antifungal agent.

- REFERENCES:
1. N.S. Ryder *Br. J. Dermatol.* 1992 **126** supp 39,2
 2. B. Favre *et al. Arch. Biochem. Biophys.* 1997 **340** 265
 3. P.Y. Ho *et al. Int. J. Cancer* 2004 **111** 51
 4. T. Rosen *et al. Int. J. Dermatol.* 1997 **36** 788
 5. C.S. Sander *et al. Mycoses* 2002 **45** 152

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.