

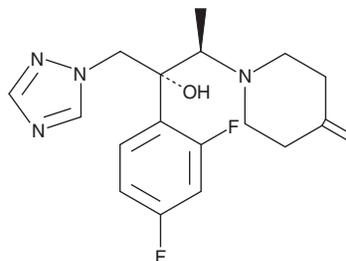
# PRODUCT INFORMATION



## Efinaconazole

Item No. 23839

**CAS Registry No.:** 164650-44-6  
**Formal Name:**  $\alpha$ R-(2,4-difluorophenyl)- $\beta$ R-methyl-4-methylene- $\alpha$ -(1H-1,2,4-triazol-1-ylmethyl)-1-piperidineethanol  
**Synonym:** KP-103  
**MF:** C<sub>18</sub>H<sub>22</sub>F<sub>2</sub>N<sub>4</sub>O  
**FW:** 348.4  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 262 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

Efinaconazole is supplied as a crystalline solid. A stock solution may be made by dissolving the efinaconazole in the solvent of choice, which should be purged with an inert gas. Efinaconazole is soluble in the organic solvent DMSO. The solubility of efinaconazole in DMSO is >14 mg/ml. It is also slightly soluble in methanol and chloroform.

### Description

Efinaconazole is a broad-spectrum triazole antifungal agent with activity against *Acremonium*, *Aspergillus*, *Candida*, *Cryptococcus*, *Epidermophyton*, *Fusarium*, *Microsporum*, *Paecilomyces*, *Pseudallescheria*, *Scopulariopsis*, *Trichophyton*, and *Trichosporon*.<sup>1</sup> It inhibits the growth of *T. rubrum* and *T. mentagrophytes* clinical isolates with MIC values ranging from  $\leq$ 2.0 to 60 ng/ml and of *C. albicans* isolates with MIC values ranging from  $\leq$ 0.5 to >250 ng/ml. Efinaconazole inhibits sterol 14 $\alpha$ -demethylase, which arrests ergosterol (Item No. 19850) biosynthesis at the fungal membrane.<sup>2</sup> It inhibits ergosterol biosynthesis in *T. mentagrophytes* and *C. albicans* with IC<sub>50</sub> values of 7.0 and 0.40 ng/ml, respectively. Topical formulations containing efinaconazole have been used for the treatment of onychomycosis.

### References

1. Jo Siu, W.J., Tatsumi, Y., Senda, H., *et al.* Comparison of *in vitro* antifungal activities of efinaconazole and currently available antifungal agents against a variety of pathogenic fungi associated with onychomycosis. *Antimicrob. Agents Chemother.* **57(4)**, 1610-1616 (2013).
2. Tatsumi, Y., Nagashima, M., Shibanushi, T., *et al.* Mechanism of action of efinaconazole, a novel triazole antifungal agent. *Antimicrob. Agents Chemother.* **57(5)**, 2405-2409 (2013).

#### 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, 11/14/2022

#### 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