

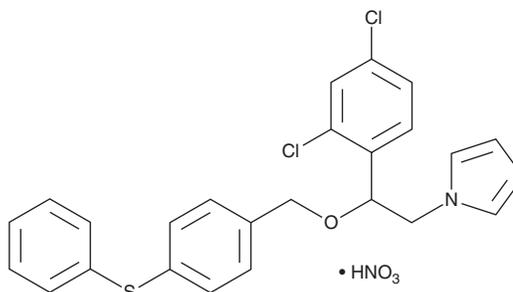
# PRODUCT INFORMATION



## Fenticonazole (nitrate)

Item No. 29002

**CAS Registry No.:** 73151-29-8  
**Formal Name:** 1-[2-(2,4-dichlorophenyl)-2-[[4-(phenylthio)phenyl]methoxy]ethyl]-1H-imidazole, mononitrate  
**Synonym:** REC 15-1476  
**MF:** C<sub>24</sub>H<sub>20</sub>Cl<sub>2</sub>N<sub>2</sub>OS • HNO<sub>3</sub>  
**FW:** 518.4  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Fenticonazole (nitrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the fenticonazole (nitrate) in the solvent of choice, which should be purged with an inert gas. Fenticonazole (nitrate) is soluble in chloroform and methanol.

### Description

Fenticonazole is an imidazole with antimicrobial activity.<sup>1,2</sup> It is active against clinical isolates of *Candida* species, including *C. albicans*, *C. parapsilosis*, *C. tropicalis*, *C. krusei*, and *C. guilliermondii* with MIC values ranging from 0.12 to 32 µg/ml.<sup>1</sup> It is also active against clinical isolates of bacteria, including *B. fragilis*, associated with bacterial vaginosis (MICs = ≤0.03-16 µg/ml), as well as bacteria, including *S. aureus* and *S. epidermidis*, associated with skin infections with MIC values ranging from 0.03 µg/ml or less to greater than 16 µg/ml.<sup>2</sup> Fenticonazole inhibits fungal growth in a guinea pig model of experimental *M. canis* dermatomycosis (ED<sub>50</sub> = 0.5%, applied topically).<sup>3</sup>

### References

1. Hernández Molina, J.M., Llosá, J., Martínez Brocal, A., et al. *In vitro* activity of cloconazole, sulconazole, butoconazole, isoconazole, fenticonazole, and five other antifungal agents against clinical isolates of *Candida albicans* and *Candida* spp. *Mycopathologia* **118**(1), 15-21 (1992).
2. Jones, B.M., Geary, I., Lee, M.E., et al. Comparison of the *in vitro* activities of fenticonazole, other imidazoles, metronidazole, and tetracycline against organisms associated with bacterial vaginosis and skin infections. *Antimicrob. Agents Chemother.* **33**(6), 970-972 (1989).
3. Veronese, M., Barzaghi, D., Bertoncini, A., et al. Fenticonazole: A new antifungal imidazole derivative *in vitro* and *in vivo* antimycotic activity. *Mykosen* **27**(4), 194-202 (1984).

#### 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, 12/04/2019

#### 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