

# PRODUCT INFORMATION



## PF-1163B

Item No. 22066

**CAS Registry No.:** 258871-60-2  
**Formal Name:** 3S-[[4-(2-hydroxyethoxy)phenyl]methyl]-4,10-dimethyl-13R-pentyl-1-oxa-4-azacyclotridecane-2,5-dione

**MF:** C<sub>27</sub>H<sub>43</sub>NO<sub>5</sub>

**FW:** 461.6

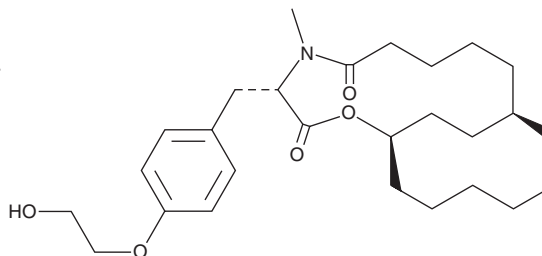
**Purity:** ≥95%

**Supplied as:** An oil

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Fungus/Unidentified sp.



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

### Laboratory Procedures

PF-1163B is supplied as an oil. A stock solution may be made by dissolving the PF-1163B in the solvent of choice, which should be purged with an inert gas. PF-1163B is soluble in acetonitrile.

### Description

PF-1163B is a depsipeptide antifungal isolated from *Penicillium* that inhibits ergosterol (ERG) synthesis (IC<sub>50</sub> = 34 ng/ml).<sup>1,2</sup> It inhibits growth of *C. albicans* (MIC = 32 µg/ml) but not of other *Candida* strains, *A. fumigatus*, or HepG2 cells.<sup>2</sup> It acts synergistically with fluconazole (Item No. 11594) to reduce growth of azole-resistant *C. albicans* (MIC = 2 and 0.0016 µg/ml alone and in combination, respectively).<sup>3</sup> PF-1163B is a hydroxy form of PF-1163A (Item No. 22065).

### References

1. Nose, H., Seki, A., Yaquchi, T., *et al.* PF1163A and B, new antifungal antibiotics produced by *Penicillium* sp. I. Taxonomy of producing strain, fermentation, isolation and biological activities. *J. Antibiot. (Tokyo)* **53(1)**, 33-37 (2000).
2. Sasaki, T., Nose, H., Hosoya, A., *et al.* PF1163A and B, new antifungal antibiotics produced by *Penicillium* sp. II. Physico-chemical properties and structure elucidation. *J. Antibiot. (Tokyo)* **53(1)**, 38-44 (2000).
3. Okabe, M., Sugita, T., Kinoshita, K., *et al.* Macrolides from a marine-derived fungus, *Penicillium meleagrinum* var. *viridiflavum*, showing synergistic effects with Fluconazole against azole-resistant *Candida albicans*. *J.Nat. Prod.* **79(4)**, 1208-1212 (2016).

#### 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, 12/20/2023

#### 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