

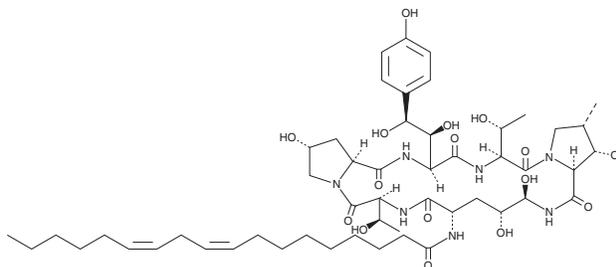
# PRODUCT INFORMATION



## Echinocandin B

Item No. 25520

CAS Registry No.: 54651-05-7  
Synonyms: Antibiotic A-22082,  
Antibiotic A-30912A,  
Antibiotic SL-7810F  
MF:  $C_{52}H_{81}N_7O_{16}$   
FW: 1,060.2  
Purity:  $\geq 85\%$   
Supplied as: A powder  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years  
Item Origin: Fungus/*Aspergillus*  
sp.



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

### Laboratory Procedures

Echinocandin B is supplied as a powder. A stock solution may be made by dissolving the echinocandin B in the solvent of choice, which should be purged with an inert gas. Echinocandin B is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Echinocandin B is an echinocandin antifungal originally isolated from *Aspergillus*.<sup>1</sup> Echinocandin B is an inhibitor of 1,3- $\beta$  glucan synthase, and it inhibits incorporation of radiolabeled glucose into alkali-insoluble glucan by *S. cerevisiae* spheroplasts in a dose-dependent manner.<sup>2,3</sup> It inhibits growth of *C. albicans* *in vitro* with an MIC value of 3.9  $\mu g/ml$  after 2 days.<sup>4</sup>

### References

1. Hector, R.F. Compounds active against cell walls of medically important fungi. *Clin. Microbiol. Rev.* **6(1)**, 1-21 (1993).
2. Baguley, C., Römmele, G., Gruner, J., et al. Papulacandin B: An inhibitor of glucan synthesis in yeast spheroplasts. *Eur. J. Biochem.* **97(2)**, 345-351 (1979).
3. Sawistowska-Schröder, E.T., Kerridge, D., and Perry, H. Echinocandin inhibition of 1,3- $\beta$ -D-glucan synthase from *Candida albicans*. *FEBS Lett.* **173(1)**, 134-138 (1984).
4. Frost, D.J., Brandt, K.D., Cugier, D., et al. A whole-cell *Candida albicans* assay for the detection of inhibitors towards fungal cell wall synthesis and assembly. *J. Antibiot. (Tokyo)* **48(4)**, 306-310 (1995).

#### 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, 08/10/2023

#### 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