

PRODUCT INFORMATION



Caspofungin-d₄ (acetate)

Item No. 29096

Formal Name: 1-[(4R,5S)-5-[(2-aminoethyl-1,1,2,2-d₄) amino]-N²-(10,12-dimethyl-1-oxotetradecyl)-4-hydroxy-L-ornithine]-5-[(3R)-3-hydroxy-L-ornithine]-pneumocandin B₀, diacetate

MF: C₅₂H₈₄D₄N₁₀O₁₅ • 2C₂H₄O₂

FW: 1,217.5

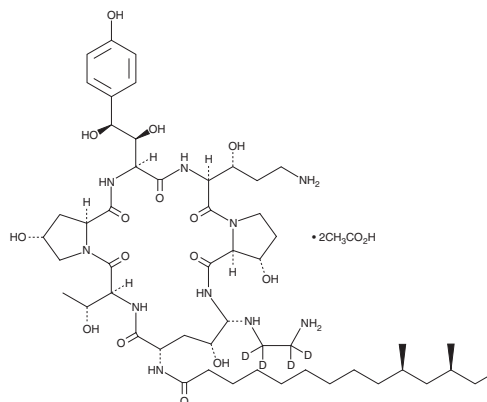
Chemical Purity: ≥95% (Caspofungin)

Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Caspofungin-d₄ (acetate) is intended for use as an internal standard for the quantification of caspofungin (Item No. 15923) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Caspofungin-d₄ (acetate) is supplied as a solid. A stock solution may be made by dissolving the caspofungin-d₄ (acetate) in the solvent of choice, which should be purged with an inert gas. Caspofungin-d₄ (acetate) is slightly soluble in the organic solvent methanol.

Description

Caspofungin is an antifungal compound that is classified as an echinocandin, as it acts by noncompetitively inhibiting the enzyme 1,3-β glucan synthase and preventing the synthesis of glucan in the fungal cell wall.¹ It is effective against disseminated or invasive aspergillosis and candidiasis.^{2,3} Caspofungin has a 50% minimum effective concentration against *Aspergillus* spp. of 0.015 μg/ml.⁴ It does not antagonize other antifungal compounds, like amphotericin B, and, in some cases, may show additive or synergistic activity.^{1,3,5}

References

1. Letscher-Bru, V. and Herbrecht, R. Caspofungin: The first representative of a new antifungal class. *J. Antimicrob. Chemother.* **51(3)**, 513-521 (2003).
2. Abruzzo, G.K., Flattery, A.M., Gill, C.J., et al. Evaluation of the echinocandin antifungal MK-0991 (L-743,872): Efficacies in mouse models of disseminated aspergillosis, candidiasis, and cryptococcosis. *Antimicrob. Agents Chemother.* **41(11)**, 2333-2338 (1997).
3. Deresinski, S.C. and Stevens, D.A. Caspofungin. *Clin. Infect. Dis.* **36(11)**, 1445-1457 (2003).
4. Pfaller, M.A., Boyken, L., Hollis, R.J., et al. In vitro susceptibility of clinical isolates of *Aspergillus* spp. to anidulafungin, caspofungin, and micafungin: A head-to-head comparison using the CLSI M38-A2 broth microdilution method. *J. Clin. Microbiol.* **47(10)**, 3323-3325 (2009).
5. Herbrecht, R., Nivoix, Y., Fohrer, C., et al. Management of systemic fungal infections: Alternatives to itraconazole. *J. Antimicrob. Chemother.* **56(Suppl 1)**, i39-i48 (2005).

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, 10/24/2022

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