

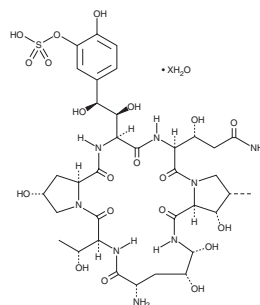
# PRODUCT INFORMATION



## FR179642 (hydrate)

Item No. 29647

**CAS Registry No.:** 1392483-15-6  
**Formal Name:** 1-[(4R,5R)-4,5-dihydroxy-L-ornithine]-4-[(4S)-4-hydroxy-4-[4-hydroxy-3-(sulfooxy)phenyl]-L-threonine]-pneumocandin A<sub>0</sub>, hydrate  
**MF:** C<sub>35</sub>H<sub>52</sub>N<sub>8</sub>O<sub>20</sub>S • XH<sub>2</sub>O  
**FW:** 936.9  
**Purity:** ≥95%  
**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

FR179642 (hydrate) is supplied as a solid. A stock solution may be made by dissolving the FR179642 (hydrate) in the solvent of choice, which should be purged with an inert gas. FR179642 (hydrate) is slightly soluble in DMSO and methanol (warmed).

### Description

FR179642 is the cyclic peptide nucleus of the lipopeptide antifungal FR901379 and an intermediate in the synthesis of the echinocandin antifungal FK463 (micafungin; Item No. 18009).<sup>1,2</sup>

### References

1. Tomishima, M., Ohki, H., Yamada, A., *et al.* FK463, a novel water-soluble echinocandin lipopeptide: Synthesis and antifungal activity. *J. Antibiot. (Tokyo)* **52(7)**, 674-676 (1999).
2. Ueda, S., Shibata, T., Ito, K., *et al.* Cloning and expression of the FR901379 acylase gene from *Streptomyces* sp. no. 6907. *J. Antibiot. (Tokyo)* **64(2)**, 169-175 (2011).

#### 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, 11/17/2022

#### 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