

# PRODUCT INFORMATION

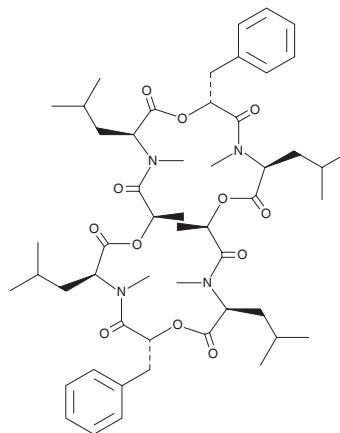


## PF-1022A

Item No. 29700

**CAS Registry No.:** 133413-70-4  
**Formal Name:** cyclo[(αR)-α-hydroxybenzenepropanoyl-N-methyl-L-leucyl-(2R)-2-hydroxypropanoyl-N-methyl-L-leucyl-(αR)-α-hydroxybenzenepropanoyl-N-methyl-L-leucyl-(2R)-2-hydroxypropanoyl-N-methyl-L-leucyl]

**MF:** C<sub>52</sub>H<sub>76</sub>N<sub>4</sub>O<sub>12</sub>  
**FW:** 949.2  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



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

### Laboratory Procedures

PF-1022A is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-1022A in the solvent of choice, which should be purged with an inert gas. PF-1022A is soluble in the organic solvent DMSO at a concentration of approximately 43 mg/ml.

### Description

PF-1022A is a cyclodepsipeptide fungal metabolite originally isolated from mycelia sterilia and has anthelmintic activity.<sup>1</sup> Oral administration of PF-1022A (0.5-2 mg/kg) induces worm excretion in a chicken model of *A. galli* infection. It induces parasite clearance in jird models of *H. contortus*, *O. ostertagi*, or *T. colubriformis* infection (ED<sub>95</sub>s = 0.11, 0.33, and 0.11 mg/animal).<sup>2</sup>

### References

1. Sasaki, T., Takagi, M., Yaguchi, T., *et al.* A new anthelmintic cyclodepsipeptide, PF1022A. *J. Antibiot. (Tokyo)* **45(5)**, 692-697 (1992).
2. Conder, G.A., Johnson, S.S., Nowakowski, D.S., *et al.* Anthelmintic profile of the cyclodepsipeptide PF1022A in *in vitro* and *in vivo* models. *J. Antibiot. (Tokyo)* **48(8)**, 820-823 (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, 11/03/2022

#### 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