

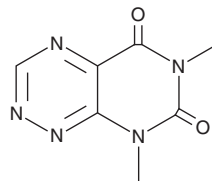
PRODUCT INFORMATION



Fervenulin

Item No. 33499

CAS Registry No.: 483-57-8
Formal Name: 6,8-dimethyl-pyrimido[5,4-e]-1,2,4-triazine-5,7(6H,8H)-dione
Synonyms: NSC 68158, Planomycin, U 7118
MF: C₇H₇N₅O₂
FW: 193.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

Fervenulin is supplied as a solid. A stock solution may be made by dissolving the fervenulin in the solvent of choice, which should be purged with an inert gas. Fervenulin is soluble in methanol and DMSO.

Description

Fervenulin is a bacterial metabolite originally isolated from *Streptomyces* that has nematocidal activity.^{1,2} It inhibits egg hatching of the root-knot nematode *M. incognita* (MIC = 30 µg/ml), as well as induces mortality in *M. incognita* second-stage juveniles when used at a concentration of 250 µg/ml.²

References

1. DeBoer, C., Dietz, A., Evans, J.S., *et al.* Fervenulin, a new crystalline antibiotic. I. Discovery and biological activities. *Antibiot Annu.* **7**, 220-226 (1960).
2. Ruanpanun, P., Laatsch, H., Tangchitsomkid, N., *et al.* Nematicidal activity of fervenulin isolated from a nematicidal actinomycete, *Streptomyces* sp. CMU-MH021, on *Meloidogyne incognita*. *World J. Microbiol. Biotechnol.* **27(6)**, 1373-1380 (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, 11/15/2022

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