

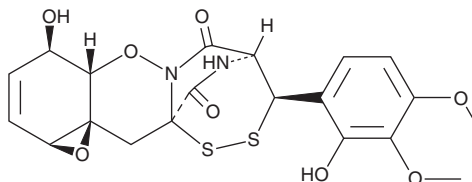
PRODUCT INFORMATION



Gliovirin

Item No. 28583

CAS Registry No.: 83912-90-7
Formal Name: (1aS,4R,4aS,8S,9R,11aR,12aS)-4,4a,8,9-tetrahydro-4-hydroxy-9-(2-hydroxy-3,4-dimethoxyphenyl)-12H-8,11a-(iminomethano)-1aH,7H-[1,2,4]dithiazepino[4,3-b]oxireno[e][1,2]benzoxazine-7,13-dione



MF: C₂₀H₂₀N₂O₈S₂
FW: 480.5
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Penicillium* sp.

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

Laboratory Procedures

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

Description

Gliovirin is a fungal metabolite that has been found in *T. harzianum* and has fungicidal, antimicrobial and anti-inflammatory activities.¹ It is active against the plant pathogenic fungus *P. ultimum* (MIC = 60 ng/ml) and the parasite *T. brucei brucei* (IC₅₀ = 90 ng/ml), but has no effect on the plant pathogenic fungi *R. solani*, *P. omnivorum*, *T. basicola*, *R. arrhizus*, and *V. dahliae* or the bacteria *B. thuringiensis*, *P. fluorescens*, and *X. malvacearum* when used at concentrations up to 1,000 ng/ml.^{2,3} Gliovirin decreases phorbol 12-myristate 13-acetate (TPA)- and ionomycin-induced increased expression of COX-2 (IC₅₀ = 1 μM) and protein levels of IL-2 in Jurkat cells (IC₅₀ = 5.2 μM).¹

References

1. Rether, J., Serwe, A., Anke, T., *et al.* Inhibition of inducible tumor necrosis factor-α expression by the fungal epipolythiodiketopiperazine gliovirin. *Biol. Chem.* **388**(6), 627-637 (2007).
2. Howell, C.R. and Stipanovic, R.D. Gliovirin, a new antibiotic from *Gliocladium virens*, and its role in the biological control of *Pythium ultimum*. *Can. J. Microbiol.* **29**(3), 321-324 (1983).
3. Iwatsuki, M., Otoguro, K., Ishiyama, A., *et al.* *In vitro* antitrypanosomal activity of 12 low-molecular-weight antibiotics and observations of structure/activity relationships. *J. Antibiot. (Tokyo)* **63**(10), 619-622 (2010).

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, 08/03/2023

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