

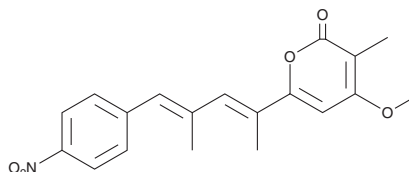
PRODUCT INFORMATION



Luteoreticulin

Item No. 25515

CAS Registry No.: 22388-89-2
Formal Name: 6-[(1E,3E)-1,3-dimethyl-4-(4-nitrophenyl)-1,3-butadien-1-yl]-4-methoxy-3-methyl-2H-pyran-2-one
Synonym: Griseulin
MF: C₁₉H₁₉NO₅
FW: 341.4
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 369 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces luteoreticuli*



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

Laboratory Procedures

Luteoreticulin is supplied as a solid. A stock solution may be made by dissolving the luteoreticulin in the solvent of choice, which should be purged with an inert gas. Luteoreticulin is soluble in organic solvents such as DMSO and dimethyl formamide.

Description

Luteoreticulin is a nitro-containing bacterial metabolite originally isolated from *S. luteoreticuli*.¹ It has mosquitocidal activity against *A. aegypti* when used at a concentration of 6.25 mg/L and nematocidal activity against *C. elegans*.²

References

1. Koyama, Y., Fukakusa, Y., Kyomura, N., *et al.* The structure of luteoreticulin, a nitro-containing metabolite of *Streptomyces luteoreticuli*. *Tetrahedron Lett.* **10(5)**, 355-358 (1969).
2. Nair, M.G., Chandra, A., Thorogod, D.L., *et al.* Nematicidal and mosquitocidal aromatic nitro compounds produced by *Streptomyces* spp. *Pesticide Science* **43(4)**, 361-365 (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, 12/07/2022

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