

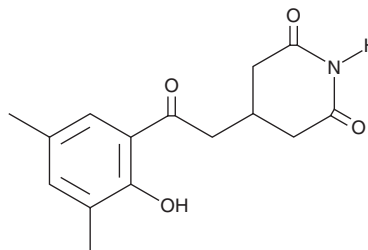
PRODUCT INFORMATION



Actiphenol

Item No. 33119

CAS Registry No.: 526-02-3
Formal Name: 4-[2-(2-hydroxy-3,5-dimethylphenyl)-2-oxoethyl]-2,6-piperidinedione
Synonyms: Actinophenol, NSC 58413
MF: C₁₅H₁₇NO₄
FW: 275.3
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

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

Description

Actiphenol is a bacterial metabolite that has been found in *Streptomyces* and has antiviral activity.^{1,2} It is active against coxsackievirus B3 and influenza A virus (IC₅₀s = 14.37 and 34.4 µg/ml, respectively).¹ Actiphenol (50 µg/disc) potentiates the antifungal activity of miconazole (Item No. 15420) against *C. albicans* in a disc assay.²

References

1. Ji, X.-Y., Zhong, Z.-J., Xue, S.-T., *et al.* Synthesis and antiviral activities of synthetic glutarimide derivatives. *Chem. Pharm. Bull. (Tokyo)* **58(11)**, 1436-1441 (2010).
2. Fukuda, T., Matsumoto, A., Takahashi, Y., *et al.* Phenatic acids A and B, new potentiators of antifungal miconazole activity produced by *Streptomyces* sp. K03-0132. *J. Antibiot. (Tokyo)* **58(4)**, 252-259 (2005).

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, 01/03/2023

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