

PRODUCT INFORMATION



Sultricin

Item No. 27976

CAS Registry No.: 131774-59-9

Formal Name: (5S,6S)-5,6-dihydro-5-hydroxy-6-[(1E,4S,5S,6S,7Z,9Z,11E)-6-hydroxy-5-methyl-4-(phosphonooxy)-1,7,9,11-heptadecatetraen-1-yl]-2H-pyran-2-one, monosodium salt

Synonyms: BU-3285T, Phostricin

MF: $C_{23}H_{34}O_8P \cdot Na$

FW: 492.5

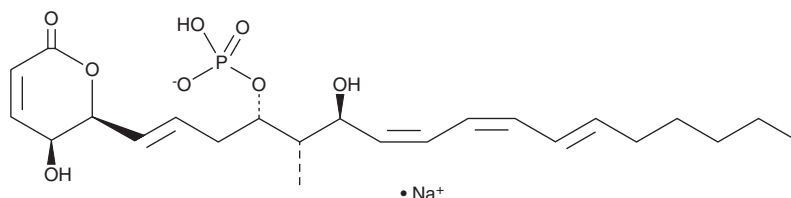
Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

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

Sultricin is supplied as a solid. A stock solution may be made by dissolving the sultricin in water. We do not recommend storing the aqueous solution for more than one day.

Description

Sultricin is a fungal metabolite originally isolated from *S. roseiscleroticus*.¹ It is active against a panel of 15 fungal strains (MICs = 0.8-50 $\mu g/ml$). Sultricin is cytotoxic to B16/F10, HCT116, Moser, P388, A549, and K562 cancer cells (IC_{50} s = 37.7, 1.92, 0.85, 3.8, 4.14, and 0.67 $\mu g/ml$, respectively). It also inhibits RNA and protein synthesis in L1210 leukemia cells (IC_{50} s = 5.5 and 2.4 $\mu g/ml$, respectively). Sultricin (0.1-10 mg/kg) increases survival time in a B16/F10 syngeneic mouse melanoma model and a P388 murine leukemia model.

Reference

1. Ohkuma, H., Naruse, N., Nishiyama, Y., *et al.* Sultricin, a new antifungal and antitumor antibiotic from *Streptomyces roseiscleroticus*. Production, isolation, structure and biological activity. *J. Antibiot. (Tokyo)* 45(8), 1239-1249 (1992).

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, 12/22/2022

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