

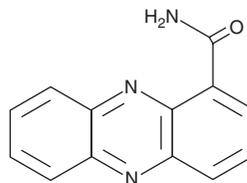
PRODUCT INFORMATION



Oxychlororaphine

Item No. 27624

CAS Registry No.: 550-89-0
Formal Name: 1-phenazinecarboxamide
MF: $C_{13}H_9N_3O$
FW: 223.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 249, 367 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Synthetic



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

Description

Oxychlororaphine is a phenazine microbial metabolite that has been found in a variety of bacteria and has anticancer, antibacterial, and antifungal activities.^{1,2} It is cytotoxic to A549, HeLa, and SW480 cancer cells (IC_{50} s = 32-40 μM).¹ Oxychlororaphine halts the cell cycle at the G_1 phase and induces apoptosis in A549 cells. It is active against *Streptomyces* bacteria and *T. mentagrophytes*, *E. floccosum*, *C. albicans*, *M. audouini*, and *B. dermatitidis* fungi when used at concentrations ranging from 12.5 to 50 ppm.²

References

1. Ali, H.M., El-Shikh, M.S., Salem, M.Z.M., *et al.* Isolation of bioactive phenazine-1-carboxamide from the soil bacterium *Pantoea agglomerans* and study of its anticancer potency on different cancer cell lines. *J. AOAC Int.* **99**(5), 1233-1239 (2016).
2. Sierra, G. and Veringa, H.A. Effect of oxy-chlororaphin on the growth *in vitro* of *Streptomyces* species and some pathogenic fungi. *Nature* **182**(4630), 265 (1958).

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, 12/19/2022

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