

PRODUCT INFORMATION

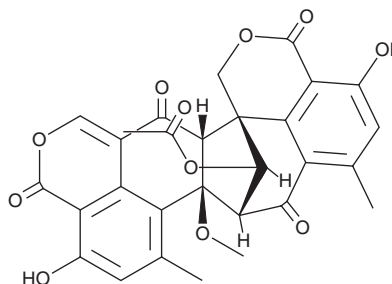


Duclauxin

Item No. 22068

CAS Registry No.: 1732-37-2
Formal Name: (8S,8aS,15aR,15bR,16S)-16-(acetyloxy)-8a,15a-dihydro-4,11-7H-8,15b-methano-1H,3H,12H-benzo[de]cyclohepta[1,2-g:3,4,5-d'e']bis[2]benzopyran-3,7,12,15(8H)-tetrone

MF: C₂₉H₂₂O₁₁
FW: 546.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Description

Duclauxin decreases proliferation of tumor cells *in vitro* and increases the lifespan of mice inoculated with Ehrlich ascitic tumor cells.^{1,2} Duclauxin (10-30 µg/ml) inhibits mitochondrial respiration of P-388 tumor cells and nucleic acid synthesis in cell culture of Ehrlich cancer, the lymphadenomas NK/LI and L 5178, and sarcoma 37.^{3,4} Duclauxin also inhibits growth of wheat coleoptile.^{4,5}

References

1. Fusková, A., Proksa, B., and Fuska, J. *In vitro* effect of duclauxin and derivatives of coumarin on nucleic acid and protein synthesis in Ehrlich's Ascites Carcinoma cells (EAC). *Pharmazie* **32(5)**, 291-293 (1977).
2. Kawai, K., Nozawa, Y., Ito, T., *et al.* Effects of xanthomegnin and duclauxin on culture cells of murine leukemia and Ehrlich ascitic tumor. *Res. Commun. Chem. Pathol. Pharmacol.* **36(3)**, 429-438 (1982).
3. Kovác, L., Böhmerová, E., and Fuska, J. Inhibition of mitochondrial functions by the antibiotics, bikaverin and duclauxine. *J. Antibiot. (Tokyo)* **31(6)**, 616-620 (1978).
4. Fuska, L.A., Ivanitskaia, L.P., Makukho, L.V., *et al.* The effects of the antibiotics vermiculin PSX-1, bicaverin and duclauxin, isolated from fungi, on nucleic acid synthesis in several tumors *Antibiotiki* **19(10)**, 33-37 (1974).
5. Bryant, F.O., Cutler, H.F., and Jacyno, J.M. Properties and cost effective method for production of the antitumor agent decauxin from sporulating *Penicillium herquei*. *J. Pharm. Sci.* **82(12)**, 1214-1217 (1993).

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, 11/11/2022

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