

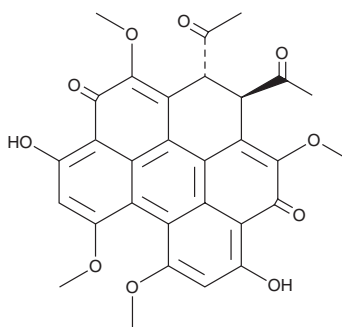
PRODUCT INFORMATION



Elsinochrome A

Item No. 33072

CAS Registry No.: 24568-67-0
Formal Name: (1R,2R)-*rel*-1,2-diacetyl-1,2-dihydro-5,10-dihydroxy-3,7,8,12-tetramethoxybenzo[ghi]perylene-4,11-dione
Synonym: NSC 623609
MF: C₃₀H₂₄O₁₀
FW: 544.5
Purity: ≥95%
UV/Vis.: λ_{max}: 261, 457 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungi/*Aspergillus* sp.



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

Laboratory Procedures

Elsinochrome A is supplied as a solid. A stock solution may be made by dissolving the elsinochrome A in the solvent of choice, which should be purged with an inert gas. Elsinochrome A is soluble in ethanol.

Description

Elsinochrome A is a perylenequinone fungal metabolite that has been found in *Hypomyces* and has anticancer activity.¹ Photoactivation of elsinochrome A induces the production of singlet oxygen and apoptosis in R366.4 rhesus monkey embryonic stem cells and HCe-8693 colorectal carcinoma cells when used at concentrations ranging from 1 to 100 μM.

Reference

1. Ma, L., Tai, H., Li, C., *et al.* Photodynamic inhibitory effects of three perylenequinones on human colorectal carcinoma cell line and primate embryonic stem cell line. *World J. Gastroenterol.* **9(3)**, 485-490 (2003).

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, 12/16/2022

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