

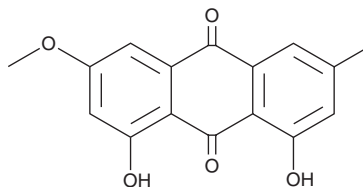
PRODUCT INFORMATION



Physcion

Item No. 19863

CAS Registry No.: 521-61-9
Formal Name: 1,8-dihydroxy-3-methoxy-6-methyl-9,10-anthracenedione
Synonyms: 6-O-methyl Emodin, NSC 251670
MF: C₁₆H₁₂O₅
FW: 284.3
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 266, 286, 434 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

Physcion is a natural anthraquinone. It has been reported to have laxative and anti-inflammatory actions, inhibit kinases and tyrosinase, and have cytotoxicity against cancer cells.¹ Physcion inhibits 6-phosphogluconate dehydrogenase (IC₅₀ = 38.5 μM), decreasing lipogenesis and RNA biosynthesis in cancer cells, leading to reduced cell proliferation and tumor growth in mice.²

References

1. Chien, S.-C., Wu, Y.-C., and Chen, Z.-W. Naturally occurring anthraquinones: Chemistry and therapeutic potential in autoimmune diabetes. *Evid. Based Complement. Alternat. Med.* **2015:35757** (2015).
2. Lin, R., Elf, S., Shan, C., et al. 6-Phosphogluconate dehydrogenase links oxidative PPP, lipogenesis and tumour growth by inhibiting LKB1-AMPK signalling. *Nat. Cell Biol.* **17(11)**, 1484-1496 (2015).

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, 11/29/2022

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