

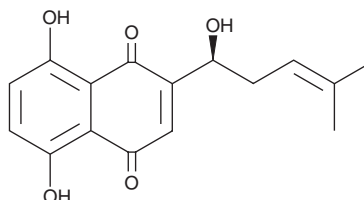
PRODUCT INFORMATION



Alkannin

Item No. 25058

CAS Registry No.: 517-88-4
Formal Name: 5,8-dihydroxy-2-[(1S)-1-hydroxy-4-methyl-3-penten-1-yl]-1,4-naphthalenedione
Synonyms: (-)-Alkannin, NSC 94524
MF: C₁₆H₁₆O₅
FW: 288.3
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 250, 279, 516, 555 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Arnebia sinensis*



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

Laboratory Procedures

Alkannin is supplied as a crystalline solid. A stock solution may be made by dissolving the alkannin in the solvent of choice. Alkannin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of alkannin in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Alkannin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, alkannin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Alkannin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Alkannin is a naphthoquinone and enantiomer of shikonin (Item No. 14751) that has been found in *A. tinctoria* and has diverse biological activities.¹⁻⁵ It inhibits tumor-specific pyruvate kinase M2 (PKM2; IC₅₀ = 0.3 μM) with 20- and 10-fold selectivity over PKM1 and pyruvate kinase L (PKL), respectively.¹ Alkannin inhibits proliferation of HCT116 and SW480 colorectal cancer cells with IC₅₀ values of 2.38 and 4.53 μM, respectively.² It halts the cell cycle at the G₁ phase and induces apoptosis in HCT116 cells when used at a concentration of 3 μM. Alkannin (1 μM) increases levels of Hsp70 in untreated and UVB-irradiated HaCaT cells as well as inhibits UVB-induced DNA fragmentation and caspase-3 activity in HaCaT cells.³ It is active against methicillin-sensitive and -resistant *S. aureus* (MICs = 6.25 μg/ml) as well as vancomycin-sensitive and -resistant *E. faecalis* (MICs = 50 and 25 μg/ml, respectively).⁴ Alkannin scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals in a cell-free assay (EC₅₀ = 22 ppm).⁵

References

1. Chen, J., Xie, J., Jiang, Z., et al. *Oncogene* **30**(42), 4297-4306 (2011).
2. Huu Tung, N., Du, G.-J., Wang, C.-Z., et al. *Phytother. Res.* **27**(1), 66-70 (2013).
3. Yoshihisa, Y., Hassan, M.A., Furusawa, Y., et al. *PLoS One* **7**(10), e47903 (2012).
4. Shen, C.-C., Syu, W.-J., Li, S.-Y., et al. *J. Nat. Prod.* **65**(12), 1857-1862 (2002).
5. Assimopoulou, A.N. and Papageorgiou, V.P. *Phytother. Res.* **19**(2), 141-147 (2005).

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, 12/15/2022

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