

PRODUCT INFORMATION



Ilimaquinone

Item No. 15124

CAS Registry No.: 71678-03-0
Formal Name: 3-[[[(1R,2S,4aS,8aS)-decahydro-1,2,4a-trimethyl-5-methylene-1-naphthalenyl]methyl]-2-hydroxy-5-methoxy-2,5-cyclohexadiene-1,4-dione

MF: C₂₂H₃₀O₄

FW: 358.5

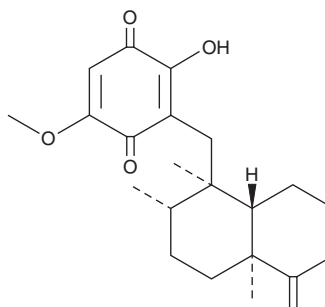
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Anthelia* sp.



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

Laboratory Procedures

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

Description

Ilimaquinone is a natural sesquiterpene quinone that has antimicrobial, anti-HIV, anti-mitotic, and anti-inflammatory properties.¹ In mammalian cells, 25 μM ilimaquinone reversibly induces vesiculation of Golgi membranes, blocking the secretory pathway.^{2,3} It inhibits the conversion of S-adenosylhomocysteine (SAH) to homocysteine by SAH hydrolase (IC₅₀ = 40 μM).^{1,4} Ilimaquinone also inhibits DNA polymerase β and dual specificity phosphatase Cdc25B (IC₅₀ = 45.2 and 92 μM, respectively) and, at 10 μM, activates gene expression through hypoxia-inducible factor-1.^{5,6}

References

1. Radeke, H.S., Digits, C.A., Casaubon, R.L., *et al.* Interactions of (-)-ilimaquinone with methylation enzymes: Implications for vesicular-mediated secretion. *Chem. Biol.* **6(9)**, 639-647 (1999).
2. Takizawa, P.A., Yucel, J.K., Veit, B., *et al.* Complete vesiculation of golgi membranes and inhibition of protein transport by a novel sea sponge metabolite, ilimaquinone. *Cell* **73(6)**, 1079-1090 (1993).
3. Cruciani, V., Leithe, E., and Mikalsen, S.O. Ilimaquinone inhibits gap-junctional communication prior to Golgi fragmentation and block in protein transport. *Exp. Cell Res.* **287(1)**, 130-142 (2003).
4. Kim, B.G., Chun, T.G., Lee, H.Y., *et al.* A new structural class of S-adenosylhomocysteine hydrolase inhibitors. *Bioorg. Med. Chem.* **17(18)**, 6707-6714 (2009).
5. Cao, S., Gao, Z., Thomas, S.J., *et al.* Marine sesquiterpenoids that inhibit the lyase activity of DNA polymerase β. *J. Nat. Prod.* **67(10)**, 1716-1718 (2004).
6. Du, L., Zhou, Y.D., and Nagle, D.G. Inducers of hypoxic response: Marine sesquiterpene quinones activate HIF-1. *J. Nat. Prod.* **76(6)**, 1175-1181 (2013).

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, 11/15/2022

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