

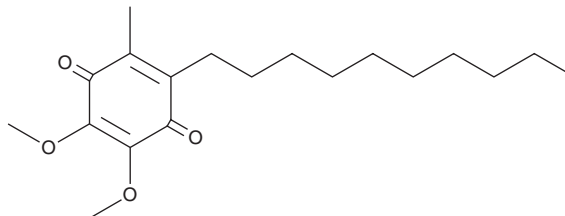
PRODUCT INFORMATION



Decylubiquinone

Item No. 21027

CAS Registry No.: 55486-00-5
Formal Name: 2-decyl-5,6-dimethoxy-3-methyl-2,5-cyclohexadiene-1,4-dione
MF: C₁₉H₃₀O₄
FW: 322.4
Purity: ≥98%
UV/Vis.: λ_{max}: 277, 403 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Decylubiquinone is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of XX in these solvents is approximately 10 mg/ml.

Decylubiquinone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of decylubiquinone should be diluted with the aqueous buffer of choice. Decylubiquinone has a solubility of approximately 0.3 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Decylubiquinone is an analog of ubiquinone (coenzyme Q₁₀; Item No. 11506). It blocks reactive oxygen species production in response to glutathione depletion and inhibits activation of the mitochondrial permeability transition.¹ Decylubiquinone inhibited opening of the mitochondrial permeability transition pore in response to calcium overload in rat liver mitochondria in different cell lines at concentrations between 50-100 μM.²

References

1. Armstrong, J.S., Whiteman, M., Rose, P., *et al.* The Coenzyme Q₁₀ analog decylubiquinone inhibits the redox-activated mitochondrial permeability transition: Role of mitochondrial complex III. *J. Biol. Chem.* **278(49)**, 49079-49084 (2003).
2. Devun, F., Walter, L., Belliere, J., *et al.* Ubiquinone analogs: A mitochondrial permeability transition pore-dependent pathway to selective cell death. *PLoS One* **5(7)**, (2010).

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, 02/09/2018

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