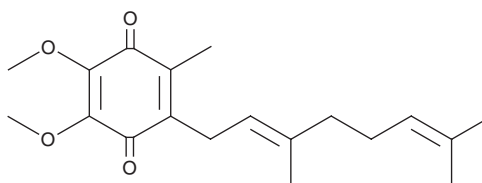


PRODUCT INFORMATION



Coenzyme Q₂ Item No. 17327

CAS Registry No.: 606-06-4
Formal Name: 2-[(2E)-3,7-dimethyl-2,6-octadien-1-yl]-5,6-dimethoxy-3-methyl-2,5-cyclohexadiene-1,4-dione
Synonyms: CoQ₂, Ubiquinone-2, Ubiquinone Q₂
MF: C₁₉H₂₆O₄
FW: 318.4
Purity: ≥95%
UV/Vis.: λ_{max}: 274 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

Coenzyme Q₂ 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 ethanol and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of coenzyme Q₂ in these solvents is approximately 10 mg/ml.

Coenzyme Q₂ is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Coenzyme Q₂ (CoQ₂) is a biosynthetic precursor to the electron transport chain cofactor CoQ₁₀ (Item No. 11506) and an inhibitor of mitochondrial complex I, also known as NADH dehydrogenase.^{1,2} It is an electron acceptor that accepts electrons from mitochondrial complex I. CoQ₂ inhibits the proliferation of BALL-1 human B cell acute lymphoblastic leukemia cells (IC₅₀ = 20 μM) and induces apoptosis in Huh7 and HepG2 hepatoma cells when used at a concentration of 100 μM.³ It increases the production of reactive oxygen species (ROS) in BALL-1 cells but inhibits ROS production in mitochondrial membranes derived from bovine heart when used in combination with the mitochondrial complex I inhibitors rotenone (Item No. 13995), piericidin A (Item No. 15379), or rolliniastatin 1 or -2.^{1,3} CoQ₂ increases the production of hydrogen peroxide in isolated rabbit heart mitochondria and decreases it in isolated rabbit liver mitochondria.⁴

References

1. Fato, R., Bergamini, C., Bortolus, M., *et al.* *Biochim. Biophys. Acta.* **1787(5)**, 384-392 (2009).
2. Ziegler, D. and Rieske, J.S. *Arch. Biochem. Biophys.* **97**, 231-235 (1962).
3. Esaka, Y., Nagahara, Y., Hasome, Y., *et al.* *Biochim. Biophys. Acta* **1724(1-2)**, 49-58 (2005).
4. Gharib, A., De Paulis, D., Li, B., *et al.* *J. Mol. Cell. Cardiol.* **52(5)**, 1091-1095 (2012).

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, 10/07/2022

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