

# PRODUCT INFORMATION



## Coenzyme Q<sub>9</sub> Item No. 16866

**CAS Registry No.:** 303-97-9  
**Formal Name:** 2,3-dimethoxy-5-methyl-6-(3,7,11,15,19,23,27,31,35-nonamethyl-2E,6E,10E,14E,18E, 22E,26E,30E,34-hexatriacontanonaen-1-yl)-2,5-cyclohexadiene-1,4-dione

**Synonyms:** CoQ<sub>9</sub>, NSC 226993, Ubiquinone-9, Ubiquinone Q<sub>9</sub>

**MF:** C<sub>54</sub>H<sub>82</sub>O<sub>4</sub>

**FW:** 795.20

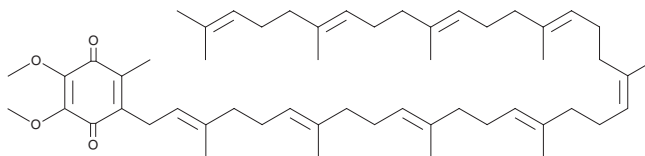
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 210, 275 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.

### Laboratory Procedures

Coenzyme Q<sub>9</sub> (CoQ<sub>9</sub>) is supplied as a crystalline solid. A stock solution may be made by dissolving the CoQ<sub>9</sub> in the solvent of choice, which should be purged with an inert gas. CoQ<sub>9</sub> is soluble in organic solvents such as ethanol, chloroform, and dimethyl formamide. The solubility of CoQ<sub>9</sub> in these solvents is approximately 0.1, 15, and 10 mg/ml, respectively.

### Description

CoQ<sub>9</sub> is a mitochondrial ubiquinone with a role in the electron transport chain.<sup>1</sup> It reduces the incidence of ischemia and reperfusion-induced arrhythmia and improves post-ischemic heart rate, coronary flow, aortic flow, and left ventricular developed pressure in isolated guinea pig hearts when used at a dose of 5 mg/kg.<sup>1</sup> CoQ<sub>9</sub> (5 mg/kg) also reduces ischemia and reperfusion-induced myocardial infarct size and the number of apoptotic cardiomyocytes in the same model.

### Reference

1. Lekli, I., Das, S., Das, S., *et al.* Coenzyme Q<sub>9</sub> provides cardioprotection after converting into coenzyme Q<sub>10</sub>. *J. Agric. Food Chem.* **56**(13), 5331-5337 (2008).

#### 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, 10/13/2022

#### 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