

# Decylubiquinone

## MPTP inhibitor

Decylubiquinone is an analogue of ubiquinone. Decylubiquinone inhibits the  $\text{Ca}^{2+}$ -dependent mitochondrial permeability transition pore (MPTP) without inhibiting respiration. Maximum effect on  $\text{Ca}^{2+}$  retention capacity of rat liver mitochondria can be achieved with decylubiquinone at  $100\mu\text{M}$ . This inhibitory effect of decylubiquinone on MPTP can be counteracted by ubiquinone-5 (Prod. No. ALX-270-294). Decylubiquinone also inhibits etoposide-induced apoptosis in L929 cells at  $5\mu\text{M}$ .

Citations: 7

[View Online »](#)

## Ordering Information

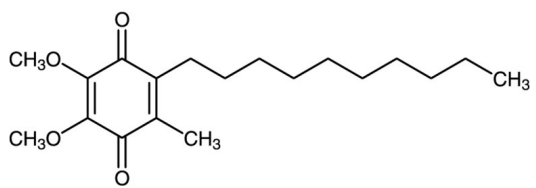
[Order Online »](#)

BML-CM115-0010	10mg
BML-CM115-0050	50mg

## Manuals, SDS & CofA

[View Online »](#)

- MPTP inhibitor used in cellular metabolism studies



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	2,3-Dimethoxy-5-methyl-6-decyl-1,4-benzoquinone
Appearance	Neat yellow to orange colored oil.
CAS	55486-00-5
Couple Target	MPTP
Couple Type	Inhibitor
Formula	$C_{19}H_{30}O_4$
MW	322.4
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (10mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)