

Ubiquinone-10

MPTP inhibitor

Inhibitor of the mitochondrial permeability transition pore (MPTP). This inhibitory effect can be counteracted by ubiquinone-5 (Prod. No. ALX-270-294).

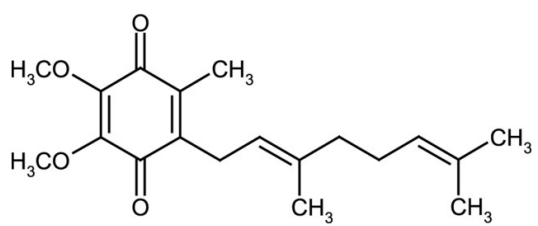
Ordering Information

[Order Online »](#)

ALX-270-295-M002	2mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Coenzyme Q2, 2,3-Dimethoxy-5-methyl-geranyl-1,4-benzoquinone

Appearance Red liquid.

CAS 606-06-4

Couple Target MPTP

Couple Type Inhibitor

Formula $C_{19}H_{26}O_4$

Identity Identity determined by 1H-NMR.

MW 318.4

Purity ≥90% (HPLC)

Solubility Soluble in acetone (50mg/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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)