

Idebenone

Quinone antioxidant

Idebenone is a quinone antioxidant chemically related to ubiquinone. It is a free radical scavenger that protects mitochondrial membranes from lipid peroxidation. It improves overall brain function in various animal models and protective effect on astrocytes is due in part *via* increased NGF production.

Citations: 3

[View Online »](#)

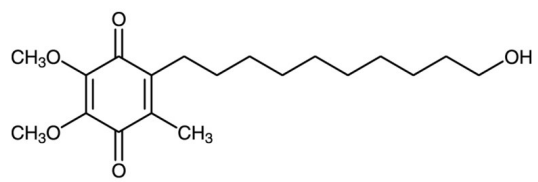
Ordering Information

[Order Online »](#)

BML-FR114-0020	20mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	2-(10-Hydroxydecyl)-5,6-dimethoxy-3-methyl-2,5-cyclohexadiene-1,4-dione
Appearance	Orange solid.
CAS	58186-27-9
Formula	$C_{19}H_{30}O_5$
MW	338.4
Purity	≥98% (HPLC)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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