

DMNQ

Inducer of intracellular ROS generation

Redox-cycling agent that induces intracellular superoxide anion formation and, depending on the concentration, induces cell proliferation, apoptosis or necrosis. DMNQ does not react with free thiol groups, is non-alkylating and adduct-forming in contrast to other quinones. Thus, DMNQ is a valuable tool for the generation of reactive oxygen species (ROS) in order to study the role of ROS in cell toxicity, apoptosis and necrosis.

Citations: 9

[View Online »](#)

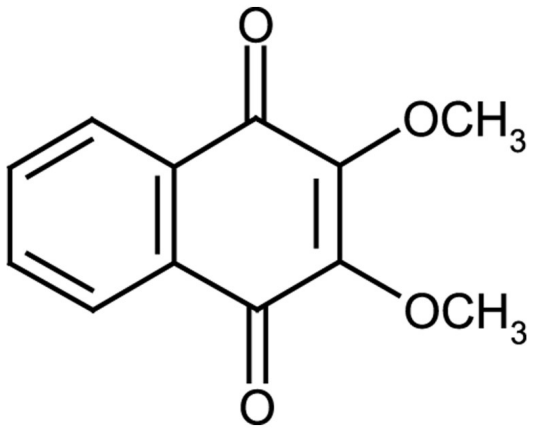
Ordering Information

[Order Online »](#)

ALX-420-027-M005	5mg
ALX-420-027-M010	10mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	2,3-Dimethoxy-1,4-naphthoquinone
Appearance	Yellow crystalline solid.
CAS	6956-96-3
Couple Target	ROS
Couple Type	Activator
Formula	$C_{12}H_{10}O_4$
Identity	Determined by NMR.
MW	218.2
Purity	≥99% (TLC)
Solubility	Soluble in acetone, DMSO or methanol.

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