

DMPO

Spin trapping reagent

Cell permeable hydrophilic spin trap for both *in vivo* and *in vitro* studies of superoxide, O-, C-, S- and N- centered free radicals.

Citations: 48

[View Online »](#)

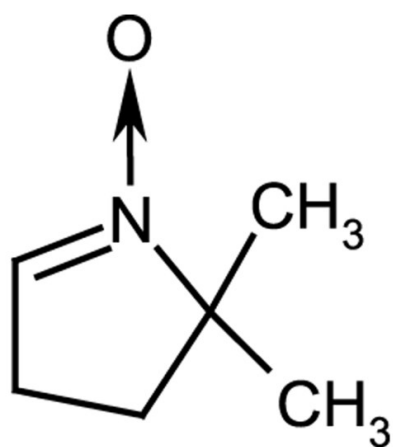
Ordering Information

[Order Online »](#)

ALX-430-090-M500	500mg
ALX-430-090-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture. Keep under inert gas. Hygroscopic.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5,5-Dimethyl-1-pyrroline-N-oxide, 2,2-Dimethyl-3,4-dihydro-2H-pyrrole 1-oxide
Appearance	Pale yellow solid or oil.
CAS	3317-61-1
Formula	$C_6H_{11}NO$
MW	113.2
Purity	>98% (HPLC)
Solubility	Soluble in DMSO (up to 25 mg/ml) or Ethanol (up to 25 mg/ml).

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