

DEPMPO

Spin trapping reagent

Most efficient spin trap for the *in vitro* and *in vivo* detection of O-, N-, S- and C-centered free radicals. Has a longer life-time than DMPO (Prod. No. ALX-430-090). Can distinguish between superoxide-dependent and independent mechanisms that lead to the hydroxyl radical. Less lipophilic ($K_p=0.16$) than DIPPMPO (Prod. No. ALX-430-119).

Citations: 36

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-430-093-M050	50mg
ALX-430-093-M500	500mg

Manuals, SDS & CofA

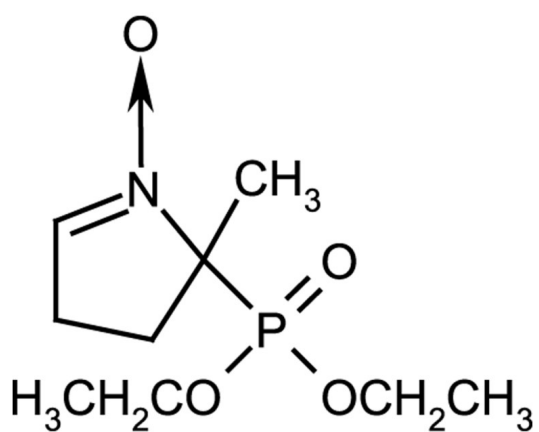
[View Online »](#)



Spectrum of the (O²⁻)-DEPMPO adduct.



Spectrum of the (HO)-DEPMPO adduct.



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Unstable in solution. Reconstitute just prior to use.
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-(Diethoxyphosphoryl)-5-methyl-1-pyrroline-N-oxide, 2-Diethylphosphono-2-methyl-3,4-dihydro-2H-pyrrole 1-oxide, (2-Methyl-3,4-dihydro-1-oxide-2H-pyrrol-2-yl)diethylphosphonate
Appearance	Colorless to light yellow oil.
CAS	157230-67-6
Formula	$C_9H_{18}NO_4P$
MW	235.2
Purity	≥99% (TLC)
Solubility	Soluble in water, methylene chloride or 100% ethanol.



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-fr@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