

EMPO

Spin trapping reagent

Beside DEPMPO (Prod. No. ALX-430-093), 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).

Citations: 16

[View Online »](#)

Ordering Information

[Order Online »](#)

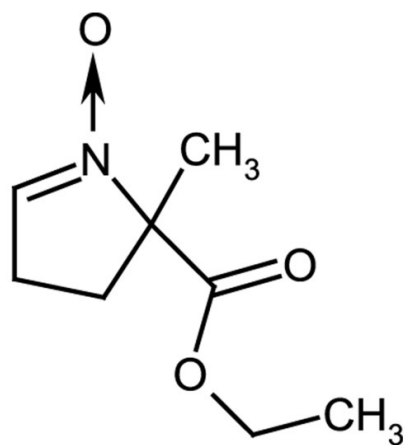
ALX-430-098-M010	10mg
ALX-430-098-M050	50mg

Manuals, SDS & CofA

[View Online »](#)



Spectrum of the (O²⁺) – EMPO adduct.



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C.
Handling	Protect from light. Hygroscopic.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-Ethoxycarbonyl-2-methyl-3,4-dihydro-2H-pyrrole-1-oxide
Appearance	Colorless to light yellow oil.
Formula	$C_8H_{13}NO_3$
Identity	Determined by NMR.
MW	171.2
Purity	≥95% (HPLC)
Solubility	Soluble in water, DMSO, ethanol or dichloromethane.

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-eu@enzolifesciences.com)

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