

TEMPONE

Spin trapping reagent

Spin trap, useful for both *in vitro* and *in vivo* experiments.

Citations: 8

[View Online »](#)

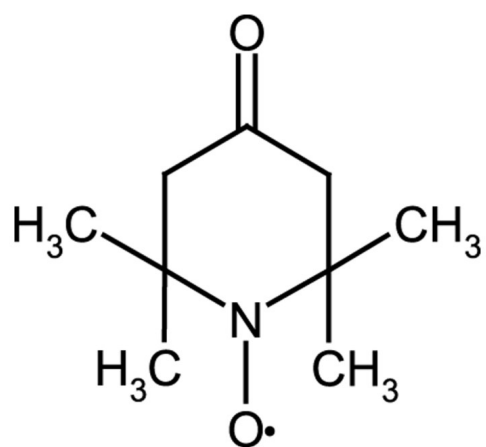
Ordering Information

[Order Online »](#)

ALX-430-079-M250	250mg
------------------	-------

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	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-Oxo-2,2,6,6-tetramethylpiperidine-1-oxyl, 4-Oxo-TEMPO
Appearance	Orange crystalline powder.
CAS	2896-70-0
Formula	$C_9H_{16}NO_2$
Identity	Determined by IR.
MW	170.2
Purity	≥98% (Assay)
Solubility	Soluble in acetone, DMSO, diethylether, 100% ethanol or water.

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-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