

mito-TEMPO-H

Spin probe

Spin probe for detection of mitochondrial superoxide.
Mitochondria-targeted antioxidant with superoxide and alkyl radical scavenging properties, which may be used for both *in vivo* and *in vitro* experiments.

Citations: 7

[View Online »](#)

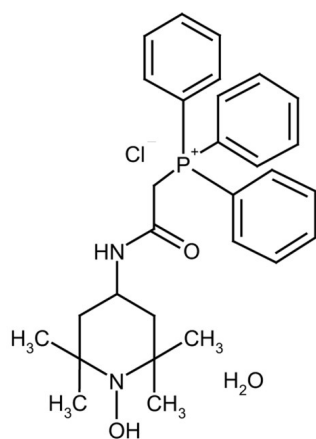
Ordering Information

[Order Online »](#)

ALX-430-171-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



H₂O

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-hydroxy-4-[2-triphenylphosphonio)-acetamido]-2,2,6,6-tetramethylpiperidine, 1-Hydroxy-2,2,6,6-tetramethyl-4-[2-(triphenylphosphonio)acetamido]piperidinium dichloride
Appearance	Colorless to pink.
Application Notes	Mito-TEMPO-H can be used for superoxide detection.
Formula	$C_{29}H_{37}N_2O_2Cl_2P \cdot H_2O$
MW	511.1 . 18.0
Purity	≥95% (EPR)
Solubility	Soluble in water, 100% ethanol or methanol.

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-
be@enzolifesciences.com](mailto:info-be@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)