

Mito-TEMPO

Mitochondria-targeted antioxidant

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

Citations: 66

[View Online »](#)

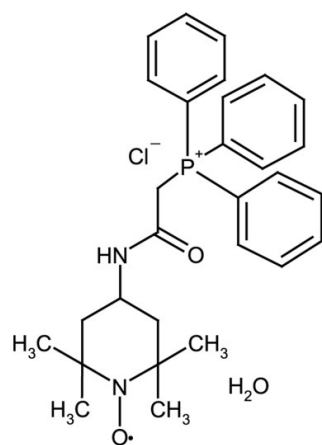
Ordering Information

[Order Online »](#)

ALX-430-150-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in water, saline, ethanol or DMSO are stable at +4°C for 1-2 weeks. DMSO stock solution can be stored also at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(2-(2,2,6,6-Tetramethylpiperidin-1-oxyl-4-ylamino)-2-oxoethyl)triphenylphosphonium chloride
Appearance	Colorless film to pale orange solid.
CAS	1334850-99-5
Formula	$C_{29}H_{35}N_2O_2ClP$
Identity	Determined by MS.
MW	510.03
Purity	≥95% (TLC)
Solubility	Soluble in water, DMSO, 100% ethanol or methanol.

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