

CMH . hydrochloride

Spin probe reagent

Analog of CPH (Prod. No. ALX-430-078) with higher cell permeability and 3-fold better reactivity with superoxide. Has been used to study intracellular production of superoxide radicals in plasma, endothelial cells and isolated hearts.

Citations: 28

[View Online »](#)

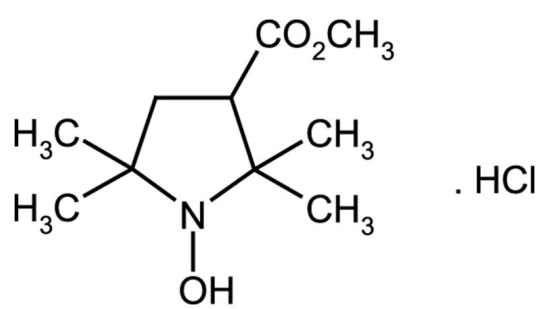
Ordering Information

[Order Online »](#)

ALX-430-117-M010	10mg
ALX-430-117-M050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Prepare fresh solutions daily before use and store 4-6 hours on ice under a nitrogen/argon flow. Oxygen will lead to oxidation and therefore it is strongly recommended to dissolve the probe in water and not in organic solvents such as DMSO, ethanol or methanol. Oxygen is 10 times more soluble in the latter solvents. Freezing solutions will not lead to better results, because mixing with air after melting the probe solutions will result in faster oxidation. Also, it is strongly recommended to store the product and its solutions in glass vials.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-Hydroxy-3-methoxycarbonyl-2,2,5,5-tetramethylpyrrolidine . HCl
Appearance	White to light-yellow crystalline powder.
Formula	$C_{10}H_{19}NO_3 \cdot HCl$
MW	201.3 . 36.5
Purity	≥99% (EPR)
Solubility	Soluble in DMSO, 100% ethanol, methanol or water.

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)