

# CPH . hydrochloride

## Spin probe reagent

Effective, cell permeable and non-toxic spin probe for the detection of superoxide radical and peroxynitrite, both *in vitro* and *in vivo*. Resistant to reduction by vitamin C and thiols.

Citations: 14

[View Online »](#)

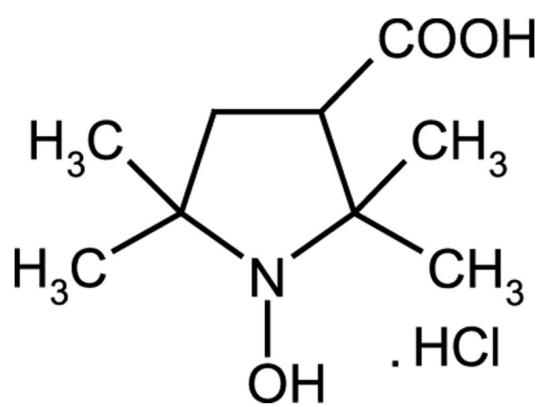
## Ordering Information

[Order Online »](#)

ALX-430-078-M050	50mg
ALX-430-078-M250	250mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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 o
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	1-Hydroxy-3-carboxy-2,2,5,5-tetramethylpyrrolidine . HCl
<b>Appearance</b>	White to off-white solid.
<b>CAS</b>	164364-89-0
<b>Formula</b>	$C_9H_{17}NO_3 \cdot HCl$
<b>MW</b>	223.7
<b>Purity</b>	≥99% (EPR)
<b>Solubility</b>	Soluble in DMSO, 100% ethanol, methanol or water.



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)