

CAT1H . hydrochloride

Spin probe

Cyclic hydroxylamine spin probe for *in vivo*, intraperitoneal and intravenous injections. Lifetime of more than 4 hours in blood plasma, in cells and tissue. Allows quantitative measurements of extracellular $O_2^{\bullet-}$ and quantification of intracellular $O_2^{\bullet-}$ in cells and tissue samples and *in vivo* $O_2^{\bullet-}$ detection.

Citations: 7

[View Online »](#)

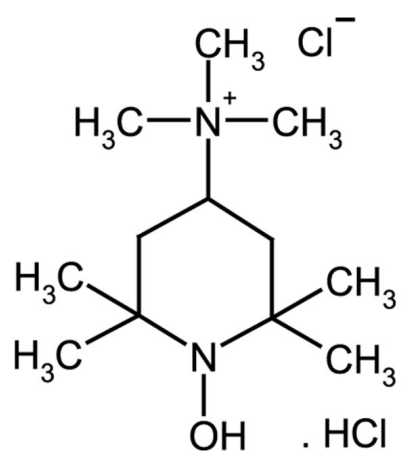
Ordering Information

[Order Online »](#)

ALX-430-131-M050	50mg
ALX-430-131-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 moisture.
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-2,2,6,6-tetramethylpiperidin-4-yl-trimethylammonium chloride . HCl
Appearance	Colorless crystalline solid.
Formula	$C_{12}H_{26}N_2O \cdot 2HCl$
MW	214.4 . 72.9
Purity Detail	Paramagnetic impurities ≤0.01% (EPR).
Solubility	Soluble in water, methanol or 100% ethanol.

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)