

PPH . hydrochloride

Spin trapping reagent

Effective, non-cell permeable and non-toxic analog of CPH for the detection of superoxide radicals. Resistant against reduction of vitamin C. For cell permeable analog see CPH (Prod. No. ALX-430-078).

Citations: 6

[View Online »](#)

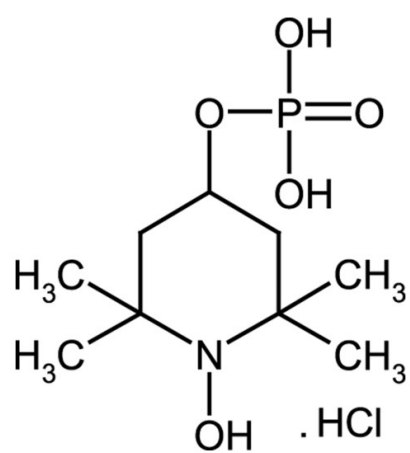
Ordering Information

[Order Online »](#)

ALX-430-080-M010	10mg
ALX-430-080-M050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-Hydroxy-4-phosphono-oxy-2,2,6,6-tetramethylpiperidine . HCl

Appearance White to off-white solid.

Formula $C_9H_{20}NO_5P \cdot HCl$

MW 253.2 . 36.5

Purity ≥99% (EPR)

Purity Detail ≤0.1% paramagnetic impurities (EPR)

Solubility Soluble in DMSO, 100% ethanol, methanol or water.

Technical Info / Product Notes The stabilized acid form of PPH is not well soluble at neutral pH (~0.1-0.01%). It is recommended to adjust pH to 7.4 for higher solubility. Dissolve PPH in oxygen-free buffers at neutral pH and use 1M NaOH to adjust pH to 7.4. Solubility is 1% in water, 0.8% in methanol or 0.5% in ethanol at pH 7.4. Solubility is 3% in DMSO at pH 7.4 (first add NaOH to PPH crystals then DMSO). Higher concentrations require further pH adjustment using NaOH.

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@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)