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 »

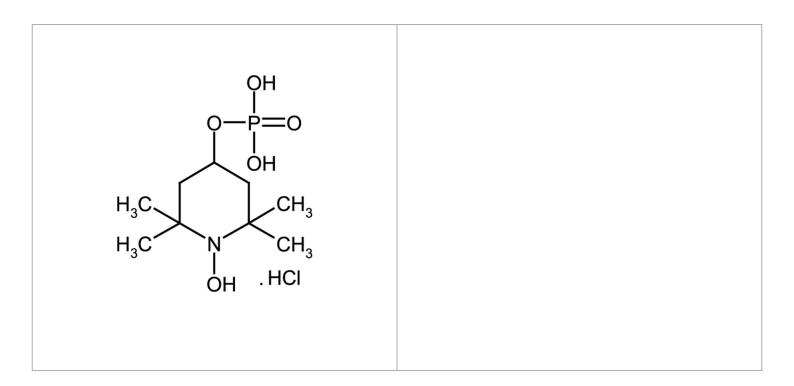
Ordering Information

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

. HCI

Appearance White to off-white solid.

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 NotesThe 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

