# **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** 

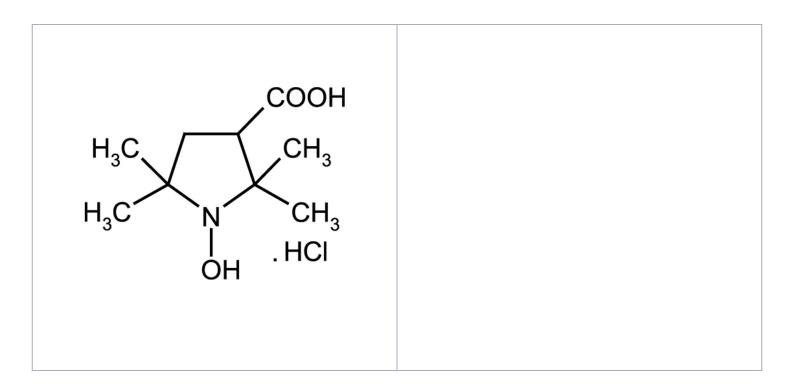
Order Online »

ALX-430-078-M050

50mg

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 o

**Long Term Storage** -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name 1-Hydroxy-3-carboxy-2,2,5,5-tetramethylpyrrolidine . HCl

**Appearance** White to off-white solid.

**CAS** 164364-89-0

Formula  $C_9H_{17}NO_3$  . HCI

**MW** 223.7

**Purity** ≥99% (EPR)

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

Last modified: August 15, 2025