

BMPO (high purity)

Spin trapping reagent

BMPO is a cyclic nitroxide spin trap for the specific *in vivo* or *in vitro* detection of short-lived hydroxyl, superoxide, and thiyl radicals. BMPO forms distinguishable adducts that can be measured by EPR spectroscopy. Unlike with DMPO (Prod. No. ALX-430-090), the superoxide adduct does not decay into a hydroxyl adduct and has a much longer half-life in cells ($t_{1/2}=23\text{min}$).

Citations: 14

[View Online »](#)

Ordering Information

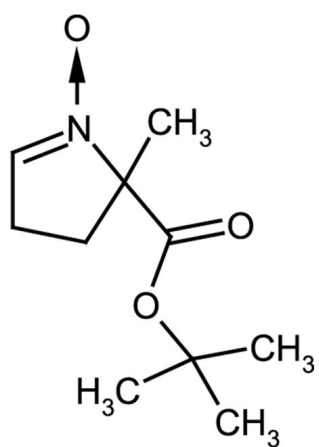
[Order Online »](#)

ALX-430-141-M010	10mg
ALX-430-141-M050	50mg

Manuals, SDS & CofA

[View Online »](#)

- Highly pure spin trapping reagent
- Contains very low paramagnetic impurities and does not require further purification
- Comes as a crystalline solid for a longer shelf-life



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-tert-Butoxycarbonyl-5-methyl-1-pyrroline-N-oxide
Appearance	White to pinkish crystalline solid.
CAS	387334-31-8
Formula	$C_{10}H_{17}NO_3$
Identity	Identity determined by MS and NMR.
MW	199.3
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (25mg/ml), 100% ethanol, 100% methanol, and water.
Source	Synthetic

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-
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)