

PBN

Spin trapping reagent

Cell permeable spin trap for both *in vivo* and *in vitro* studies.

Citations: 21

[View Online »](#)

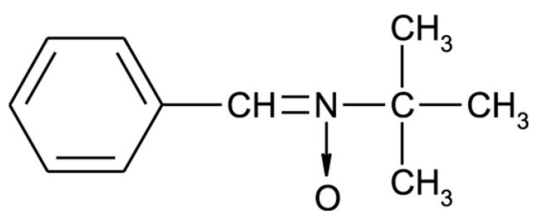
Ordering Information

[Order Online »](#)

ALX-430-082-G001	1g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-tert-Butyl- α -phenylnitrone
Appearance	White to off-white crystalline solid.
CAS	3376-24-7
Formula	$C_{11}H_{15}NO$
MI	14: 7056
MW	177.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO or 100% ethanol; slightly soluble in water.
Technical Info / Product Notes	For selected literature references on spin trapping and beer please click here.

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)