

POBN (high purity)

Spin trapping reagent

Cell permeable hydrophilic spin trap for both *in vivo* and *in vitro* studies. Water soluble analog of PBN (Prod. No. ALX-430-082). Low paramagnetic impurities. No further purification required.

Citations: 8

[View Online »](#)

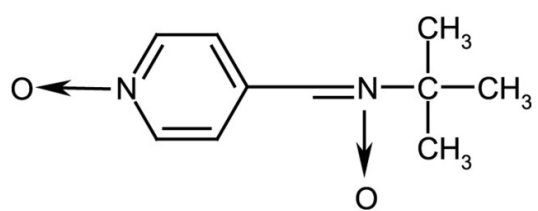
Ordering Information

[Order Online »](#)

ALX-430-091-M500	500mg
ALX-430-091-G001	1g

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 α -(4-Pyridyl 1-oxide)-N-tert-butylnitron

Appearance White to off-white or colorless crystalline solid.

CAS 66893-81-0

Formula $C_{10}H_{14}N_2O_2$

MW 194.2

Purity $\geq 99\%$ (HPLC)

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

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