

MCPIO

Spin trapping reagent

Cell permeable, stable spin trap. Alternative to PBN (Prod. No. ALX-430-082).

Citations: 2

[View Online »](#)

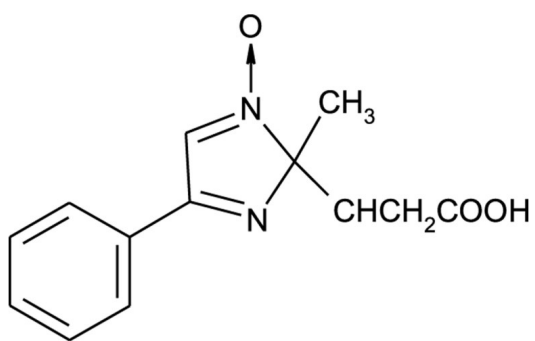
Ordering Information

[Order Online »](#)

ALX-430-083-M010	10mg
------------------	------

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 2-(2-Carboxyethyl)-2-methyl-4-phenyl-2H-imidazole-1-oxide

Appearance Off-white powder.

Formula $C_{13}H_{14}N_2O_3$

MW 246.3

Purity ≥98% (HPLC)

Solubility Soluble in water, 100% ethanol or DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com