

MnTBAP chloride

SOD mimetic. Peroxynitrite scavenger.

Cell permeable superoxide dismutase (SOD) mimetic. Potent inhibitor of peroxynitrite-induced oxidative reactions (peroxynitrite scavenger), but not a scavenger of nitric oxide (NO). Inhibits neuronal apoptosis.

Citations: 10

[View Online »](#)

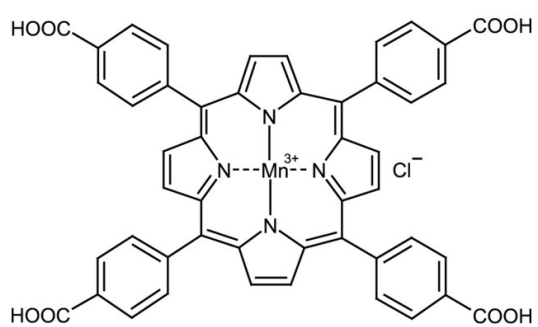
Ordering Information

[Order Online »](#)

ALX-430-069-M010	10mg
ALX-430-069-M050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 2 years after receipt when stored at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Black crystalline powder.
Formula	$C_{48}H_{28}MnN_4O_8Cl$
MW	879.2
Purity	≥95% (TLC)
Solubility	Dissolve at 0.1M in aqueous base, then add buffer to lower pH to pH 7.0 or higher as needed.

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)