

TPEN

Heavy metal chelator

A cell permeable and extremely high-affinity chelator of heavy metals, yet with low affinity for Mg^{2+} and Ca^{2+} . Lipid-soluble, crossing both natural and artificial membranes. TPEN can be used in conjunction with QUIN 2 (Prod. No. ENZ-52013) or QUIN 2/AM to reduce heavy-metal quenching of QUIN 2 fluorescence in cells. Ion specificity: $Zn^{2+} > Fe^{2+} > Mn^{2+} >> Ca^{2+} = Mg^{2+}$. Complex with Fe^{2+} catalyzes superoxide dismutation. Induces DNA fragmentation when chelated with Zn^{2+} and induces apoptosis in T cells.

Citations: 12

[View Online »](#)

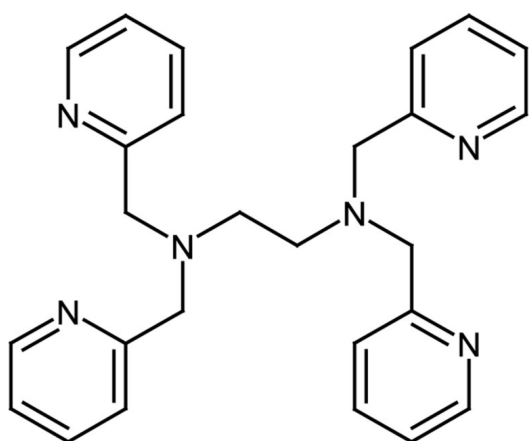
Ordering Information

[Order Online »](#)

ALX-450-011-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N,N,N',N'-Tetrakis-(2-pyridylmethyl)ethylenediamine

Appearance White to light yellow to brown solid.

CAS 16858-02-9

Formula $C_{26}H_{28}N_6$

MW 424.5

Purity ≥97% (Nonaqueous Titration)

Solubility Soluble in dimethyl formamide, DMSO, 100% ethanol or methanol.

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)