

Zinc(II) Protoporphyrin IX

Heme oxygenase inhibitor

Potent and selective inhibitor of heme oxygenase ($K_i=3\text{nM}$), the enzyme that generates carbon monoxide (CO). Inhibits soluble guanylyl cyclase. Produces a time- and concentration-dependent inactivation of all three isoforms of nitric oxide synthase ($\text{IC}_{50}=0.8\mu\text{M}$, $4.0\mu\text{M}$, and $5.0\mu\text{M}$ for nNOS (NOS I), iNOS (NOS II) and eNOS (NOS III), respectively). Does not cross the blood-brain barrier. An useful tool to elucidate the role of CO in signal transduction and as a neurotransmitter.

Citations: 24

[View Online »](#)

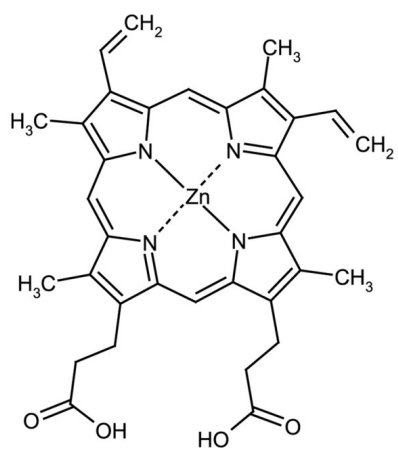
Ordering Information

[Order Online »](#)

ALX-430-049-M025	25mg
ALX-430-049-M100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 3 years after receipt when stored at -20°C. Make solutions fresh daily. Product may be unstable in solution if exposed to light.
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Dark solid.
CAS	15442-64-5
Couple Target	Heme oxygenase
Couple Type	Inhibitor
Formula	$C_{34}H_{32}N_4O_4Zn$
Identity	Identity determined by UV and NMR.
MW	626.0
Purity	≥95% (TLC)
Solubility	Soluble in pyridine, dimethyl formamide or DMSO. Colloidal in aqueous base; slightly soluble in aqueous base/organic solvent mixtures.



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)