

# 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 »](#)

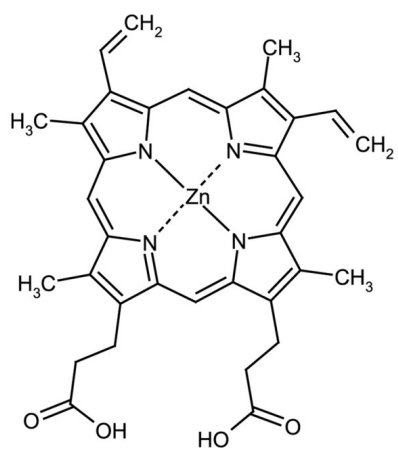
## Ordering Information

[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.

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-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-fr@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)