

Tin Protoporphyrin IX . dichloride

Heme oxygenase inhibitor

Potent inhibitor of heme oxygenase.

Citations: 7

[View Online »](#)

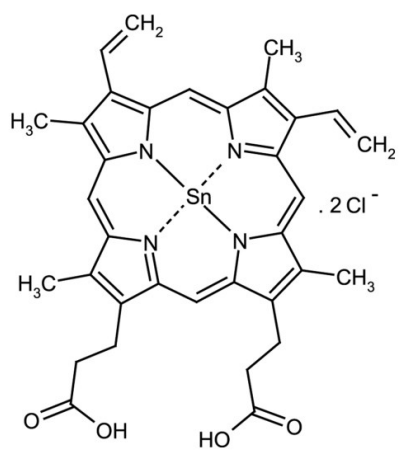
Ordering Information

[Order Online »](#)

ALX-430-051-M005	5mg
ALX-430-051-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Dark red to violet solid.
CAS	14325-05-4
Couple Target	Heme oxygenase
Couple Type	Inhibitor
Formula	$C_{34}H_{32}N_4O_4Sn \cdot 2Cl$
Identity	Identity determined by NMR and UV VIS.
MW	750.3
Purity	≥95% (HPLC)
Solubility	Soluble in aqueous base.

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-
fr@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)