

Copper(II) Protoporphyrin IX (free acid)

Negative control for the heme oxygenase inhibitor zinc protoporphyrin IX (Prod. No. ALX-430-049).

Citations: 3

[View Online »](#)

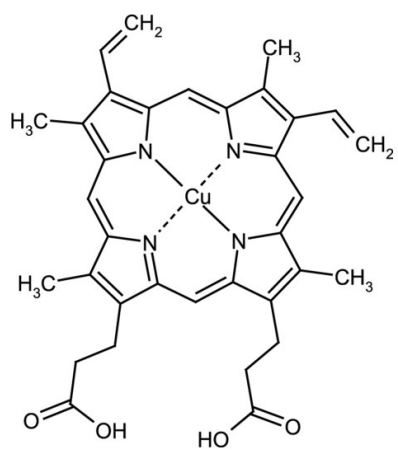
Ordering Information

[Order Online »](#)

ALX-430-048-M025	25mg
------------------	------

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. 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	Red to brown solid.
Formula	$C_{34}H_{32}CuN_4O_4$
MW	624.2
Purity	≥95%
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-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)