

Cobalt (III) Protoporphyrin IX chloride

Heme oxygenase activator

The ability of cobalt ions to induce heme oxygenase (HO) activity in liver cells was first reported almost 30 years ago. Cobalt protoporphyrin IX (CoPP) was later shown to be a substrate for HO, now identified as the inducible isoform HO-1 (heat shock protein 32; hsp32). CoPP has been shown to down-regulate various cytochrome P450 isoforms and various mechanisms of action have been attributed to CoPP and other metalloporphyrins.

Citations: 11

[View Online »](#)

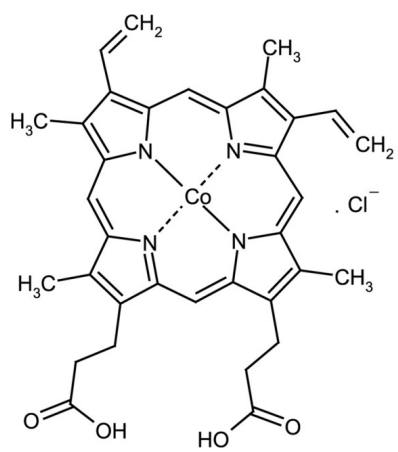
Ordering Information

[Order Online »](#)

ALX-430-076-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 and moisture.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Dark purple solid.

CAS 102601-60-5

Couple Target Heme oxygenase

Couple Type Activator

Formula $C_{34}H_{32}CoN_4O_4 \cdot Cl$

Identity Identity determined by UV/VIS and NMR.

MW 654.5

Purity ≥95% (TLC)

Solubility Soluble in basic aqueous solutions at pH between 8.5-9.0. Once in solution pH can be adjusted to physiological pH.



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)